



Helios Gene Gun Diffusion Screen (165-2475)

The purpose of using a screen is to reduce tissue damage in the center of the shot. Use of the screen is beneficial when bombarding attached tissue culture cells, some internal organs, and tissues that may be damaged by the high concentration of gold particles in the center of the shot. Results should be compared with and without a screen to determine whether it is beneficial in the system you are using.

Cleaning and Sterilizing

The screen may be cleaned with a soft cloth. It may be sterilized by autoclaving or by soaking in 70% or 100% ethanol.

Installation

1. Place the gene gun on a table with the barrel facing upward.
2. Place the screen over the opening of the barrel and between the spacer arms so that the support around the perimeter of the screen rests on the base of the barrel.
3. Grasp the knurled edges of the screen and turn it to lock it in place against the spacer arms. The screen has a self-centering feature so that it is positioned over the center of the barrel opening.

