




Biolistic Particle Delivery Systems

Factors Affecting Transformation	Helios Gene Gun System	PDS-1000/He System	PDS-1000/He System With Hepta™ Adaptor
			
Experimental conditions	In situ, in vitro, in vivo, ex vivo	In vitro, ex vivo, in vivo (plants)	In vitro, ex vivo, in vivo (plants)
Target area	Small (2 cm ²)	Large (40 cm ²)	Largest (~75 cm ²)
Pressure range	100–600 psi	450–2,200 psi	450–2,200 psi, reduced by 7-way
Target type	Animals: Any tissue exposed to barrel (skin, organs); cell, explant, and organ culture Plants: Field and greenhouse use, plant cell culture, explants Yeast, bacteria, other microbes	Animals: Cell and organ culture Plants: Small intact plants, plant cell culture, explants Yeast, bacteria, other microbes Organelles (chloroplasts, mitochondria, etc.)	Animals: Cell and organ culture Plants: Cells with thin cell walls Yeast, bacteria, other microbes



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Web site www.bio-rad.com **USA** 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 01 877 89 01 **Belgium** 09 385 55 11 **Brazil** 55 11 5044 5699
Canada 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 420 241 430 532 **Denmark** 44 52 10 00 **Finland** 09 804 22 00
France 01 47 95 69 65 **Germany** 089 31 884 0 **Greece** 30 210 9532 220 **Hong Kong** 852 2789 3300 **Hungary** 36 1 459 6100 **India** 91 124 4029300
Israel 03 963 6050 **Italy** 39 02 216091 **Japan** 03 6361 7000 **Korea** 82 2 3473 4460 **Mexico** 52 555 488 7670 **The Netherlands** 0318 540666
New Zealand 64 9 415 2280 **Norway** 23 38 41 30 **Poland** 48 22 331 99 99 **Portugal** 351 21 472 7700 **Russia** 7 495 721 14 04
Singapore 65 6415 3188 **South Africa** 27 861 246 723 **Spain** 34 91 590 5200 **Sweden** 08 555 12700 **Switzerland** 061 717 95 55
Taiwan 886 2 2578 7189 **Thailand** 800 88 22 88 **United Kingdom** 020 8328 2000