



Biolistic® PDS-1000/He Particle Delivery System

**Catalog Numbers
165-2257 and
165-2250LEASE
to 165-2255LEASE**

BIO-RAD

Warranty and Regulatory Notices

Warranty Statement

This warranty may vary outside of the continental United States. Contact your local Bio-Rad office for the exact terms of your warranty.

Bio-Rad Laboratories warrants that the Biolistic PDS-1000/He system (catalog numbers 165-2257 and 165-2250LEASE to 165-2255LEASE) will be free from defects in material and workmanship, and will meet all performance specifications for the period of 1 year from the date of shipment. This warranty covers all parts and labor.

In the event that the instrument must be returned to the factory for repair under warranty, the instrument must be packed for return in the original packaging.

Bio-Rad shall not be liable for any incidental, special, or consequential loss, damage, or expense directly or indirectly arising from the use of the Biolistic PDS-1000/He system. Bio-Rad makes no warranty whatsoever in regard to products or parts furnished by third parties, such being subject to the warranty of their respective manufacturers. Service under this warranty shall be requested by contacting your nearest Bio-Rad office.

The following items are considered customer-installed consumables: fuses, microcarriers, macrocarriers, and rupture disks. These parts are not covered by this warranty. All customer-installed parts are warranted only to be free from defects in workmanship.

This warranty does not extend to any instruments or parts thereof that have been subject to misuse, neglect, or accident, or that have been modified by anyone other than Bio-Rad or that have been used in violation of Bio-Rad instructions.

The foregoing obligations are in lieu of all other obligations and liabilities including negligence and all warranties, of merchantability, fitness for a particular purpose or otherwise, expressed or implied in fact or by law, and state Bio-Rad's entire and exclusive liability and buyer's exclusive remedy for any claims or damages in connection with the furnishing of goods or parts, their design, suitability for use, installation, or operation. Bio-Rad will in no event be liable for any special, incidental, or consequential damages whatsoever, and Bio-Rad's liability under no circumstances will exceed the contract price for the goods for which liability is claimed.

Regulatory Notices

Important: This Bio-Rad instrument is designed and certified to meet EN55011, EN50082-1, and IEC 1010-1 requirements, which are internationally accepted electrical safety standards. Certified products are safe to use when operated in accordance with the instruction manual. This instrument should not be modified or altered in any way. Alteration of this instrument will:

- Void the manufacturer's warranty.

- Void the regulatory certifications.

- Create a potential safety hazard.

Note: This equipment has been tested and found to comply with the limits for a Class A digital device, pursuant to Part 15 of the FCC rules. These limits are designed to provide reasonable protection against harmful interference when the equipment is operated in a commercial environment. This equipment generates, uses, and can radiate radio frequency energy and, if not installed and used in accordance with the instruction manual, may cause harmful interference to radio communications. Operation of this equipment in a residential area is likely to cause harmful interference in which case the user will be required to correct the interference at his own expense.

Table of Contents

Section 1	Biolistic PDS-1000/He Particle Delivery System.....	1
1.1	Particle Delivery Technology	1
1.2	Overview of PDS-1000/He Particle Delivery System	1
1.3	Important Safety Information.....	3
1.4	Requirements for System Operation.....	3
Section 2	Product Description	5
2.1	Packing List.....	5
2.2	Identification of Unit Controls and Components	10
Section 3	Installation	13
3.1	Connecting the PDS-1000/He System to a Helium Source	13
3.2	Connecting the PDS-1000/He to a Vacuum Source.....	15
3.3	Power Cord/Voltage Regulator.....	16
Section 4	Operation of the PDS-1000/He Instrument	17
4.1	Preparation of System Components Prior to Bombardment.....	17
4.2	Performing a Bombardment.....	22
4.3	Removal of Residual Helium Pressure–Shut Down	30
Section 5	Selection and Adjustment of System Bombardment Parameters.....	30
5.1	Overview -Matrix of Variables: Cell Types, Settings and Conditions	30
5.2	Vacuum Level in Bombardment Chamber.....	31
5.3	Helium Pressure / Rupture Disk Selection	31
5.4	Solenoid Valve Adjustment	32
5.5	Vacuum Flow Rate Control Valves	32
5.6	Distance Between Rupture Disk and Macrocarrier	32
5.7	Distance Between Macrocarrier and Stopping Screen	32
5.8	Distance Between Stopping Screen and Target Shelf (microcarrier flight distance).....	32
5.9	Microcarrier Selection.....	33
5.10	Preparation of Biological Material for Bombardment	34
Section 6	Troubleshooting.....	37
6.1	Rupture Disk Bursts at Incorrect Pressure	37
6.2	Stopping Screen Forced Through Screen Support Ring	37
6.3	Excessive Gas Usage.....	38
6.4	Chamber Will Not Hold Vacuum	38
6.5	Sample Damage From Gas Pressure Wave	38
6.6	Unit Will Not Pressurize Gas Acceleration Tube.....	39
Section 7	Product Information.....	40
7.1	Biolistic System.....	40
7.2	Spare Parts	41
Section 8	Appendices	43
8.1	Cleaning the PDS-1000/He Device	43
8.2	Metal Case Version	43
8.3	Specifications	46
8.4	Performing a Bombardment—Quick Guide Tear-Out.....	47

Section 1

Introduction to Particle Delivery

1.1 Particle Delivery Technology

Biolistic particle delivery is a method of transformation that uses helium pressure to introduce DNA-coated microcarriers into cells. Microprojectile bombardment can transform such diverse targets as bacterial, fungal, insect, plant, and animal cells and intracellular organelles. Particle delivery is a convenient method for transforming intact cells in culture since minimal pre- or post-bombardment manipulation is necessary. In addition, this technique is much easier and faster to perform than the tedious task of micro-injection. Both stable and transient transformation are possible with the Biolistic particle delivery system.

1.2 Overview of PDS-1000/He Particle Delivery System

The Biolistic System

The Biolistic PDS-1000/He instrument uses pressurized helium to accelerate sub-cellular sized microprojectiles coated with DNA (or other biological material) over a range of velocities necessary to optimally transform many different cell types. The system consists of the bombardment chamber (main unit), connective tubing for attachment to vacuum source, and all components necessary for attachment and delivery of high pressure helium to the main unit (helium regulator, solenoid valve, and connective tubing).

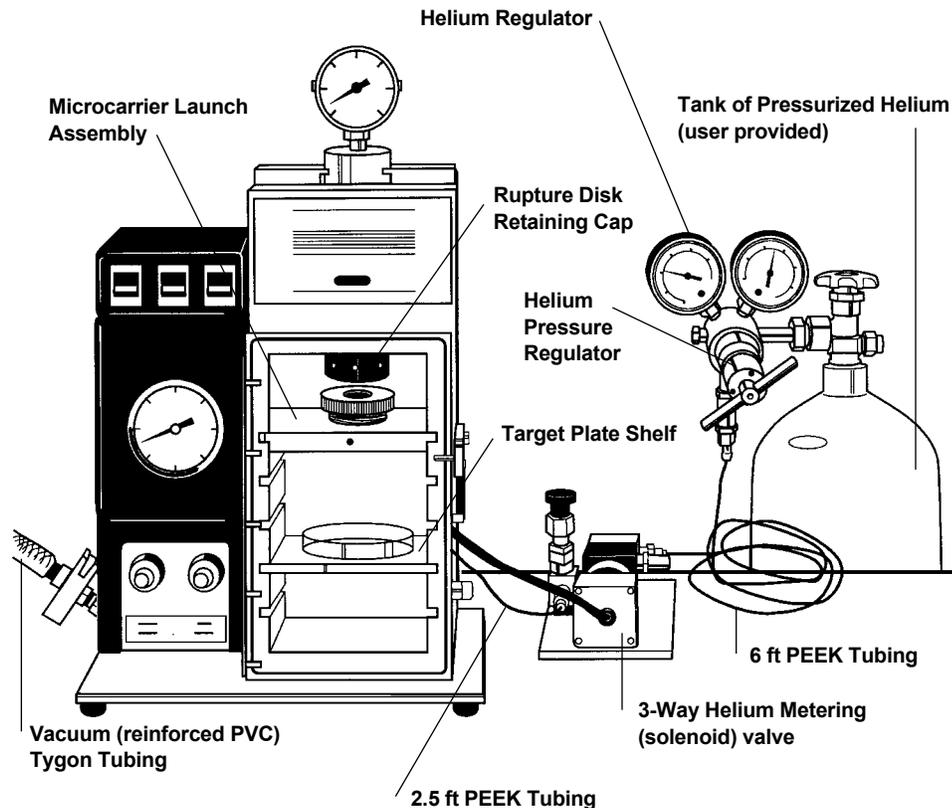


Fig. 1.1. Unit components, front view.

The Biolistic Process

The Biolistic PDS-1000/He system uses high pressure helium, released by a rupture disk, and partial vacuum to propel a macrocarrier sheet loaded with millions of microscopic tungsten or gold microcarriers toward target cells at high velocity. The microcarriers are coated with DNA or other biological material for transformation. The macrocarrier is halted after a short distance by a stopping screen. The DNA-coated microcarriers continue traveling toward the target to penetrate and transform the cells.

The launch velocity of microcarriers for each bombardment is dependent upon the helium pressure (rupture disk selection), the amount of vacuum in the bombardment chamber, the distance from the rupture disk to the macrocarrier (A), the macrocarrier travel distance to the stopping screen (B), and the distance between the stopping screen and target cells (C).

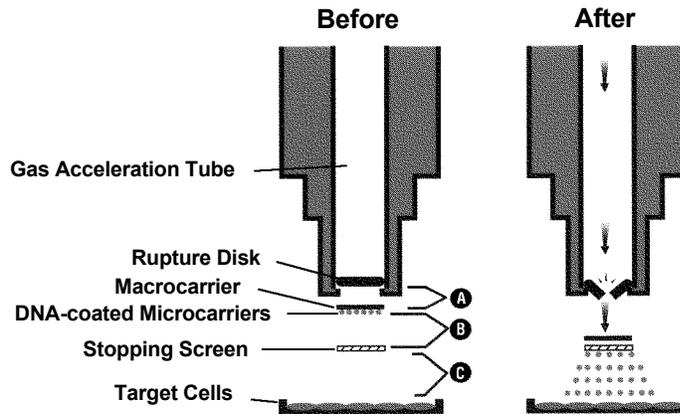


Fig. 1.2. The Biolistic bombardment process.

Design Improvements

The original Biolistic device used a gunpowder explosion to accelerate DNA-coated microcarriers into target cells. The helium technology used in the current PDS-1000/He system has primary advantages of providing cleaner, safer, and more reproducible particle acceleration. This stems from the use of rupture disks that burst at a defined pressure. In addition, helium inflicts less tissue damage. Bulletin 1689 offers a comparative analysis of gunpowder and helium target patterns and transformation efficiencies.

The PDS-1000/He device was updated in March 1995 to improve the quality of key components. We removed parts originally designed to operate with the gunpowder acceleration method. The most notable change was the conversion of the material used for the bombardment chamber from metal to a strong, lightweight plastic. This makes the instrument easier to transport and clean, with no change in bombardment performance (identical internal chamber dimensions). An over-pressure relief valve and a particle filter on the vacuum vent supply were also added.

The actual steps for performing a particle bombardment of a biological sample are unchanged, and the consumables are also the same with the plastic case version. Extensive testing involving the genetic transformation of yeast, plant, and animal cells by both Bio-Rad and independent researchers demonstrated that the gene transfer results obtained with the new plastic chamber design are equivalent to those of the previous metal chamber model. See Appendix 8.2 for a description of parts unique to the metal-chamber design.

1.3 Important Safety Information



Pressurized Helium Safety Information

Caution: Although helium is neither toxic nor flammable, all gases under pressure are potentially dangerous if used improperly. Never use a helium tank with, or attach a tank to, the PDS-1000/He system unless the tank is properly secured. Follow the instructions provided with the helium cylinder from the supplier and those that are applicable for your institution (site safety officer). Bio-Rad has supplied tubing, fittings, a control valve, and a pressure regulator capable of safely handling the high pressure helium gas used in the Biolistic bombardment process. These components have been carefully selected and are the only parts to be used with the PDS-1000/He system.

Power Safety Information

Figure 1.3 shows the serial number certification label which is found at the rear of the Biolistic PDS-1000/He unit. This label provides the manufacturing data about the instrument, its voltage settings, and CDRH standards for electrical safety. This instrument and its accessories conform to the IEC and CDRH standards for electrical safety.

BIO-RAD	Made in U.S.A.
Model No.	Serial No.
<input type="text"/>	<input type="text"/>
Voltage	
<input type="text"/>	

Fig. 1.3. Instrument serial number label on the rear of the instrument.

1.4 Requirements for System Operation

Selecting Site for Operation

Prepare a space 61 cm wide x 46 cm long x 61 cm high (24 inches x 18 inches x 24 inches) preferably in a bio-containment hood or other tissue preparation area, near a standard electrical outlet (110 V/60 Hz in the U.S.). Also, allow for placement of the vacuum pump near the site of operation if house vacuum is not used (see Vacuum Supply).

User Supplied Components

Helium Supply

Only helium gas is to be used with PDS-1000/He system. The low atomic weight of helium permits maximum gas expansion into the bombardment chamber. Thus, sufficient acceleration of the DNA-coated microcarriers is generated for penetration of the target cell membrane.

Obtain a high pressure (2,400 to 2,600 psi) tank of high purity helium for optimization of bombardment conditions for the biological system of choice. This allows use of all of the rupture disks (the highest disk has a 2,200 psi rating). Only grade 5 (99.999%) or grade 4.5 (99.995%) helium is to be used, since helium of a lesser grade contains contaminating material which may obstruct gas flow within the PDS-1000/He system, as well as contaminate the biological sample. Follow all safety instructions provided by helium supplier for helium tank installation.

The helium pressure regulator (supplied) has a CGA 580, female fitting (standard in the United States) for attachment to the user-supplied helium tank. An adaptor to this fitting may be required outside the United States. Contact your local Bio-Rad office for information on the helium pressure regulator adaptor requirements in your location.

A user-supplied, 1 inch adjustable wrench is required for attachment of the regulator to a helium tank having a capacity of 55 cubic feet or greater.

Vacuum Source

The main unit of the PDS-1000/He system must be connected to a vacuum source capable of evacuating the bombardment chamber to a minimum of 5 inches of mercury for operation. This minimum vacuum requirement is part of the instrument safety system (the chamber door must be sealed for helium pressure to be delivered into the main unit). Tubing is supplied to connect the PDS-1000/He system to vacuum source.

For maximum evacuation capacity, we recommend connecting the PDS-1000/He device to an oil-filled, rotary vane vacuum pump, either single or dual stage, with an pumping speed of 90–150 liters/minute (3–5 cubic feet/minute). This pumping rate minimizes the time target cells are exposed to vacuum. Oil for the vacuum pump must be supplied by the user. An exhaust mist eliminator on the vacuum pump is also recommended.

The level of the vacuum required in the bombardment chamber depends on the biological system being targeted for transformation. Higher vacuum reduces drag forces on microcarriers during helium-driven acceleration. During the brief bombardment process (less than 1 minute), typical protocols require a vacuum level within the bombardment chamber between 15–29 inches of mercury. Some cells, tissues, and intact plant cells require a high vacuum (up to 28 inches of mercury) for efficient transformation.

House vacuum may be sufficient for bombardment of certain cell types (mammalian cells), but house vacuum pumping rates can fluctuate and vary greatly in overall evacuation capacity (typically 20 inches of mercury, maximum).

Consumables

The 500 Optimization Kit (catalog number 165-2278) provides the consumables needed for 500 bombardments. It is recommended for users who have yet to determine the optimal conditions for the bombardment of the biological system of interest. The kit contains 0.25 g each of 0.6 μ , 1.0 μ , and 1.6 μ gold microcarriers, 100 each of the nine different rupture disks (ranging from 450 psi to 2,200 psi), 500 macrocarriers, and 500 stopping screens. After optimal conditions are determined, Standard Pressure Kits are available (see Section 7 or the current Bio-Rad catalog for a complete listing).

Additional Laboratory Supplies and Equipment

Vortex mixer is needed for microcarrier preparation.

Common laboratory supplies, such as 95% ethanol, pipettes, etc. are required, as cited in standard protocols.

Additional Items Available from Bio-Rad

The Yeast Optimization Kit (catalog number 170-3100) allows first-time users to become familiar with the Biolistic instrument. It is also helpful for experienced users wishing to periodically standardize their bombardment conditions. This kit provides all of the biological material needed to transform yeast. Yeast provides a system that is quickly and easily assayed. The kit demonstrates the effect of varying the DNA concentration, rupture disk pressure, and the target distance (from the stopping screen), and includes *Saccharomyces cerevisiae* strain

948, YEp352 DNA, CaCl₂, spermidine, culture medium, and plating medium. Enough material for 60 bombardments is provided.

Macrocarrier Holders, set of 5 (catalog number 165-2322), are included with the Biolistic PDS-1000/He system. Additional holders are desirable to facilitate a series of bombardments in one experiment.

The Disk-Vac (catalog number 165-2323) is a small pen-shaped device capable of generating a suction for efficient handling of rupture disks and macrocarriers. Use of the Disc-Vac reduces static generated during manipulation and prevents contamination with glove powder or oil from skin.

Customers outside the continental USA and Canada will require a Voltage Converter (catalog number 165-2259). To use the system, you must locally obtain a cord set which has an IEC/320/CEE 22 connector on one end. This connects to the Voltage Converter and is the type commonly found on computers or televisions. The other end of the power cord will terminate in a plug which will fit the receptacle used in your location. Contact your local Bio-Rad office for more information on this Voltage Converter.

Section 2 Product Description

2.1 Packing List

Check the items received with your PDS-1000/He unit against the list below. If items are missing, contact your local Bio-Rad office.

Instruction Manual

Pressure Regulator for Helium Cylinder (with in-line 0.45 μ filter)

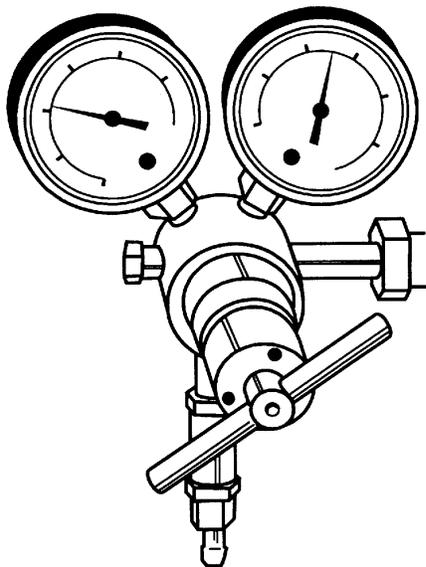


Fig. 2.1. Helium pressure regulator.

6.0 ft PEEK plastic tubing (1/16" OD x .010" ID tubing & fittings), used to connect 3-way helium metering (solenoid) valve to helium pressure regulator

2.5 ft PEEK plastic tubing (1/16" OD x .010" ID tubing and fittings), used to connect 3-way helium metering (solenoid) valve to rear of main unit

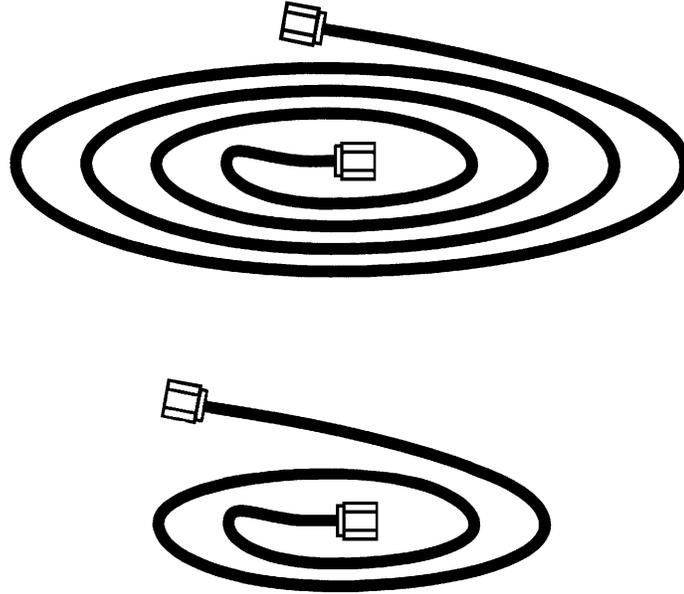


Fig. 2.2. 6 ft and 2.5 ft PEEK tubing.

3-Way Helium Metering (solenoid) Valve, with attached cord for power connection to main unit

Power cord (US and Canada only, 120 V three prong plug)

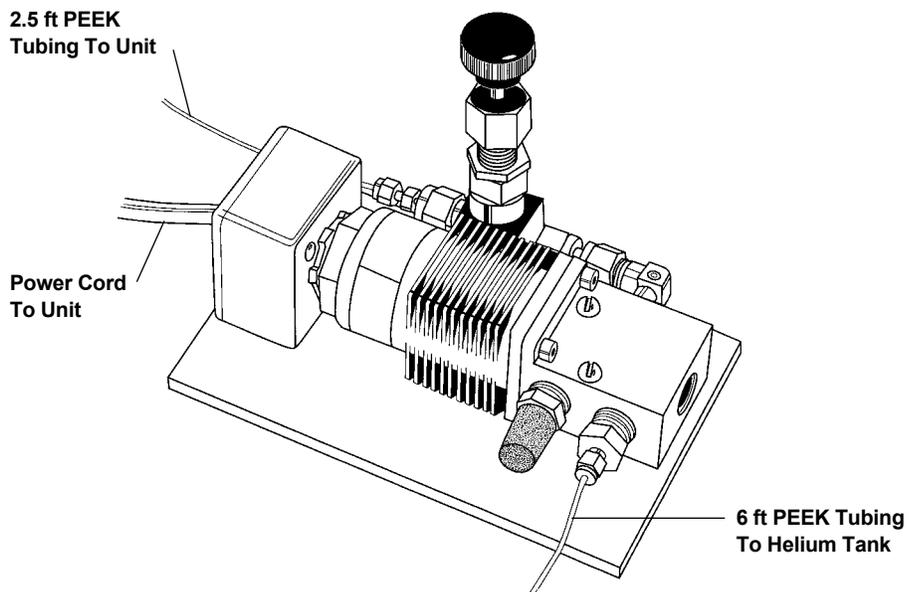


Fig. 2.3. 3-way helium metering (solenoid) valve.

Tools

- 3 Hexagonal gap setting tools (1 each- 1/8", 1/4", and 3/8"), used to set gap between the bottom of the rupture disk retaining cap and the lid of the microcarrier launch assembly
- 1 3/16" hex key (Allen) wrench, used for removal of gas acceleration tube (service only)
- 1 1/8" hex key (Allen) wrench, used for releasing set screw in microcarrier launch assembly shelf
- 2 1/4" x 5/16" open-end wrenches, used for connecting small Swagelock fittings of plastic PEEK tubing to rear of instrument, external solenoid valve, and helium pressure regulator
- 1 torque wrench for rupture disk retaining cap
- 1 seating tool for macrocarriers

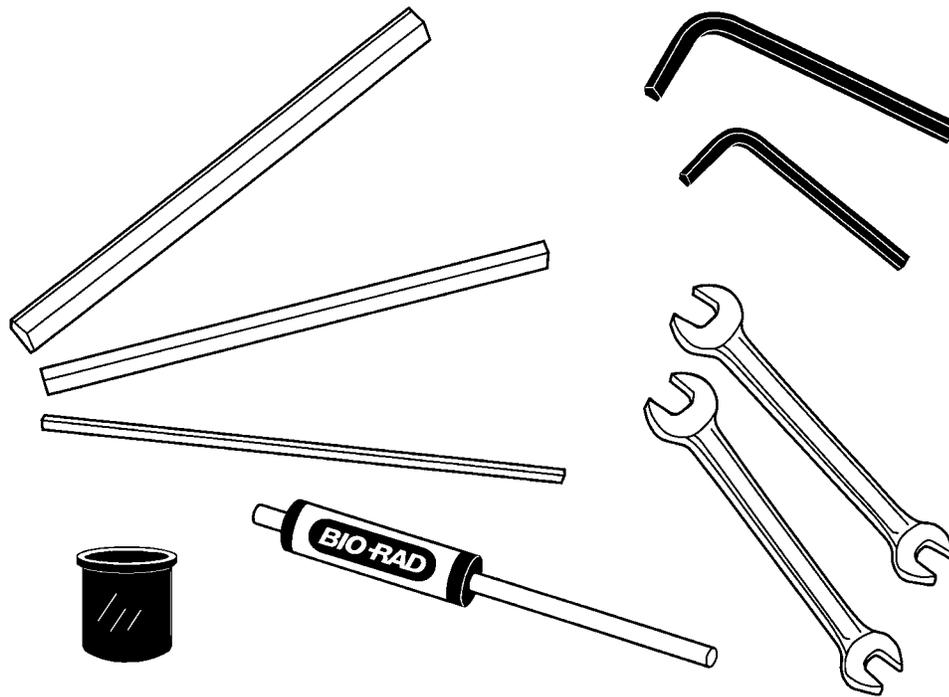


Fig. 2.4. PDS-1000/He tools.

PDS-1000/He Main unit (bombardment chamber with control panel and gauges; shipped fully assembled)

Reinforced PVC vacuum tubing (1/2" ID x 3/4" OD, 5 ft. length & fitting) attached to rear of unit by clamp assembly (centering ring, vacuum hose clamping ring and nozzle adaptor; see Section 3.2 for individual components)

Rupture disk retaining cap (with torque wrench placement holes; Figure 10), attached to gas acceleration tube within bombardment chamber

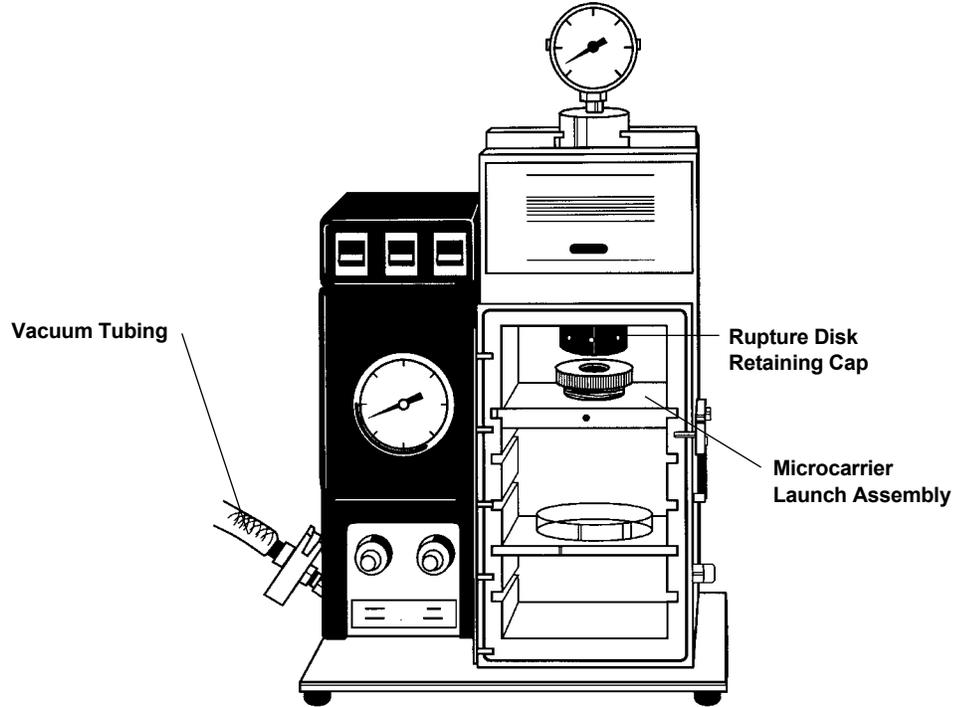


Fig. 2.5. PDS-1000/He main unit.

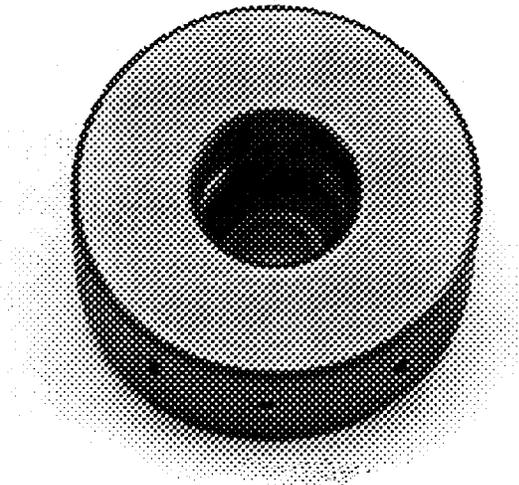


Fig. 2.6. Rupture disk retaining cap.

Microcarrier launch assembly (shipped fully assembled) consists of the following:

- Launch Assembly Shelf with Recessed Set Screw
- Macrocarrier Cover Lid
- Adjustable Nest
- Fixed Nest with Retaining Spring
- Stopping screen Support Ring
- Spacer Rings, 5 mm height, 2
- Macrocarrier Holders, 5, for use within microcarrier launch assembly, after macrocarrier is inserted using macrocarrier insertion tool

Target Plate Shelf

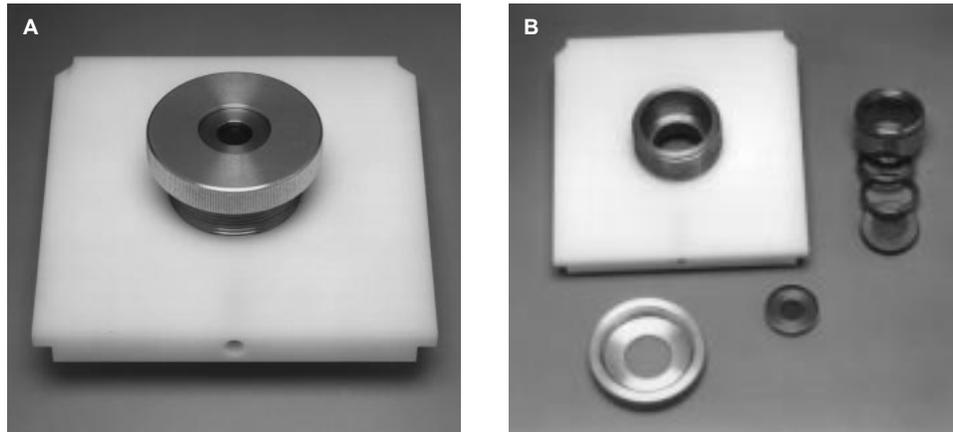


Fig. 2.7. Microcarrier launch assembly (A) and disassembled components (B).

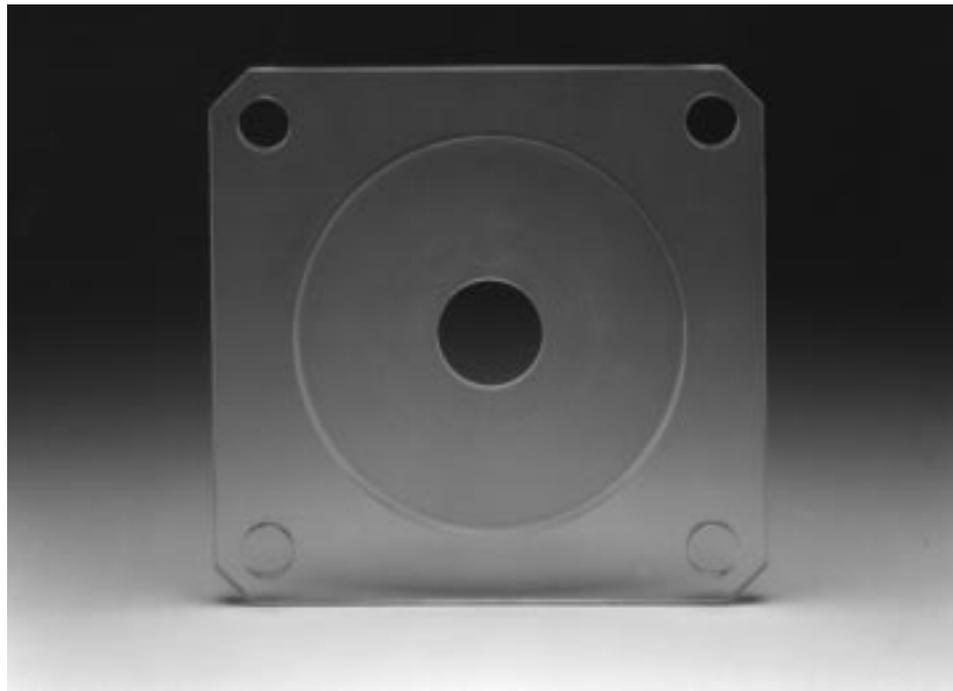


Fig. 2.8. Target plate shelf.

2.2 Identification of Unit Controls and Components

The following is a brief description of the operation controls for the PDS-1000/He unit (Figure 2.9)

Table 2.1. Unit Controls and Components

Controls and Components	Description
Front View- Exterior	
Power Switch, ON/OFF	Controls supply of line electrical power to the instrument.
Fire Switch	<p>Controls flow of helium into Gas Acceleration Tube by activating Solenoid Valve. Illuminated red when enabled, <i>i.e.</i> when safety interlock is satisfied that at least 5 inches Hg. vacuum is present in chamber.</p> <p>Fire Switch must be held ON continuously until Rupture Disk bursts; then release Fire Switch to stop flow of helium.</p> <p>If the Fire Switch is released before the disk ruptures, the helium is vented via a safety vent in the external three-way metering (solenoid) valve.</p>
Vac/Vent/Hold Switch	<p>Controls application of vacuum to bombardment chamber. Vac applies vacuum from line source. Vent releases vacuum using filtered air. Hold maintains vacuum by isolating chamber.</p> <p><i>Bombardments should be performed with this switch in "Hold" position.</i></p>
Vacuum Gauge	Indicates level of vacuum in bombardment chamber, in inches of mercury where zero equals ambient atmospheric pressure.
Vacuum/Vent Rate Control Valves	Regulate rate of application and relief of vacuum in bombardment chamber. Clockwise rotation closes valves.
Helium Pressure Gauge	Indicates helium pressure in Gas Acceleration Tube, in psi. When solenoid valve is activated by Fire Switch, the needle in this oil-filled gauge rotates clockwise until rupture disk bursts. <i>Watch this gauge carefully during a bombardment and note the actual rupture pressure.</i>
Gas Acceleration Tube	Helium accumulates within this tube when it is sealed by rupture disk at chamber-end of Tube. Helium PEEK tubing connects at top of Gas Acceleration Tube outside, rear of chamber.
Bombardment Chamber	Holds Rupture Disk, Microcarrier Launch Assembly, and biological target under vacuum during a bombardment.

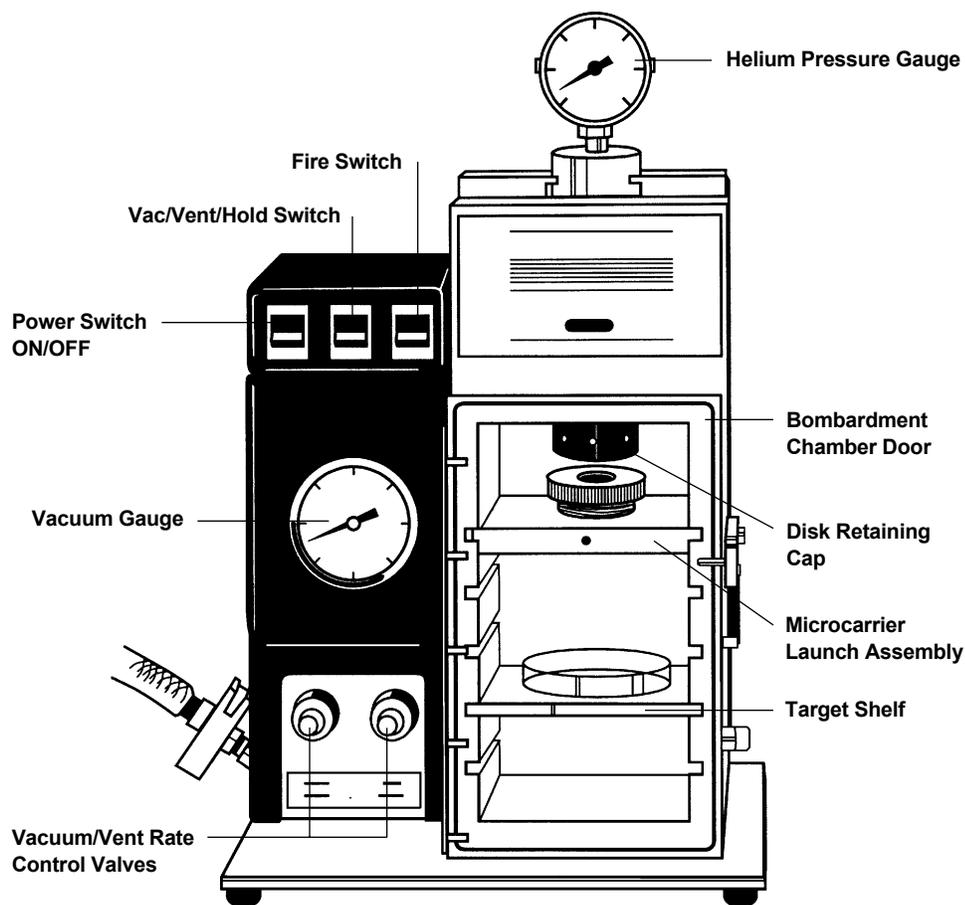


Fig. 2.9. Front view of PDS-1000/He unit.

Table 2.2. Front View-Interior of Bombardment Chamber

Bombardment Chamber Door (with Brace)	Closes chamber with a solid piece of polycarbonate plastic. Note the single, large o-ring which seals vacuum in the chamber, and the self-positioning brace which eliminates flex of chamber walls during bombardment cycle.
Rupture Disk Retaining Cap	Seals Rupture Disk against chamber end of Gas Acceleration Tube. This must be tightened securely. The Torque Wrench is used in the holes in the cap which are visible in the photo.
Microcarrier Launch Assembly	Holds the DNA/microcarrier preparation on a Macrocarrier sheet over the Stopping Screen in the path of the helium shock wave.
Target Shelf	Holds the biological target in a Petri plate in the path of the accelerated DNA/microcarrier preparation. Particle flight distance is determined by positioning the shelf at one of four levels using slots in the chamber walls.

Table 2.3. Rear Connections

Refer to Figure 2.10 for a rear perspective view of the unit.

Helium Connection to Gas Tube	Connects top of Gas Acceleration Tube to plastic tubing from Solenoid Valve, supplying high pressure helium.
Over-Pressure Relief Valve	Opens at 0.5 psi chamber pressure to relieve accumulation of gas. A new safety feature. Automatically resets after activation.
Vacuum Line, Chamber to Controls	Supplies vacuum to chamber from control valves.
Helium Metering (Solenoid) Valve Electrical Connection	Supplies electric power to 3-way Solenoid Valve in helium line. Plug the three-pin connector from the Solenoid Valve into this receptacle.
Line Cord Electrical Connection	Supplies electric power to Biolistic unit from house line. Plug connector from Line Cord into this receptacle.
Vacuum Line, Connection to Source	Connects unit to house vacuum supply or vacuum pump via fittings included. Feeds into vacuum flow controls.

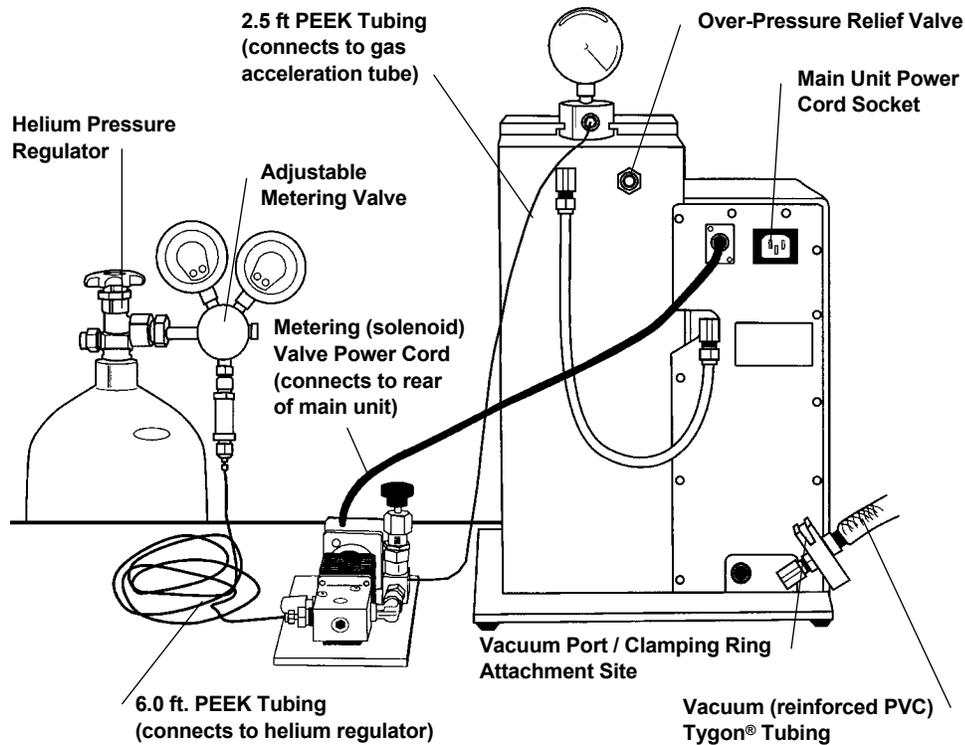


Fig. 2.10. Rear view of main unit, component connection points.

Section 3 Installation

3.1 Connecting the PDS-1000/He System to a Helium Source

Refer to Section 2.2, Identification of Unit Controls and Components, prior to system installation.

Helium Pressure Regulator Installation

Connecting the helium pressure regulator to a tank of pressurized helium.

Components needed:

- Pressure regulator for helium cylinder (with 0.45 micron in-line filter), provided with unit (Figure 3.1).
- Cylinder of grade 4.5 to 5.0 helium (minimum 99.995% pure); maximum pressure of 2,600 psi, user supplied.
- 1 1/8" open-end wrench or a 10" or 12" adjustable wrench, user supplied.

Note: The regulator is intended for use only with helium gas under a maximum of 2,600 psi of pressure. The outlet on pressurized helium cylinders used in the United States (maximum 2,600 psi) is compatible with the fitting supplied on the pressure regulator with the PDS-1000/He unit (CGA 580, female fitting). Outside of the US, contact your local Bio-Rad office for information regarding the proper cylinder/regulator fitting in your area.

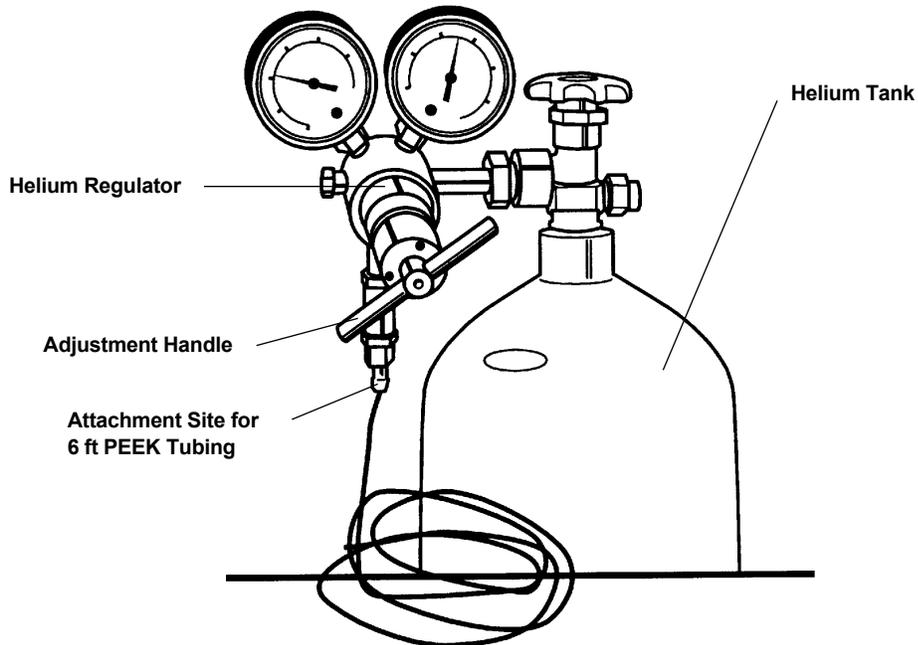


Fig. 3.1. Helium regulator/tank attachment.

1. Secure the cylinder to a wall, post, or other anchored fixture so it will not tip or fall during use.
2. Inspect the cylinder valve for dirt, dust, oil, grease, or damaged threads. Remove dust and dirt with a clean cloth. Do not attach the regulator if you determine that the valve port is damaged or cannot be cleaned. Inform your gas supplier of this condition and request a replacement cylinder.
3. Clear the valve port of any foreign matter by standing to the side of the cylinder and quickly opening and closing the cylinder valve.
4. Attach the regulator to the cylinder valve and tighten securely with a 1 1/8" open-end wrench or a 10" or 12" adjustable wrench.

Three-Way Metering Valve (Solenoid) Installation

Connecting the helium pressure regulator to the 3-way helium metering (solenoid) valve

Components needed:

- Two 5/16" open end wrenches (Figure 2.4)
- PEEK tubing, 6 feet (Figure 2.2)
- Solenoid valve (Figure 2.3)

Note: The Swagelock fittings used on the flexible tubing are easily damaged by over tightening. Turn finger-tight, then use wrench for 1/6 additional turn (60°).

1. Connect one end of the 6 foot tubing assembly to helium pressure regulator (Figure 3.1). Attach the tubing to the in-line, 0.45 micron filter at the end of the regulator. Use two 5/16" open end wrenches to tighten the connection.
2. Connect the other end of the 6 foot tubing assembly to the inlet port of the 3-way helium (solenoid) valve assembly. This is the port nearest the metal exhaust filter on the 3-way metering valve. Use two 5/16" open end wrench to tighten the connection (Figures 2.3 and 2.10).

Connecting the 3-way helium metering (solenoid) valve to the PDS-1000/He main unit

Components needed:

- Two 5/16" open end wrenches (Figure 2.4)
- PEEK tubing, 2.5 feet (Figure 2.2)
- Solenoid valve (Figure 2.3)

Note: The Swagelock fittings used on the flexible tubing are easily damaged by over tightening. Turn finger-tight, then use the wrench for 1/6 additional turn (60°).

1. Connect one end of the 2.5 foot PEEK tubing assembly to the outlet port of the 3-way helium metering (solenoid) valve assembly (Figures 2.3 and 2.10). This is the port nearest the adjustable metering valve (black knob). Use two 5/16" open end wrenches to tighten the connection. Do not alter the setting of the metering valve at this time.
2. Connect one end of the 2.5 foot long PEEK tubing assembly to the PDS-1000/He unit. This port is located on the gas acceleration tube, below the helium pressure gauge (top, rear of unit). Use a 5/16" open end wrench to tighten the connection (Figure 2.10).
3. Plug the power cord of the 3-way helium metering (solenoid) valve into the rear of the PDS-1000/He unit (Figure 2.10).

3.2 Connecting the PDS-1000/He Unit to a Vacuum Source

The main unit is shipped with the vacuum hose assembly attached to the port on the rear of the unit, via the clamping assembly. Only the connection of the free end of the tubing to a vacuum source is required for connecting the vacuum source to the system.

Choice of Vacuum Source

A vacuum system with a 3 cfm (cubic feet per minute) or 100 l/min pumping capacity is recommended for use with the PDS-1000/He unit. This minimizes the total time the target cells/tissue are exposed to reduced atmospheric pressure when inside the bombardment chamber. It is also recommended that the pump have an exhaust mist eliminator.

House or shared source vacuum systems commonly have variable evacuation capacities, from 20 inches to 23 inches Hg. When multiple users access a common vacuum system, the actual vacuum pressure will vary, often decreasing to less than 20 inches Hg. per outlet. This is an inadequate vacuum level for yeast and plant cell/tissue bombardment.

Most rotary vacuum pumps will achieve up to 26–28 inches of mercury (Hg) vacuum. However, the length of time required to attain the desired vacuum depends on the pumping capacity of the vacuum source. The higher the pumping capacity, the less time is required to achieve the desired vacuum. Variation in evacuation time in the bombardment chamber has an unknown effect on target cells and tissue transformation efficiency.

Table 3.2. Pumping rates of various vacuum systems with the PDS-1000/He system

Vacuum Source Pumping Capacity	Vacuum Level (inches of Hg)	Evacuation Time†
House/Common Source	20	10–15 seconds
	29	Typically not attainable
Pump - 1 cfm or 28 l/min	20	12 seconds
	29	45 seconds
Pump - 3 cfm or 100 l/min	20	7 seconds
	29	25 seconds

† Vacuum Flow Rate Valve fully open.

Replacement and Re-assembly of connecting components

See Figure 3.2 for a diagram outlining component assembly. The clamp consists of an O-ring and a vacuum hose centering ring with male nozzle fitting screwed to it.

1. Place the clamping assembly at one end of the vacuum tubing. The vacuum tubing is PVC-reinforced Tygon tubing (1/2" ID x 3/4" OD; 5 feet length).
2. The vacuum tubing is shipped with the clamping ring assembly (Figure 3.2) attached to the PDS-1000/He unit at the brass vacuum port located on the rear of the unit.
3. Attach the free end of the vacuum hose assembly to the vacuum source.

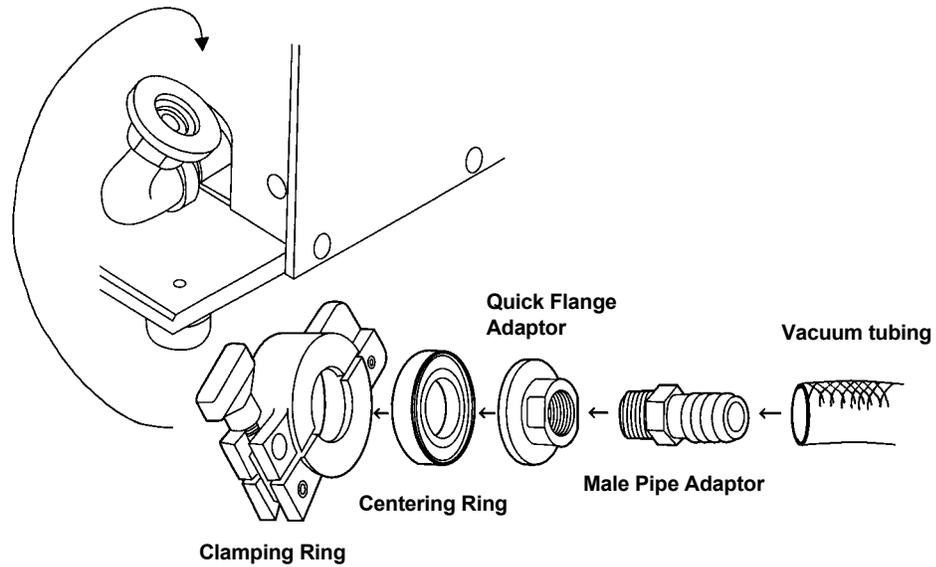


Fig. 3.2. Components used to connect vacuum source to main unit.

3.3 Power Cord / Voltage Regulator

The power cord must be plugged directly into a 100 V-120 V outlet. Plug other end into unit. For use with a 220 V or 240 V line voltage (users outside the US or Canada), connect a voltage converter to the power cord prior to use (catalog number 165-2259, see Figure 3.3). You must supply a cord set that has an IEC-320/CEE-22 connector on the end that attaches to the voltage converter; the other end of the cord set will terminate with a plug that is compatible with the electrical outlets in your area.



Fig. 3.3. Voltage converter.

Section 4 Operation of the PDS-1000/He Instrument

4.1 Preparation of System Components Prior to Bombardment

Instrument Preparation

1. Verify that helium tank has 200 psi in excess of desired rupture disk pressure for bombardment.
2. Set gap distance between rupture disk retaining cap and microcarrier launch assembly.

When the rupture disk retaining cap is in place, insert the fully assembled microcarrier launch assembly (with cover lid) inside the bombardment chamber on the highest possible bombardment chamber wall slot (Figure 4.1). The set screw in the white plastic shelf should face outward. Release the set screw on the front of the microcarrier launch assembly with the smaller of the two hex key wrenches provided (Figure 2.4).

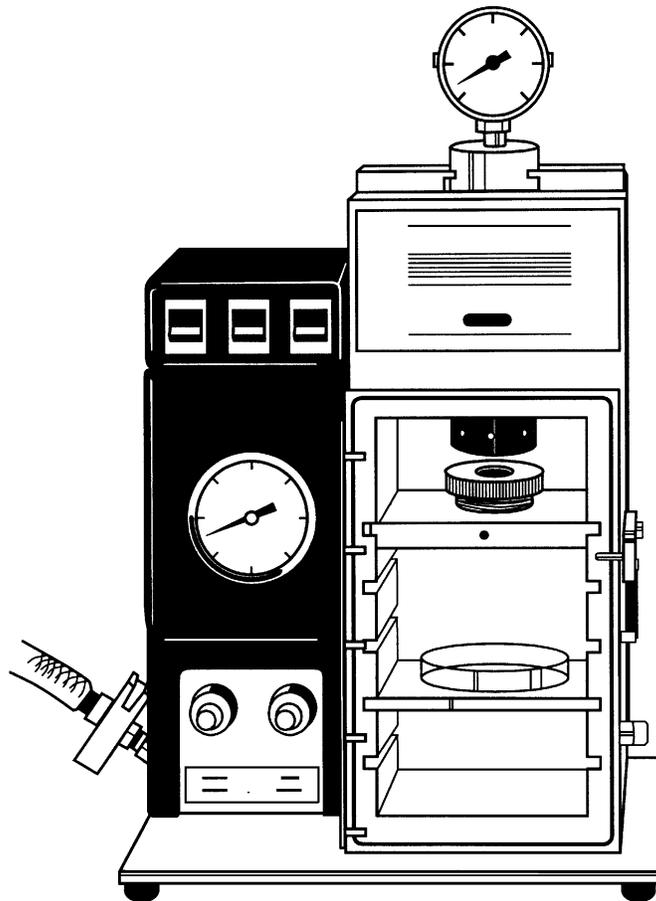


Fig. 4.1. Microcarrier launch assembly and target shelf inside bombardment chamber.

Three hexagonal gap adjustment tools of 1/8", 1/4", and 3/8" have been provided to reproducibly set the gap distance (Figure 2.4). A 1/4" distance between the rupture disk retaining cap and the macrocarrier cover lid is recommended for initial optimization bombardments.

While holding the hexagonal gap adjustment tool against the bottom of the rupture disk retaining cap, turn the adjustable nest until the macrocarrier cover lid touches the gap adjustment tool. The position of the adjustable nest is fixed by tightening the set screw in the white plastic shelf with the hex key wrench.

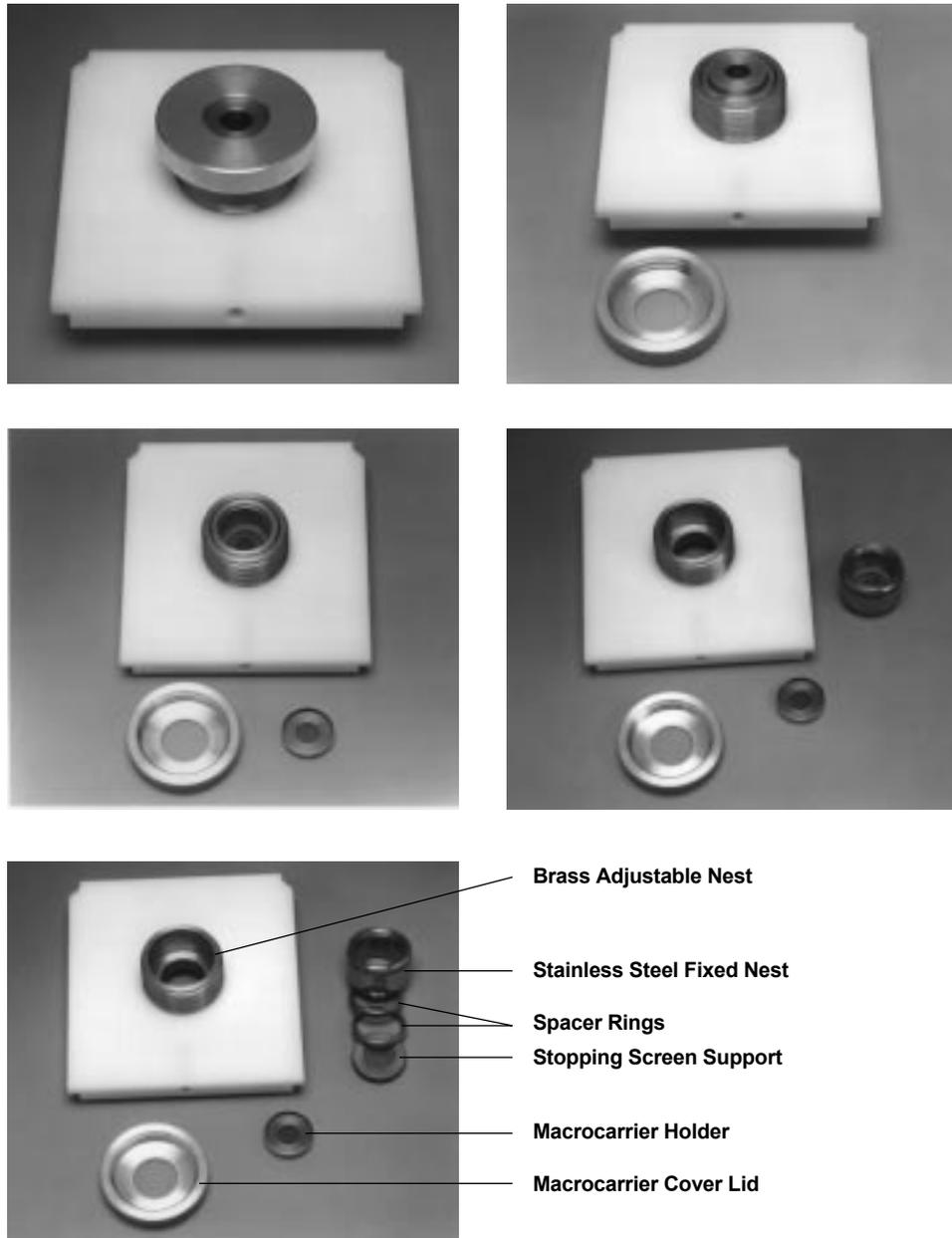


Fig. 4.2. Microcarrier Launch Assembly adjustment.

3. Prepare the Rupture Disk Retaining Cap. After setting the gap between cap and microcarrier launch assembly, wrap the Rupture Disk Retaining Cap in aluminum foil and sterilize by autoclaving.

4. Prepare the Microcarrier Launch Assembly.

The effect of the gas shock wave on the microcarrier velocities is determined in part by the gap between the rupture disk and the macrocarrier.

The macrocarrier flight distance can be adjusted by varying the positions of the stopping screen support and the spacer rings inside the fixed nest (Figure 4.2). The Stopping Screen Support is placed in the middle position at the factory. This is the recommended position when initially optimizing bombardment parameters.

To make this adjustment, remove the microcarrier launch assembly from the PDS-1000/He unit. Unscrew and remove the macrocarrier cover lid. Disassemble the components of the microcarrier launch assembly by placing the macrocarrier insertion tool into the bottom of the assembly and pushing up. This releases the stainless steel fixed nest from the brass adjustable nest. The two spacer rings and the stopping screen support will fall out from within the fixed nest.

To change the factory-set position to the position for the minimum macrocarrier travel distance (6 mm), invert the fixed nest and insert the stopping screen support inside the fixed nest so that the conical side of the stopping screen support faces down in the final orientation; then insert the two spacer rings (5 mm thickness, each). With the fixed nest still inverted, place the macrocarrier launch assembly over the fixed nest and seat the fixed nest within the brass adjustable nest.

If greater macrocarrier travel is desired, rearrange the spacer rings and stopping screen support accordingly. The macrocarrier travel distance can be increased in two 5 mm steps, to a maximum of 16 mm.

Sterilize the microcarrier launch assembly by wrapping in aluminum foil and autoclaving. Alternatively, this assembly can be sterilized by wiping with 70% ethanol, followed by drying in a sterile environment.

5. Target Shelf

Sterilize the target shelf by wiping with 70% ethanol, followed by drying in a sterile environment just prior to use. This part may not be autoclaved.

Consumable Preparation

Several consumables are available for use with the PDS-1000/He system (Figure 4.3).

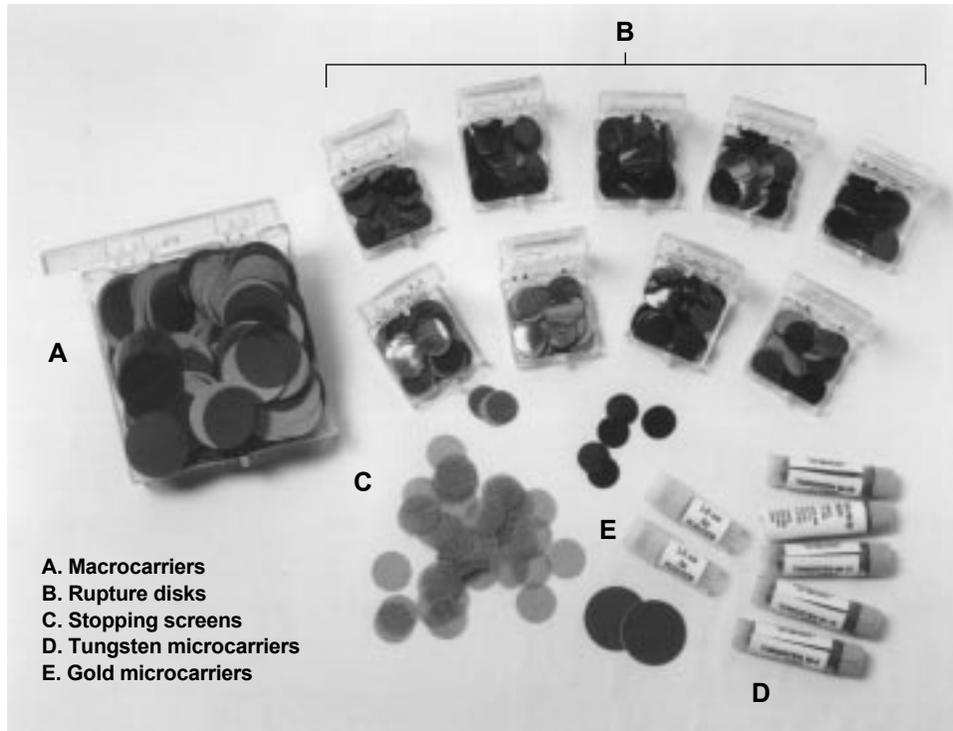


Fig. 4.3. Consumables for the PDS-1000/He instrument. The 0.6 micron gold is not pictured.

1. Macrocarriers

Pre-assemble and pre-sterilize the macrocarrier set in a macrocarrier holder prior to performing sample cell/tissue bombardments. The Disk-Vac (catalog number 165-2323) provides ease in handling rupture disks and macrocarriers.

Macrocarriers are shipped in quantities of 500/box, with paper linings between disks. Transfer selected macrocarriers to individual Petri dishes for easier handling. Remove the paper lining from between the macrocarriers. Place the macrocarrier into the macrocarrier holder using the seating tool (Figure 4.4). The edge of the macrocarrier should be securely inserted under the lip of the macrocarrier holder. The macrocarrier holders, with macrocarriers already in place, should be sterilized by autoclaving.

2. Rupture disks

Rupture disks (Figure 4.3) are packaged and shipped in quantities of 100/box. Transfer selected rupture disks to individual Petri dishes for easier handling. Sterilize rupture disks by briefly dipping them in 70% isopropanol just prior to insertion in the Retaining Cap. Do not soak for more than a few seconds. Extensive soaking may delaminate the disks, resulting in premature rupture. All disks, with the exception of those rated at 450, 650, and 1,100 psi, are laminated. Autoclaving is not recommended because of potential delamination.

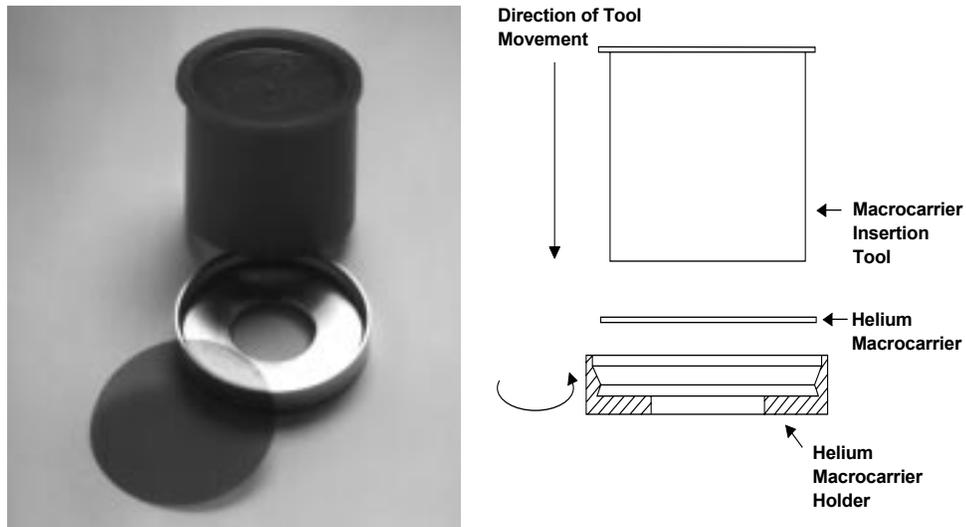


Fig. 4.4A and 4.4B. Insertion of macrocarrier into macrocarrier holder with plastic insertion tool.

3. Stopping screens

Transfer selected stopping screens (Figure 4.3) to individual Petri dishes for easier handling. Sterilization by autoclaving is recommended. Alternatively, these parts can be sterilized by soaking in 70% ethanol, followed by drying in a sterile environment.

4. Microcarriers

The following procedure prepares tungsten or gold microcarriers for 120 bombardments using 500 μg of the microcarrier per bombardment, based on the method of Sanford, *et al.* [*Methods in Enzymology*, **217**, 482-509 (1993)].

Weigh out 30 mg of microparticles into a 1.5 ml microfuge tube.

Add 1 ml of 70% ethanol (v/v).

Vortex vigorously for 3–5 minutes (a platform vortexer is useful).

Allow the particles to soak in 70% ethanol for 15 minutes.

Pellet the microparticles by spinning for 5 seconds in a microfuge.

Remove and discard the supernatant.

Repeat the following wash steps three times:

- Add 1 ml of sterile water.
- Vortex vigorously for 1 minute.
- Allow the particles to settle for 1 minute.
- Pellet the microparticles by briefly spinning in a microfuge.
- Remove the liquid and discard.

After the third wash, add 500 μl sterile 50% glycerol to bring the microparticle concentration to 60 mg/ml (assume no loss during preparation).

The microparticles can be stored at room temperature for up to two weeks. Tungsten aliquots should be stored at $-20\text{ }^{\circ}\text{C}$ to prevent oxidation. Gold aliquots can be stored at $4\text{ }^{\circ}\text{C}$ or room temperature.

Store dry tungsten and gold microcarriers in a dry, non-oxidizing environment to minimize agglomeration.

Coating Washed Microcarriers with DNA

The following procedure is sufficient for six bombardments; if fewer bombardments are needed, adjust the quantities accordingly.

Vortex the microcarriers prepared in 50% glycerol (30 mg/ml) for 5 minutes on a platform vortexer to resuspend and disrupt agglomerated particles.

When removing aliquots of microcarriers, it is important to continuously vortex the tube containing the microcarriers to maximize uniform sampling. When pipetting aliquots, hold the microcentrifuge tube firmly at the top while continually vortexing the base of the tube.

Remove 50 μ l (3 mg) of microcarriers to a 1.5 ml microcentrifuge tube.

Continuous agitation of the microcarriers is needed for uniform DNA precipitation onto microcarriers. For added convenience and/or multiple samples, use a platform attachment on your vortex mixer for holding microcentrifuge tubes.

While vortexing vigorously, add in order:

- 5 μ l DNA (1 μ g/ μ l)
- 50 μ l 2.5 M CaCl_2
- 20 μ l 0.1 M spermidine (free base, tissue culture grade)

Continue vortexing for 2–3 minutes.

Allow the microcarriers to settle for 1 minute.

Pellet microcarriers by spinning for 2 seconds in a microfuge.

Remove the liquid and discard.

Add 140 μ l of 70% ethanol (HPLC or spectrophotometric grade).

Remove the liquid and discard.

Add 140 μ l of 100% ethanol.

Remove the liquid and discard.

Add 48 μ l of 100% ethanol.

Gently resuspend the pellet by tapping the side of the tube several times, and then by vortexing at low speed for 2–3 seconds.

4.2 Performing a Bombardment

Quick Guide (This summary is repeated in Section 8.4 as a tear-out sheet.)

Before the Bombardment

1. Select/adjust bombardment parameters for gap distance between rupture disk retaining cap and microcarrier launch assembly. Placement of stopping screen support in proper position inside fixed nest of microcarrier launch assembly
2. Check helium supply (200 psi in excess of desired rupture pressure).
3. Clean/sterilize:
Equipment: rupture disk retaining cap, microcarrier launch assembly
Consumables: macrocarriers/macrocarrier holders
4. Wash microcarriers and resuspend in 50% glycerol
5. Coat microcarriers with DNA and load onto sterile macrocarrier/macrocarrier holder the day of experiment

Firing the Device

1. Plug in power cord from main unit to electrical outlet.
2. Power ON.
3. Sterilize chamber walls with 70% ethanol.
4. Load sterile rupture disk into sterile retaining cap.
5. Secure retaining cap to end of gas acceleration tube (inside, top of bombardment chamber) and tighten with torque wrench.
6. Load macrocarrier and stopping screen into microcarrier launch assembly.
7. Place microcarrier launch assembly and target cells in chamber and close door.
8. Evacuate chamber, hold vacuum at desired level (minimum 5 inches of mercury).
9. Bombard sample: Fire button continuously depressed until rupture disk bursts and helium pressure gauge drops to zero.
10. Release Fire button.

After the Bombardment

1. Release vacuum from chamber.
2. Target cells removed from chamber.
3. Unload macrocarrier and stopping screen from microcarrier launch assembly.
4. Unload spent rupture disk.
5. Remove helium pressure from the system (after all experiments completed for the day).

Detailed Operation Procedure

Note: We recommend that the first bombardment each day be a “dry run” with no target cells or microcarriers to ensure that the system is set up properly and the gas pathway is filled with helium, not air.

1. Power On

With the unit plugged in to the appropriate electrical outlet or voltage converter, turn on the unit by pressing the ON switch. This is the left-most red control panel switch (Figure 2.9).

2. Helium Pressure

Confirm that the helium tank pressure regulator is set to 200 psi over the selected rupture disk burst pressure (*e.g.*, set the regulator to 1,100 psi when working with a 900 psi rupture disk).

3. Pressurizing the System with Helium

Make certain that the helium pressure regulator is properly installed on the helium tank (see Set-up).

Close the helium pressure regulator by turning the regulator adjustment screw counter-clockwise until the adjusting spring pressure is released.

Release helium into the pressure regulator by carefully and slowly opening the cylinder valve on the helium tank. The cylinder pressure in the tank is indicated on the high pressure gauge (the gauge closest to cylinder).

Set the desired helium delivery pressure for the rupture disk you are using (measured on gauge on the outside, farthest from cylinder) by turning the regulator adjustment handle clockwise. The pressure should be set to 200 psi above the rupture disk rating.

4. Coating microcarriers with DNA

The day of the scheduled bombardment, coat the microcarriers with DNA. To obtain the best results, use the DNA-coated microcarriers as soon as possible.

5. Loading DNA-coated microcarriers onto a macrocarrier/macrocarrier holder

Each macrocarrier is placed inside a macrocarrier holder and sterilized, as described above. Prior to the application of the DNA-coated microcarriers onto a macrocarrier, prepare a small desiccating chamber for each macrocarrier/macrocarrier holder and place away from excessive vibration.

A small desiccating chamber consists of a sterile 60 mm tissue culture Petri dish (with lid) containing CaCl_2 as desiccant in the base of the dish (Figures 4.5 and 4.6). The desiccant is covered with a small piece of filter paper to provide a stable platform for the macrocarrier/macrocarrier holder. The sterile macrocarrier/macrocarrier holder is placed atop the filter paper, with the macrocarrier facing up and the stainless steel holder touching the filter paper.

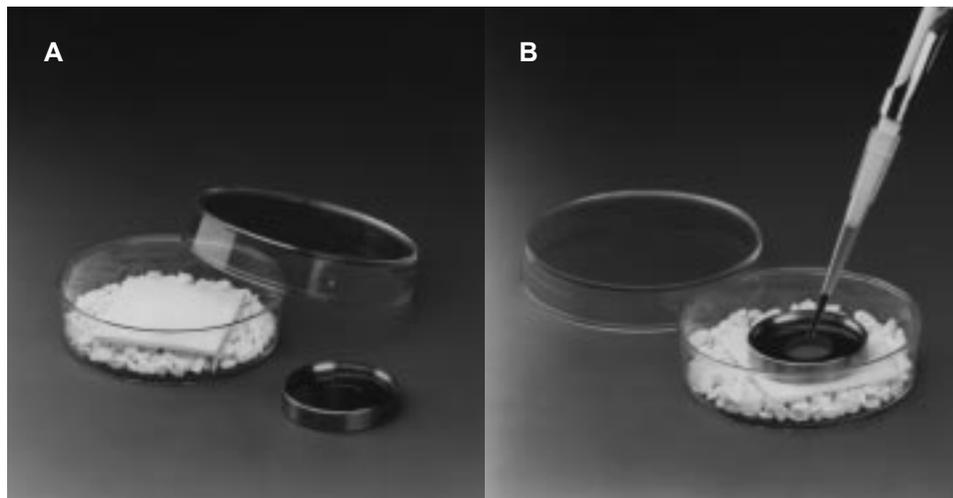


Fig. 4.5A and 4.5B. Loading DNA-coated microcarriers onto a macrocarrier/macrocarrier holder, positioned in small desiccating chamber.

This environment permits rapid access to each macrocarrier and provides a low humidity environment for the ethanol to uniformly evaporate from the microcarriers. This low humidity, along with a minimum of vibration during evaporation, minimizes microcarrier agglomeration.

For each macrocarrier, remove 6 μl aliquots (approximately 500 μg) of microcarriers and spread evenly over the central 1 cm of the macrocarrier using a pipette tip. Pipet from a continuously vortexed tube and rapidly apply suspended microcarriers to the macrocarrier, as microcarriers quickly settle out from the ethanol solution in the tube or even in the pipette tip.

Immediately cover the culture dish after application of the microcarrier suspension to the macrocarrier. The ethanol should evaporate within 10 minutes to leave the DNA-coated microcarriers adhering to the macrocarrier. The loaded macrocarriers should be used within 2 hours.

6. Cleaning chamber walls

Clean the chamber of the PDS-1000/He as desired with 70% ethanol. Allow time for drying. Do not autoclave or flame sterilize the PDS-1000/He unit.

7. Loading the rupture disk

Unscrew the rupture disk retaining cap from the gas acceleration tube from within the bombardment chamber (Figure 4.6) or unwrap cap from sterile wrapping.

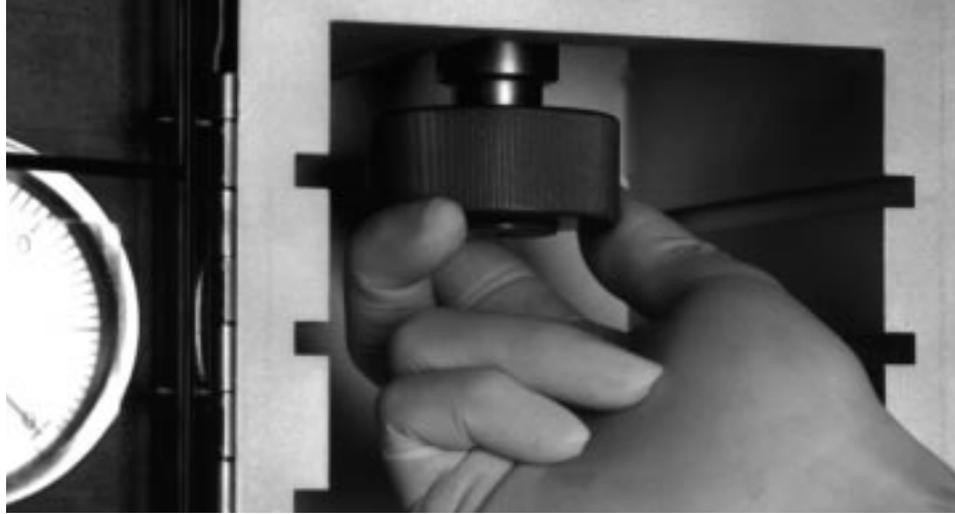


Fig. 4.6. Removal/mounting of rupture disk retaining cap onto end of gas acceleration tube inside bombardment chamber.

Select rupture disk of desired burst pressure. Handle all rupture disks with sterile forceps or Disk-Vac (catalog number 165-2323). Grease, fingerprints, or even powder from plastic gloves on the rupture disk may prevent a tight seal from forming within the retaining cap.



Fig. 4.7. Rupture disk insertion into recess of retaining cap.

Immediately before placing a rupture disk in the retaining cap, briefly wet the rupture disk in isopropanol (do not soak the disk for an extended period of time or the disk may delaminate). Loading the rupture disk while wet forms a liquid gasket when the cap is screwed onto the gas acceleration tube, and thereby reduces the failure rate of the rupture disk. Place the rupture disk in the recess of the rupture disk retaining cap (Figure 4.7).

Screw the rupture disk retaining cap onto the gas acceleration tube using a left-to-right motion. Never tighten the rupture disk retaining cap without a rupture disk in place or scratching and deformation of the two metal surfaces will occur and cause helium to leak when a rupture disk is pressurized.

The retaining cap is tightened to a torque of 60 inch/pounds with the retaining cap torque wrench. To use the torque wrench, insert the short end of the metal rod into an accessible hole in the retaining cap. Push the long end of the rod to the right only until it touches the inner surface of the black tube (Figure 4.8). If the retaining cap is not tightened sufficiently, the rupture disk may slip out of place as the gas acceleration tube fills with helium. Test fire once to fill gas tubing with helium.



Fig. 4.8. Proper torque applied to retaining cap with torque wrench.

8. Microcarrier Launch Assembly

Unscrew the macrocarrier cover lid from the assembly. Place a sterile stopping screen on the stopping screen support (Figure 4.9). **Note:** Never operate the PDS-1000/He instrument without a stopping screen in place. The target sample will be destroyed from the uninterrupted acceleration of the macrocarrier by the helium shock wave.

Install the macrocarrier/macrocarrier holder on the top rim of the fixed nest (Figure 4.10). The dried microcarriers should be facing down, toward the stopping screen. Replace the macrocarrier cover lid on the assembly and turn clockwise until snug. Do not over-tighten.

Place the microcarrier launch assembly in the top slot inside the bombardment chamber (Figures 4.11 and 5.1).

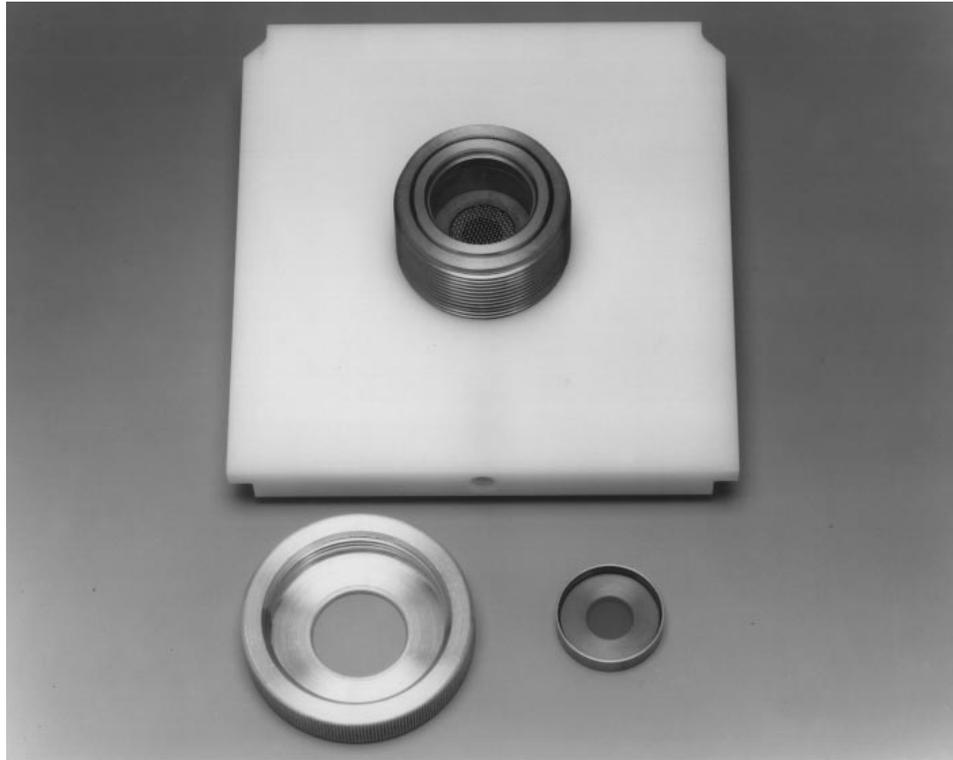


Fig. 4.9. Placement of stopping screen inside fixed nest with macrocarrier and cover lid removed.



Fig. 4.10. Removal / replacement of macrocarrier cover lid with assembled fixed nest. Macrocarrier holder (with macrocarrier properly inserted) is inverted and placed atop fixed nest.

9. **Target cells/tissue placement in chamber**

Place the Target Shelf at the desired level inside the bombardment chamber. Place the sample (usually contained within a Petri dish) on the Target Shelf. Close and latch the sample chamber door.

10. **Chamber evacuation/hold**

Turn on the vacuum source. Set the vacuum switch on the PDS-1000/He (middle red control switch, Figure 4.11) to the **VAC** position. Evacuate the sample chamber to the desired level, at least 5 inches of mercury. The rightmost red control switch (the **FIRE** switch) will be illuminated when the minimum vacuum is achieved.

When the desired vacuum level is reached, hold the chamber vacuum at that level by quickly pressing the vacuum control switch through the middle **VENT** position to the bottom **HOLD** position.



Fig. 4.11. Sample bombardment: power switch ON, vacuum switch on HOLD position, and user continuously pressing FIRE button.

11. **Bombard the sample**

With the vacuum level in the bombardment chamber stabilized, press and hold the **FIRE** switch to allow helium pressure to build inside the gas acceleration tube that is sealed by a selected rupture disk (Figure 4.11).

Estimate rupture disk burst pressure by observing the helium pressure gauge at the top of the acceleration tube. A small pop will be heard when the rupture disk bursts. The rupture disk should burst within 10% of the indicated rupture pressure and within 11–13 seconds.

Release the **FIRE** switch immediately after the disk ruptures to avoid wasting helium. Releasing the **FIRE** switch prior to disk rupture will vent the gas acceleration tube via the 3-way helium metering (solenoid) valve.

Note: Variation in the burst pressure indicated on the helium pressure gauge (on the top of the unit) from the rated rupture disk pressure may be observed if the gas acceleration tube fill rate is improperly set. See Section 5.4 for solenoid valve adjustment procedure.

12. Release vacuum from chamber

Release the vacuum in the sample chamber by setting the **VACUUM** switch to the middle **VENT** position.

13. Target cells removal from chamber

After vacuum is released, the vacuum gauge should read 0 inches of mercury (Hg) of vacuum. Open the sample chamber door. Remove the sample and treat as appropriate.

14. Macrocarrier and stopping screen removal from microcarrier launch assembly

Remove the microcarrier launch assembly. Unscrew the lid and remove the macrocarrier holder. Discard the used macrocarrier and stopping screen (Figure 4.12).

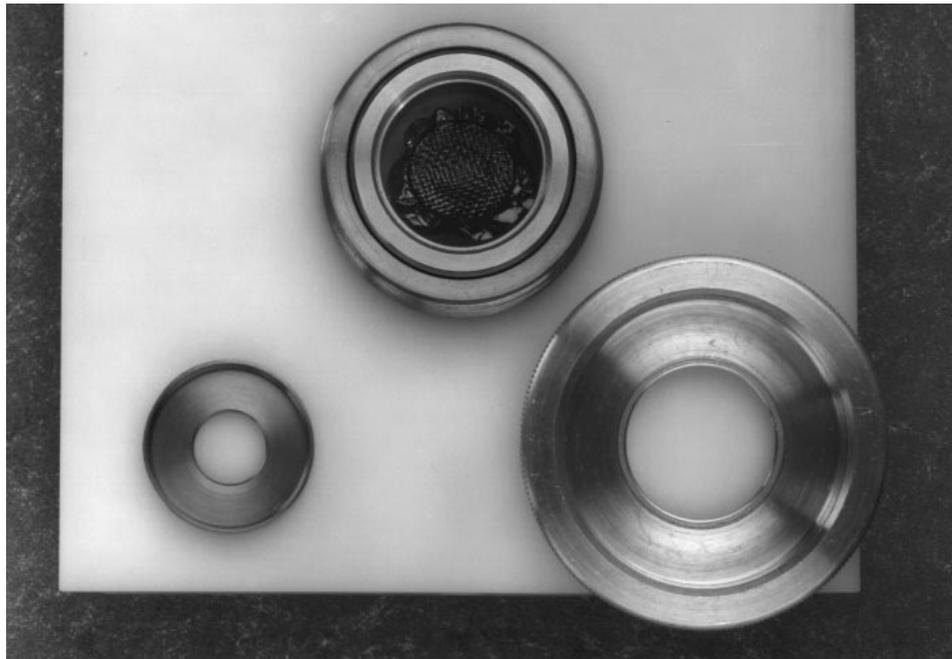


Fig. 4.12. View of used macrocarrier and stopping screen within disassembled microcarrier launch assembly after a bombardment.

15. Removal of spent rupture disk

Unscrew the rupture disk retaining cap from the gas acceleration tube. Remove the remains of the rupture disk (Figure 4.13). The next bombardment may now be performed (from step 7).

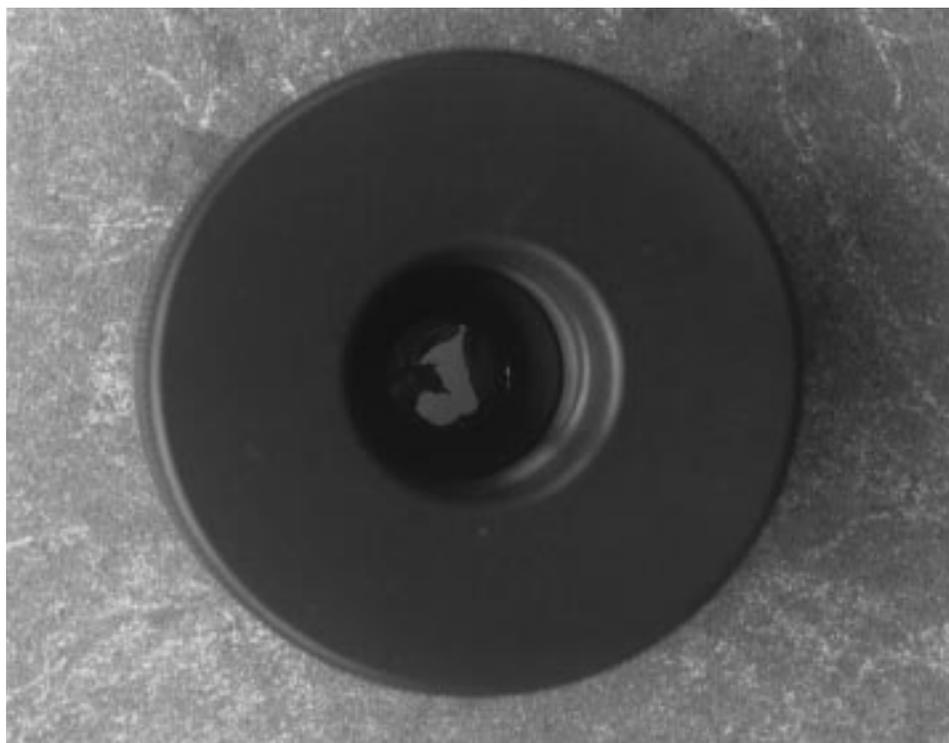


Fig. 4.13. View of spent rupture disk within the retaining cap.

4.3 Removal of Residual Helium Pressure—Shut Down

After completing all bombardment(s), remove the helium pressure from the PDS-1000/He system and close the helium cylinder valve. Perform the following steps to remove helium pressure from the system.

1. Close the helium cylinder valve and chamber door.
2. With at least 5 inches of Hg of vacuum in bombardment chamber, remove residual line pressure from the regulator, solenoid and PEEK tubing by activating the **FIRE** button.
3. Release the **FIRE** button on the apparatus, and remove all tension on the pressure adjustment screw of the helium regulator, turning counter-clockwise, until it turns freely.
4. Vent any residual vacuum from the bombardment chamber by setting the vacuum switch to the **VENT** position.

Section 5 Selection and Adjustment of System Bombardment Parameters

5.1 Overview—Matrix of Variables

Factors affecting bombardment efficiency are numerous, and interact in complex ways. All possible variables cannot be addressed in a single large factorial experiment. Instead, most users will find it sufficient to prioritize variables by the magnitude of their effects, optimizing them individually, and then testing their interactions on a more limited scale. Table 5.1 shows some examples of conditions used in published protocols for a variety of target materials.

Table 5.1. Cell Types, Settings, and Conditions

Cell Type	Growth Phase	Cell Density	Osmoticum	Vacuum (inches Hg)	Target Distance (cm)	Helium Pressure (psi)	Particle Size
Bacteria	Late log to early stationary	10 ⁸ –10 ⁹ per 100 mm plate	0.75 M sorbitol	29	6	1,100	M5 tungsten
Yeast	Early stationary	10 ⁸ –10 ⁹ per 100 mm plate	0.75 M sorbitol and 0.75 M manitol	28	6	1,300	0.6 μ gold
Algae	Log	10 ⁸ –10 ⁹ per 100 mm plate		29	6	1,300	0.6 μ gold
Plant							
• embryos	–	10 explants per 100 mm plate	None	28	6	1,300	1.0 μ gold
• callus or cell culture	Log	0.75 ml packed cell volume	None	28	9	1,100	1.0 μ gold
Subcellular Organelles	Mid-log	5 x 10 ⁷ per 100 mm plate	None	28	6	1,300	0.6 μ gold
Animal							
• tissue culture	Log	50–80% confluent on 35 mm plates	None	15	3	1,100	1.6 μ gold
• tissue sections	1 hr–4 days post-excision	400 μm sections	None	25	9	1,100	1.6 μ gold

5.2 Vacuum Level in Bombardment Chamber

The vacuum in the bombardment chamber reduces the frictional drag of the microcarriers as they are accelerated toward the target cells. The unit should be connected to a vacuum system (see previous section) that can evacuate the bombardment chamber to 28–29 inches Hg in less than 30 seconds. This level of vacuum is useful for most plant and microbial cells/tissues. Mammalian cells/tissue should be bombarded at approximately 15 inches Hg.

Helium gas enters the evacuated bombardment chamber once the rupture disk bursts. The use of low molecular weight helium minimizes the deceleration of the microcarriers as they pass through helium and also reduces the force of the gas shock wave that hits the target cells. This reduced impact will help minimize target tissue damage.

5.3 Helium Pressure / Rupture Disk Selection

Each of the nine different rupture disks available ruptures at a specific pressure, ranging in rating from 450 to 2,200 psi. The rupture pressure determines the power of the shock wave entering the bombardment chamber. Increasing helium pressure will increase particle acceleration and subsequent target tissue penetration by the DNA-coated microcarriers. Since the shock wave or resulting acoustic wave may cause damage to the target cells or tissue, use the lowest helium pressure used that gives high transformation efficiency.

Suggested starting helium pressure conditions for optimizing various biological systems are:

Cell type	Rupture disk pressure
Bacteria	1,100 psi
Fungi	1,300 psi
Yeast	1,300 psi
Plant cells/tissue	1,100 psi
Mammalian cells	1,100 psi

5.4 Solenoid Valve Adjustment

A factory pre-set metering valve (black knob) on the 3-way helium metering (solenoid) valve assembly controls the fill rate of the gas acceleration tube. The proper fill rate is set at the factory for a 1,550 psi rupture disk to burst within 12–15 seconds. A more rapid fill rate may result in what appears to be a lower burst pressure, due to a lag of needle movement in the oil-filled gauge. The metering valve (black knob) may be adjusted if desired. Adjust knob in small increments: a clockwise rotation of the black knob will lengthen the fill rate.

5.5 Vacuum Flow Rate Control Valves

Both valves are set at the factory to be fully open (counter clock-wise), for maximum flow rates. To decrease the rates, turn the valves (knob) clockwise until the desired evacuation and/or vent rate is achieved (see Figure 2.9).

5.6 Distance Between Rupture Disk and Macrocarrier

The effect of the gas shock wave on microcarrier velocities is determined in part by the gap between the rupture disk and the macrocarrier. The smaller the distance, the more powerful the effect of the gas shock wave on macrocarrier acceleration.

Three hexagonal gap adjustment tools of 1/8", 1/4", and 3/8" have been provided to reproducibly set the gap distance (Figure 2.4). A 1/4" distance between the rupture disk retaining cap and the macrocarrier cover lid is recommended when initially optimizing bombardment parameters.

5.7 Distance Between Macrocarrier and Stopping Screen

Macrocarrier flight instability increases with greater travel distance, therefore, the shortest travel distance is recommended. This is achieved by having the stopping screen support placed above both spacer rings inside the fixed nest when initially optimizing bombardment conditions. The macrocarrier flight distance can be adjusted by varying the position of the stopping screen support, as positioned by the spacer rings inside the fixed nest of the microcarrier launch assembly. The travel can be increased to a maximum of 16 mm in steps of 5 mm by varying spacer ring placement.

5.8 Distance Between Stopping Screen and Target Shelf (microcarrier flight distance)

One of the most important parameters to optimize is target shelf placement within the bombardment chamber. This placement directly affects the distance that the microcarriers travel to the target cells for microcarrier penetration and transformation.

Four target shelf levels are available in the bombardment chamber: level 1= 3 cm, level 2 = 6 cm, level 3 = 9 cm, and level 4 =12 cm below the stopping screen (Figure 5.1).

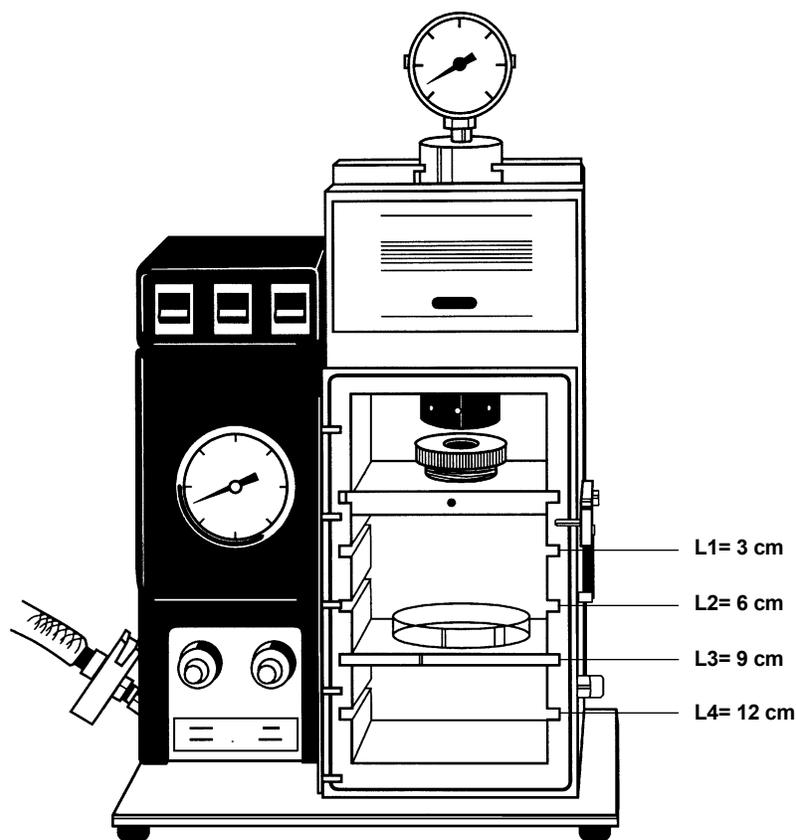


Fig. 5.1. Target shelf placement and corresponding target distances.

5.9 Microcarrier Selection

Types of Microcarrier

Two types of microcarriers are available, gold and tungsten. The density of these microcarriers is sufficient to penetrate a wide variety of cell and tissue types using the Biolistic PDS-1000/He acceleration system.

Five tungsten particle sizes are available. The tungsten particles are less expensive, but more irregular in shape and the size variation within a given particle size is much wider than gold. This size variation makes optimization using a certain particle size more difficult. Also, tungsten may be toxic to certain cell types and may oxidize to alter DNA binding.

Gold microcarriers are available in three sizes. Gold particles are more spherical and are more uniform in size within a given sample than tungsten microcarriers. Gold is biologically inert, but the DNA coating of this type of microcarrier is subject to more variation during precipitation than tungsten.

Microcarrier Size

The velocity of the chosen microcarrier increases with larger particle size. Increased microcarrier velocity has a major effect on particle penetration into target cells/tissue and thus transformation efficiency.

Recommended starting particle size/type for bombardment of various cell types is

Bacteria	0.7 μm (M5) tungsten
Yeast	0.6 μm gold
Algae	0.6 μm gold
Plant cells/tissue	1.0 μm gold
Animal cell cultures	1.6 μm gold
Sub-cellular organelles	0.6 μm gold

Table 5.1 and bulletins 1688 and 2015 give a detailed discussion of optimization parameters.

5.10 Preparation of Biological Material for Bombardment

Plant Cells

Many factors contribute to optimum performance in any system for plant gene transfer. Transformation procedures employing microcarrier bombardment are no exception, but the improved PDS-1000/He apparatus lends itself to rapid and easy modification, to tailor its performance to a particular application. Microcarrier bombardment also allows exploitation of a very broad range of explant tissue types. This permits optimization of a crucial aspect of the transformation system which usually is difficult to alter when transforming by other means. The following is a discussion of some of the critical parameters in the microcarrier transformation system, together with a brief examination of explant characteristics which affect transformation efficiency. This is not intended as a "how to" manual. Because of the extreme diversity of applications possible with the system, and the various types of tissues which can be employed, it is not possible to produce a universal protocol for all applications. Rather, this discussion is intended to aid the new user to design experiments.

Note: If you are a first-time user, it is important to be certain that your PDS-1000/He apparatus has been assembled correctly and is operating according to specifications. We suggest that you use the Yeast Optimization Kit (catalog number 170-3100) to be sure that microcarrier preparation has been carried out properly, and that microcarriers are being delivered uniformly. This also provides an opportunity to examine firsthand the pattern of particle delivery and gene expression in treated cells. This may save substantial time as you begin to optimize the system for your own purposes.

Choice of Explant Tissue

Unlike most other procedures for DNA transfer, microcarrier bombardment places few constraints on the types of tissues which can be treated. DNA can be delivered into essentially any cell or tissue which can be exposed sufficiently to allow particle penetration. Thus more emphasis usually can be given to issues such as the tissue's ability to regenerate plants, or its physiological suitability for gene regulation studies. There are, however, several aspects of the explant which contribute to the efficiency of the system.

An ideal tissue for microcarrier bombardment could perhaps be described as a cell monolayer, all cells of which are capable of expressing introduced genes, and all of which could, independently, divide and differentiate into functional plants. Microcarrier bombardment of such a tissue will likely be efficient because cells can be spread over a large area, thus efficiently capturing a large proportion of the particles delivered in each bombardment. Particles also have a high probability of penetrating such cells because they are not covered by overlying cell layers which would very likely intercept some of the particles delivered by each bombardment. Absolute transient expression levels are likely to be high, because of the number of cells affected. The low profile of cell clusters reduces the probability that the tissue will be moved about by the gas released in the shockwave during treatment. Such a tissue would also lend itself well to selection with antibiotics or herbicides, in part, at least, because it would have a favorable geometry for the establishment of a uniform concentration of the selective agent throughout the culture.

Some kinds of embryogenic suspension cultures approach this ideal. In fact, it is this type of tissue which has allowed great success in the microcarrier-mediated transformation of recalcitrant species such as maize [see, for example, Gordon-Kamm *et al.*, *The Plant Cell*, **2**, 603-618 (1990)]. Most explants, however, differ from this ideal in substantive ways. It is useful, though, to keep these characteristics in mind when choosing and preparing tissues for bombardment, whether taken from established cultures, or recently removed from donor plants. For example, tissue damage is typically most severe within the central area of bombardment where particle density is greatest. Cultured materials can usually be arranged to minimize this effect by placing them in a circle around the central area of bombardment. Even though some of the DNA-bearing particles fail to enter cells, this will often increase the absolute number of transformation events recovered from a single bombardment by reducing tissue damage. Similarly, explants from donor plants should be arrayed so as to capture as large a portion of particles as possible. Newly acquired explants can be dissected so that the cell types of interest are exposed as much as possible. The height of the tissue above the surface of its support medium can be minimized to reduce losses during treatment, which result from tissues with high profiles being blown about during bombardment. Finally, it may be useful to reduce the size of each explant or tissue mass to more closely approximate the geometry of suspension culture cells, in order to enhance the uniformity of selection.

Most tissues appear to vary with regard to their ability to express introduced genes following bombardment. For example, this effect is frequently observed as fluctuations in the frequency or intensity of transient expression of β -glucuronidase (GUS) constructions among different tissues, or even among cells in different areas of a single explant. It can be caused in part by unevenness in particle distribution during preparation of macrocarriers, but apparently can also be caused by heterogeneity of the tissue itself. In intact explants it is not uncommon to see variation in expression associated with the physiological age of the explant. Although the causes of these phenomena are unknown, it is important to realize that such effects exist, and can be a source of serious confounding in experiments designed to choose tissues for use in transformation experiments, and to optimize bombardment parameters.

Optimization of Bombardment Parameters—Transient expression vs stable transformation

The goal of many microcarrier bombardment experiments is to develop protocols for efficient production of transgenic plants. The frequency at which stable transformation events are recovered is the ultimate criterion by which these protocols must be measured. However, measurement of transient expression (*i.e.*, expression of newly introduced DNA sequences after a relatively brief period following bombardment) provides rapid feedback, and can be an invaluable aid in the determination of effective bombardment parameters for a specific tissue.

A precise quantitative relationship between the level or frequency of transient expression observed in a bombarded tissue, and the frequency at which stable transformation of cells occurs within that tissue may not exist. Transient expression is, however, a very useful indicator of the efficiency of DNA delivery, and can help to define conditions required to deliver DNA into specific cell layers. In this way, measurement of transient expression allows rapid determination of a set of parameters which permit delivery of DNA into tissues of interest, and provides critical information about which parameters have the greatest effect.

Choice of reporter gene and assay procedures for optimization of bombardment must be determined by the particular question to be addressed. Histochemical assay of β -glucuronidase expression in bombarded cells is currently the most commonly used measurement of the frequency of transient expression events in bombarded tissues. It is also ideal for determining which cell types in a heterogeneous tissue are capable of expressing the introduced sequence. This procedure is not useful, however, for determining the absolute expression levels produced by a particular set of bombardment parameters. For such measurements, it is

more appropriate to use an extractive assay procedure in which expression level can be measured, for example, as a function of enzyme activity per unit of total protein. Several excellent markers (and associated assay procedures) are available for this type of study. These include luciferase and β -glucuronidase measured by a fluorometric process, and neomycin phosphotransferase, for which a commercial ELISA assay has recently been marketed.

Yeast/Bacteria/Microbes

Cells which grow in suspension should be grown to late log to early stationary phase in media under the appropriate conditions. Pellet the cells and resuspend in sterile water. Estimate the cell density by determining the absorbance (for yeast, 1 O.D.₆₀₀ is approximately 2.8×10^9 cells/ml; for *E. coli*, 1 O.D.₆₀₀ is approximately 1.0×10^9 cells/ml). Spread 1×10^8 yeast or 2×10^9 *E. coli* on 100 mm Petri plates containing the appropriate media. In addition to the necessary supplements, the media for growing yeast should contain 0.75 M mannitol and 0.75 M sorbitol, while media for growing bacteria should contain 0.75 M sorbitol. The plates should be allowed to air dry briefly, but used within 1 or 2 hours. Transformation efficiency decreases if the plates are allowed to sit even for several hours.

Mammalian Cells

Cells in culture

Both primary and established animal cell lines are transformable using particle bombardment. Because optimum bombardment conditions are usually found when mammalian cells are placed at the top level of the bombardment chamber, the cells can be inoculated into 35 mm tissue culture plates. Only if cells are to be bombarded at Target Level 2 or farther should larger plates be used.

Cells which grow as attached cultures should be inoculated into tissue culture plates one day prior to bombardment so that they are in log phase and 50–80% confluent at the time of transformation. Cells which grow as suspension cultures should be inoculated from log phase cultures onto plates which have been aseptically coated by adsorption with Cell-Tak (Collaborative Biomedical Products, Bedford, MA) according to the manufacturer's instructions. The cells should be allowed to attach for 1 or 2 hours and should be approximately 75% confluent at the time of bombardment.

Immediately prior to bombardment, tissue culture media should be aspirated so that the cells are directly exposed to air. Cells to be bombarded are placed in the bombardment chamber centered under the stopping screen, the chamber door closed, the chamber evacuated to the proper vacuum, and the cells bombarded. After bombardment, add fresh tissue culture media to the cells and incubate for the appropriate time.

Organ cultures

Animal organs may be transformed by particle bombardment 1 hour to several days after preparation. Tissues should be approximately 400 μ m thick sections and no more than 1 cm square. The tissue may be grown in cell culture inserts containing a permeable membrane. The tissue section (in the cell culture insert) is placed in a Petri plate containing 1% agar, then transferred to the bombardment chamber. After bombardment, the insert containing the tissue can be transferred back to the growth chamber.

Section 6 Troubleshooting

6.1 Rupture Disk Bursts at Incorrect Pressure

Probable Cause	Action
1. The helium flow rate is too fast.	Note the helium burst pressure by observing the helium pressure gauge at the top of the main unit. A factory pre-set metering valve (black knob) on the external 3-way helium metering (solenoid) valve assembly controls the fill rate of the gas acceleration tube. The proper fill rate is 12–15 seconds to burst a rupture disk when the helium pressure is set 200 psi over the rated rupture pressure. A more rapid fill rate may result in what appears to be a lower burst pressure, due to a lag of needle movement in the oil-filled gauge. The metering valve (black knob) may be adjusted if desired. Adjust knob in small increments: a clockwise rotation of the black knob will lengthen the fill rate.
2. The improper disk was inserted or two disks are sandwiched together.	Replace with single, desired psi rupture disk.
3. The retaining cap was not securely tightened.	This is frequently the case when the disk slips out of the retaining cap. It is important to screw the cap onto the end of the gas acceleration tube by hand and then use the torque wrench.
4. The rupture disks are not clean.	Use tweezers or disc-vac to handle the disks after briefly rinsing in isopropanol. Do not handle with bare or gloved hands (grease and powder source).
5. The sealing edge of the Gas Acceleration Tube is not clean.	Occasionally adhesive from the rupture disks (laminated disks) may build up on the edge of this part. This can be cleaned by using a mild, non-abrasive detergent, like dish soap, to remove the accumulated adhesive.
6. The sealing edge (inside bombardment chamber) of the Gas Acceleration Tube is damaged.	The edge of this part is susceptible to physical damage if the retaining cap is tightened without a rupture disk in place. If damage has occurred such that the disk will not rupture, you must replace this part.
7. Rupture disk is defective.	Replace.

6.2 Stopping Screen Forced Through Screen Support Ring

Probable Cause	Action
1. The screen has been “unseated” from support	Carefully insert the Microcarrier Launch Assembly into bombardment chamber to prevent movement of stopping screen. This screen movement can also occur if you close the bombardment chamber door too forcefully.
2. The stopping screen support ring is damaged.	Examine the support ring for evidence of scratches or dents on the support ring that might prevent the screen from seating properly.

6.3 Excessive Gas Usage

Probable Cause	Action
1. Leaking helium fitting in the system	Isolate the components from the He cylinder to the solenoid assembly by completely closing the metering valve (turning the knob counterclockwise until it stops.) Ensure that the regulator gauge maintains pressure over time. “Soap Test” the fittings by spraying a film of soapy water at the site of the fittings and looking for bubble formation. Tighten any fittings that appear to leak. Remember to close main cylinder valve of helium tank when not in use.
2. Damaged sealing edge on Gas Acceleration Tube (scratch or dent in metal).	Replace Gas Acceleration Tube.

6.4 Chamber Will Not Hold Vacuum

Probable Cause	Action
1. Door gasket does not seal	Test the door gasket (the large rubber O-ring around the interior perimeter of the bombardment chamber door) by spraying the outer edge of the closed door with soapy water while attempting to hold a vacuum. If the soap bubbles, replace O-ring.
2. Vacuum connection at the back of the bombardment chamber is not sealed	Check for leakage (poorly fitting components) from the clamp assembly located at the rear of the unit.

6.5 Sample Damage from Gas Pressure Wave

Probable Cause	Action
1. Macrocarrier not inserted	Examine and replace
2. Microcarrier Launch Assembly not inserted into bombardment chamber	Examine and replace
3. Incorrect sample shelf position	Reposition
4. Stopping screen absent, improperly positioned or fails to function within the Microcarrier Launch Assembly	Examine and replace
5. Incorrect assembly of Microcarrier Launch Assembly components	Examine and reassemble

6.6 Unit Will Not Pressurize Gas Acceleration Tube

Probable Cause	Action
1. Rupture disk not in position or not sealed by retaining cap.	Insert disk properly
2. Main cylinder valve of Helium tank not open	Adjust according to Section 3.1
3. Insufficient vacuum	Examine vacuum connections (see Section 3.2)
4. Helium system leak	<p>Isolate leaks in the helium pressure regulator by closing the helium cylinder valve and turning the pressure-adjusting screw one turn counter-clockwise, then note the following:</p> <ul style="list-style-type: none"> • If the high pressure gauge reading drops, there is a leak in the cylinder valve, inlet fitting, or high pressure gauge. • If the low pressure gauge drops, there is a leak in the valve, hose, hose fitting, outlet fitting or low pressure gauge. • If the high pressure gauge drops, and at the same time the low pressure gauge rises, there is a leak in the regulator seat. <p>Defective regulators should be returned to Bio-Rad Laboratories for repair. Use of a damaged regulator, or one that has been repaired improperly is a safety hazard.</p>
5. Helium tank regulator not adjusted correctly	Adjust according to Section 3.1
6. Solenoid 3-way metering valve assembly not properly connected to main unit	Connect according to Section 3.1 and 5.4
7. Solenoid 3-way metering valve closed	Adjust according to Section 3.1 and 5.4

Section 7 Product Information

7.1 Biolistic Particle Delivery System

Catalog Number	Product Description
165-2250LEASE	PDS-1000/He Lease - 1st Year - Academic
165-2251LEASE	PDS-1000/He Lease - Additional Year - Academic
165-2252LEASE	PDS-1000/He Lease - 5 Year - Academic
165-2253LEASE	PDS-1000/He Lease - 1st Year - Industrial
165-2254LEASE	PDS-1000/He Lease - Additional Year - Industrial
165-2255LEASE	PDS-1000/He Lease - 5 Year - Industrial
165-2248LEASE	PDS-1000/He Expired Lease-to-Purchase Conversion
165-2256	Particle Delivery Research License
165-2257	PDS-1000/He , for nonprofit organizations
165-2258LEASE	PDS-1000 Helium Retrofit Kit , for converting the PDS-1000 to the PDS-1000/He
165-2259	Voltage Converter , for countries with 220 V or 240 V
165-2262	0.6 μ Gold Microcarriers , 0.25 g
165-2263	1.0 μ Gold Microcarriers , 0.25 g
165-2264	1.6 μ Gold Microcarriers , 0.25 g
165-2265	Tungsten M-5 Microcarriers , ~ 0.4 μ , 6 g
165-2266	Tungsten M-10 Microcarriers , ~ 0.7 μ , 6 g
165-2267	Tungsten M-17 Microcarriers , ~ 1.1 μ , 6 g
165-2268	Tungsten M-20 Microcarriers , ~ 1.3 μ , 6 g
165-2269	Tungsten M-25 Microcarriers , ~ 1.7 μ , 6 g
165-2278	500 Optimization Kit , includes 0.25 g each of 0.6 μ , 1.0 μ , and 1.6 μ gold microcarriers, 100 each of 9 rupture disks, 500 macrocarriers, 500 stopping screens
165-2280	500 Standard Pressure Kit , with 1.0 μ gold + 450 psi disks, includes 0.5 g 1.0 μ gold, 500 450 psi rupture disks, 500 macrocarriers, 500 stopping screens
165-2281	1,500 Standard Pressure Kit , with 1.0 μ gold + 650 psi disks
165-2282	500 Standard Pressure Kit , with 1.0 μ gold + 900 psi disks
165-2283	500 Standard Pressure Kit , with 1.0 μ gold + 1,100 psi disks
165-2284	500 Standard Pressure Kit , with 1.0 μ gold + 1,350 psi disks
165-2285	500 Standard Pressure Kit , with 1.0 μ gold + 1,550 psi disks
165-2286	500 Standard Pressure Kit , with 1.0 μ gold + 1,800 psi disks
165-2287	500 Standard Pressure Kit , with 1.0 μ gold + 2,000 psi disks
165-2288	500 Standard Pressure Kit , with 1.0 μ gold + 2,200 psi disks
165-2290	500 Standard Pressure Kit , with 1.6 μ gold + 450 psi disks, includes 0.5 g 1.6 μ gold, 500 450 psi rupture disks, 500 macrocarriers, 500 stopping screens
165-2291	500 Standard Pressure Kit , with 1.6 μ gold + 650 psi disks
165-2292	500 Standard Pressure Kit , with 1.6 μ gold + 900 psi disks
165-2293	500 Standard Pressure Kit , with 1.6 μ gold + 1,100 psi disks
165-2294	500 Standard Pressure Kit , with 1.6 μ gold + 1,350 psi disks
165-2295	500 Standard Pressure Kit , with 1.6 μ gold + 1,550 psi disks

Catalog Number	Product Description
165-2296	500 Standard Pressure Kit , with 1.6 μ gold + 1,800 psi disks
165-2297	500 Standard Pressure Kit , with 1.6 μ gold + 2,000 psi disks
165-2298	500 Standard Pressure Kit , with 1.6 μ gold + 2,200 psi disks
165-2322	Macrocarrier Holders , set of 5
165-2323	Disk-Vac
165-2326	450 psi Rupture Disks , 100
165-2327	650 psi Rupture Disks , 100
165-2328	900 psi Rupture Disks , 100
165-2329	1,100 psi Rupture Disks , 100
165-2330	1,350 psi Rupture Disks , 100
165-2331	1,550 psi Rupture Disks , 100
165-2332	1,800 psi Rupture Disks , 100
165-2333	2,000 psi Rupture Disks , 100
165-2334	2,200 psi Rupture Disks , 100
165-2335	Macrocarriers , 500
165-2336	Stopping Screens , 500
170-3100	Yeast Optimization Kit , includes <i>S. cerevisiae</i> , YEp352 DNA, CaCl ₂ , spermidine, glucose, culture medium, plating medium†

Consumables for the Biolistic PDS-1000 System (gunpowder-discontinued)

165-2275	500 Accelerator Kit , with 1.0 μ gold, includes 0.5 g gold, 500 accelerators, 500 macrocarriers, 500 stopping plates
165-2276	500 Accelerator Kit , with 1.6 μ gold
165-2277	Cleaning Patches, 60

7.2 Spare Parts for Helium Biolistic System Assembly (PDS-1000/He)

Spare Part Number	Product Description
Plastic Case Version	
100-9110	Macrocarrier Cover Lid
165-2322	Microcarriers , 5
165-2323	Disk-Vac
800-3007	Torque Lever Assembly
800-3013	Solenoid Valve Cord Assembly
800-3018	Gas Acceleration Tube Assembly
800-3030	Solenoid Valve
800-3036	2.5 ft PEEK Tubing Assembly
800-3039	6 ft PEEK Tubing Assembly
900-5156	Vacuum Switch Replacement Kit
910-0074	Acceleration Tube O-Ring
910-0757	Retaining Spring
910-1201	Set Screw
910-4192	Flow Control Needle Valve
910-9500	Cylinder Pressure Regulator

Spare Part

Spare Part Number	Product Description
910-9511	Helium Pressure Gauge
910-9515	Vacuum Gauge
910-9529	Wrench, Open End, 1/4" x 5/16"

Metal Case Version (Discontinued)

800-3014	Complete Door Assembly
910-9549	Complete Filter Housing
910-9526	Filter Housing O-Ring

Gunpowder Version (Discontinued)

910-9538	Accelerators, 100
920-3009	Locking Plug
910-0019	Perimeter Seal Door O-Ring
920-3061	Pin Loader
910-9522	Tornado Brush
910-9530	Cleaning Rod
910-9531	Dud Remover Rod
920-3022	Large Rubber Stopping Washer
920-3023	Steel Blue Stopping Disk
920-3026	Small Rubber Stopping Washer
920-3035	Latch, Door 2
920-3058	Mid Seal Door O-Ring
920-3007	Rupture Disk Retainer Cap
920-3008	Microcarrier Launch Shelf
920-3010	Adjustable Nest
920-3011	Fixed Nest
920-3013	Stopping Screen Holder
920-3014	5 mm Spacer Ring
920-3015	Microcarrier Cover Lid
920-3016	1/8 inch Gap Adjustment Tool
920-3017	1/4 inch Gap Adjustment Tool
920-3018	3/8 inch Gap Adjustment Tool
920-3060	Petri Dish Holder
920-3068	18 inch Vacuum Hose
920-3085	Red Seals, 500

Refer to the Bio-Rad catalog for more details or contact your local Bio-Rad office.

† Shipped on dry ice; perishable freight charges not included.

Biolistic is a registered trademark of E. I. du Pont de Nemours and Company. The Biolistic technology is exclusively licensed by Du Pont to Bio-Rad Laboratories.

Section 8 Appendices

8.1 Cleaning the PDS-1000/He Device

1. Chamber
Clean the chamber with 70% ethanol. Allow time for drying.
2. Gas Acceleration tube
 - A. Unscrew the rupture disk retaining cap from the gas acceleration tube

Do not autoclave or flame sterilize the PDS-1000/He. Autoclaving will wet electrical components, thus creating a potential electrical shock hazard.
 - B. Disconnect the tubing assembly from the gas inlet on the rear of the acceleration tube (See Figure 1.2). Remove the screws from the hold-down brackets on the top of the PDS-1000/He. Two sets are located on either side of the acceleration tube, and CAREFULLY remove the tube from the top of the sample chamber.
3. The acceleration tube can be sterilized by partially immersing in 70% ethanol followed by drying in a sterile environment. Do not autoclave the gas acceleration tube. Do not immerse the pressure gauge in the sterilizing solution.

8.2 Metal Case Version

The metal chamber design has been offered by Bio-Rad Laboratories since the product line was acquired from E.I du Pont de Nemours in 1991. The original metal case design is capable of operating either the gunpowder acceleration mechanism or the helium acceleration mechanism (Figure 8.1). The gunpowder acceleration delivery system has been discontinued. The current plastic case version of the PDS-1000/He unit was introduced in March 1995 and only operates with helium. The target distances and internal dimensions of the plastic case are equivalent to those found in the metal case when used with helium acceleration mechanism.



Fig. 8.1. Front view of Metal Case version of PDS-1000/He instrument.

Aspects unique to the metal chamber design

- Metal bombardment chamber.
- Top slots of the metal case. These are used only for the target shelf placement when using gunpowder acceleration system.
- Door construction:
Gasket design changed. Portions removed that were dedicated to use with gunpowder acceleration.
Door brace added in plastic case design for additional safety.
- External vacuum filter assembly components. Located on rear and side of metal case unit. These components are used to trap spent gunpowder in thin oil film inside trap after bombardment. See rear view (Figure 8.2) for identification of components.
- Helium tubing-stainless steel in metal case version. Composed of PEEK plastic in the plastic case version.
- Power cord is permanently attached in the metal case design. Power cord is detachable in the plastic case design.

Attachment/removal of vacuum filter assembly for the metal case PDS-1000/He unit

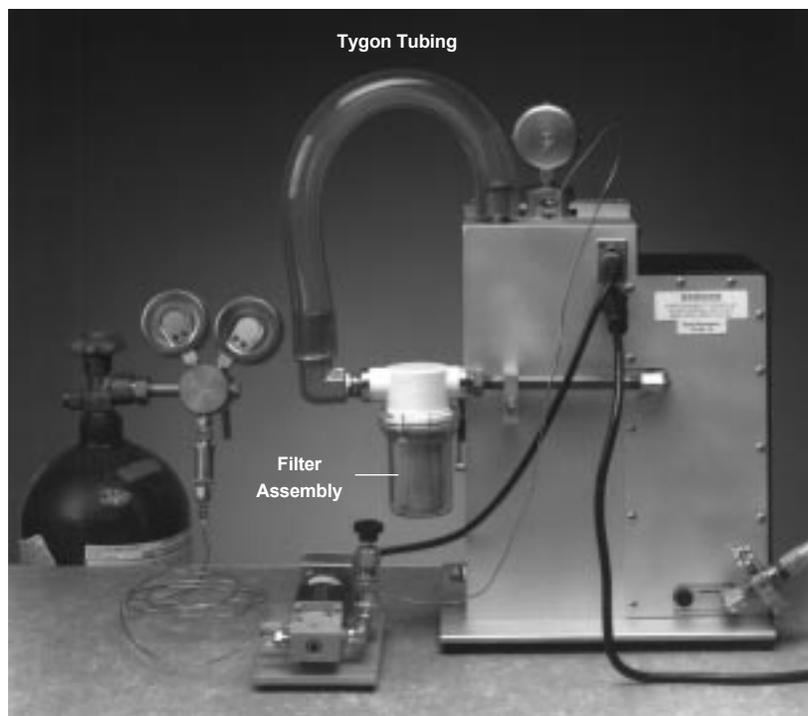


Fig. 8.2. Rear view of metal case version of PDS-1000/He system.

Wrap PTFE tape on the threaded brass vacuum filter fitting on the rear of the PDS-1000/He unit and attach the filter assembly (Figure 8.2). The vacuum filter assembly includes a plastic filter housing and brass filter hose fitting, plus the Tygon tubing. Turn the filter assembly approximately four complete turns, ending so the filter housing is parallel to the unit and the Tygon tubing is facing upward.

Push the open end of the Tygon tubing onto the baffle located on top of the PDS-1000/He unit.

Note: The filter assembly must be firmly supported as the Tygon tubing is being pushed onto the baffle on top of the unit (or onto the brass filter hose fitting) to avoid cracking the plastic filter housing.

Table 8.1 Sterilization and Decontamination Compatibility

Method	Material						
	Aluminum	Bunan	Delrin®	Neoprene	Nylon	Polycarbonate	Polyethylene
Sterilization							
Autoclave (121 °C)	U	M	M	U	U	M	M
UV Irradiation	S	*	S	*	M	U	U
Ethylene Oxide	S	U	S	*	S	S	S
Formaldehyde (gas)	S	S	S	S	S	S	S
Gluteraldehyde (2%)	S	S	S	S	S	S	S
Biological Disinfection							
Alcohol (70%)	S	S	M	S	S	M	S
Hypochlorite (5%)	U	M	U	M	S	S	S
Gluteraldehyde (2%)	S	S	S	S	S	S	S
Phenolic Derivatives	U	U	U	U	U	U	S
Formaldehyde	M	M	S	S	S	S	S

Key: S = Satisfactory; M = Moderate attack, may be satisfactory, suggest testing;
 U = Unsatisfactory, not recommended; *Performance unknown, suggest testing.

Method	Material					
	Polyvinyl Chloride	PTFE	Stainless Steel	Tygon	Acrylic	Painted Surfaces
Sterilization						
Autoclave (121 °C)	M	S	S	M	U	U
UV Irradiation	S	S	S	S	U	S
Ethylene Oxide	S	S	S	S	S	S
Formaldehyde (gas)	S	S	S	S	S	S
Gluteraldehyde (2%)	S	S	S	S	S	S
Biological Disinfection						
Alcohol (70%)	M	S	M	M	M	S
Hypochlorite (5%)	U	S	M	M	S	S
Gluteraldehyde (2%)	S	S	S	S	S	S
Phenolic Derivatives	U	S	U	U	U	S
Formaldehyde	M	S	S	M	S	S

Key: S = Satisfactory; M = Moderate attack, may be satisfactory, suggest testing;
 U = Unsatisfactory, not recommended; *Performance unknown, suggest testing.

8.3 Specifications

General Specifications, PDS-1000/He Biolistic Particle Delivery System

Dimensions	29 (width) x 25.5 (depth) x 47.5 (height) cm
Construction	Aluminum, ABS plastic and acrylic chassis
Weight	15 kg
Electrical input voltage	100-120 VAC, 50- 60 Hz
Maximum current	<5 Amps
Mechanical	
Fuse	6.3 A, 250 V, 5 x 20 mm
Vacuum	<0.4 inches Mercury/minute leakage
Over-Pressure	0.5 psi relief valve, self-resetting
Environmental	
Operating	32 °F (0 °C) to 95 °F (35 °C) temperature, 0-95% non-condensing humidity
Storage	32 °F (0 °C) to 158 °F (70 °C) temperature, 0-95% non-condensing humidity

Sole Source Specifications, PDS-1000/He Biolistic Particle Delivery System

The following are the sole source specifications for the Biolistic PDS-1000/He unit:

1. The system is covered by patents numbers 161807, 670771, 877619, 074652, IP-0844, 07/303/503, 07/529989, 437848, 4945050, 5036006, 5100792, and 621561 together with all continuations, divisionals, and continuations-in-part applications, any patents that issue on applications and reissues thereof. The Biolistic PDS-1000/He instrument is available for sale or lease accordingly. Biolistic is a registered trademark of E. I. du Pont de Nemours and Company. The Biolistic technology is exclusively licensed to Bio-Rad Laboratories.
2. A system capable of generating high pressure gas delivery with a mechanism for delivering an instantaneous cold gas shock wave into an enclosure where the gas shock is released, contained and vented.
3. The unit contains a throat region in the high pressure gas delivery system that allows for interchangeable inserts (rupture disks) which translates the cold gas shock into acceleration of microprojectiles coated with biological molecules for delivery into diverse target cells/tissue without killing the cells and/or tissue. Nine interchangeable disks are available in 450, 650, 900, 1,100, 1,350, 1,550, 1,800, 2,000 and 2,200 pounds per square inch (psi) pressure.
4. Employs a method for launching the biologically coated microcarriers from the target side of a planar surface of a flexible plastic sheet that is positioned to be accelerated by the gas shock wave such that it will move freely in the direction of the target cells until restrained by a barrier. The barrier permits the microcarriers to move toward the target cells and also effectively baffles the gas shock wave to deflect some of the force of the cold gas shock wave away from the target cells with sufficient acceleration of microprojectiles coated with biological molecules for delivery into diverse target cells/tissue without killing the cells and/or tissue.

8.4 Performing a Bombardment

Quick Guide

Before the Bombardment

1. Select/adjust bombardment parameters for Gap distance between rupture disk retaining cap and microcarrier launch assembly. Placement of stopping screen support in proper position inside fixed nest of microcarrier launch assembly
2. Check helium supply (200 psi in excess of desired rupture pressure).
3. Clean/sterilize:
Equipment: rupture disk retaining cap, microcarrier launch assembly
Consumables: macrocarriers/macrocarrier holders
4. Wash microcarriers and resuspend in 50% glycerol
5. Coat microcarriers with DNA and load onto sterile macrocarrier/macrocarrier holder the day of experiment

Firing the Device

1. Plug in power cord from main unit to electrical outlet.
2. Power ON.
3. Sterilize chamber walls with 70% ethanol.
4. Load sterile rupture disk into sterile retaining cap.
5. Secure retaining cap to end of gas acceleration tube (inside, top of bombardment chamber) and tighten with torque wrench.
6. Load macrocarrier and stopping screen into microcarrier launch assembly.
7. Place microcarrier launch assembly and target cells in chamber and close door.
8. Evacuate chamber, hold vacuum at desired level (minimum 5 inches of mercury).
9. Bombard sample: Fire button continuously depressed until rupture disk bursts and helium pressure gauge drops to zero.
10. Release Fire button.

After the Bombardment

1. Release vacuum from chamber.
2. Target cells removed from chamber.
3. Unload macrocarrier and stopping screen from microcarrier launch assembly.
4. Unload spent rupture disk.
5. Remove helium pressure from the system (after all experiments completed for the day)



**Bio-Rad
Laboratories**

**Molecular
Bioscience Group**

2000 Alfred Nobel Drive
Hercules, California 94547
Telephone (510) 741-1000
Fax: (510) 741-5800

Australia, Bio-Rad Laboratories Pty Limited, Block Y Unit 1, Regents Park Industrial Estate, 391 Park Road, Regents Park, NSW 2143 • Phone 02-9914-2800 • Fax 02-9914-2888
Austria, Bio-Rad Laboratories Ges.m.b.H., Auhofstrasse 78D, 1130 Wien • Phone (1) 877 89 01 • Fax (1) 876 56 29
Belgium, Bio-Rad Laboratories S.A./N.V., Begoniastraat 5, 9810 Nazareth Eke • Phone 09-385 55 11 • Fax 09-385 65 54
Canada, Bio-Rad Laboratories (Canada) Ltd., 5671 McAdam Road, Mississauga, Ontario L4Z 1N9 • Phone (905) 712-2771 • Fax (905) 712-2990
China, Bio-Rad Laboratories, 14, Zhi Chun Road, Hai Dian District, Beijing 100088 • Phone (01) 2046622 • Fax (01) 2051876
Denmark, Bio-Rad Laboratories, Symbion Science Park, Fruebjergvej 3, DK-2100 Copenhagen • Phone 39 17 9947 • Fax 39 27 1698
Finland, Bio-Rad Laboratories, Business Center Länsikeskus, Pihatörmä 1A SF-02240, Espoo, • Phone 90 804 2200 • Fax 90 804 1100
France, Bio-Rad S.A., 94/96 rue Victor Hugo, B.P. 220, 94 203 Ivry Sur Seine Cedex • Phone (1) 49 60 68 34 • Fax (1) 46 71 24 67
Germany, Bio-Rad Laboratories GmbH, Heidemannstraße 164, D-80939 München/Postfach 450133, D-80901 München • Phone 089 31884-0 • Fax 089 31884-100
India, Bio-Rad Laboratories, C-248 Defence Colony, New Delhi 110 024 • Phone 91-11-461-0103 • Fax 91-11-461-0765
Italy, Bio-Rad Laboratories S.r.l., Via Cellini, 18/A, 20090 Segrate Milano • Phone 02-21609 1 • Fax 02-21609-399
Japan, Nippon Bio-Rad Laboratories, 7-18, Higashi-Nippori 5-Chome, Arakawa-ku, Tokyo 116 • Phone 03-5811-6270 • Fax 03-5811-6272
The Netherlands, Bio-Rad Laboratories B. V., Fokkerstraat 10, 3905 KV Veenendaal • Phone 0318-540666 • Fax 0318-542216
New Zealand, Bio-Rad Laboratories Pty Ltd., P. O. Box 100-051, North Shore Mail Centre, Auckland 10 • Phone 09-443 3099 • Fax 09-443 3097
Pacific, Bio-Rad Laboratories, Unit 1111, 11/F., New Kowloon Plaza, 38, Tai Kok Tsui Road, Tai Kok Tsui, Kowloon, Hong Kong • Phone 7893300 • Fax 7891257
Singapore, Bio-Rad Laboratories (Singapore) Ltd., 221 Henderson Rd #05-19, Henderson Building, Singapore 0315 • Phone (65) 272-9877 • Fax (65) 273-4835
Spain, Bio-Rad Laboratories, S. A. Avda Valdelaparra 3, Pol. Ind. Alcobendas, E-28100 Alcobendas, Madrid • Phone (91) 661 70 85 • Fax (91) 661 96 98
Sweden, Bio-Rad Laboratories AB, Gärdsvägen 7D, Box 1276, S-171 24 Solna • Phone 46-(0)8-735 83 00 • Fax 46-(0)8-735 54 60
Switzerland, Bio-Rad Laboratories AG, Kanalstrasse 17, Postfach, CH-8152 Glattbrugg • Phone 01-809 55 55 • Fax 01-809 55 00
United Kingdom, Bio-Rad Laboratories Ltd., Bio-Rad House, Maylands Avenue, Hemel Hempstead, Herts HP2 7TD • Free Phone 0800 181134 • Fax 01442 259118