



Gene Shot Control Cartridges

**Catalog Number
165-2244**



Specifications

Contents Twelve cartridges, for use in the Helios Gene Gun, packed with a desiccant capsule in a cartridge storage vial. Each cartridge contains 0.5 mg of 1.6 μ gold (catalog number 165-2264) coated with 1 μ g (~0.20 pmole) of a plasmid encoding a β -galactosidase reporter gene and with 0.3 μ g (~0.076 pmole) of a plasmid encoding a luciferase reporter gene. These reporter genes are expressed from a constitutive high-level promoter that is active in all animal cells.

Storage Store tightly capped, with parafilm around the top, at 4 °C with a desiccant pellet (supplied). **Before opening, allow vial to come to room temperature.**

Quality control

The biological efficiency of the cartridges was determined by expression of β -galactosidase and luciferase in cell extracts of CHO cells 24 hr post-delivery using the Helios Gene Gun with the screen (catalog number 165-2425). CHO cells attached to 35 mm tissue culture plates bombarded at 200 psi with these cartridges express $\geq 5 \times 10^6$ RLU of luciferase in a 10 sec counting period per 10^5 cells and produce ≥ 20 nmole of 4-methylumbelliferone (from 4-methylumbelliferyl- β -D-galactopyranoside) per 10^5 cells when assayed with the FluorAce™ β -galactosidase Assay Kit (catalog number 170-3150).