

## Recommended Biolistic System by Cell Types

Transfection Technologies	Plants										Bacteria	Fungi			Animals					
	Plant Cell Suspension Cultures	Plant Cell Adherent Cultures	Plant Cell Callus or Explants	Pollen	Leaves in Situ	Detached Leaves	Meristems in Situ	Detached Meristems	Gram-Positive Bacteria	Gram-Negative Bacteria		Yeast	Other Fungi	Skin or Organs in Vivo	Organ Cultures	Ex Vivo Cells or Tissues (e.g., blood)	Suspension Cell Cultures	Mammalian Cells — Gene Knockdowns	Adherent Cells, Trypsinized	Adherent Cells, Attached
<b>Biolistic Particle Delivery</b>																				
Helios® gene gun system	•	•	•	•	•	•	•	•	•	•	•	•	•	•	•	•	•	•	•	•
PDS-1000/He™ system	•	•	•	•	•	•	•	•	•	•	•	•	•	•	•	•	•	•	•	•
PDS-1000/He system with Hepta™ adaptor	•	•	•	•	•	•	•	•	•	•	•	•	•	•	•	•	•	•	•	•

- First choice, highly recommended.
- Recommended, but other technologies may be better suited to that application.
- Possible, but little data to support its effectiveness.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Web site** [www.bio-rad.com](http://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