

# SAFETY ADDENDUM

## Biolistic® PDS-1000/He Particle Delivery System



**Warning:** Disconnect cord at power source before servicing.  
(refer to section 1.3)



**Warning:** Use the voltage converter only with the PDS-1000/He, do not use for other equipment. No user serviceable parts are inside. Do not attempt to remove cover.  
(refer to section 3.3)



**Warning:** Eye protection must be worn, especially during the steps for chamber evacuation, bombardment, and through venting.  
(refer to section 4.2)