

## Cleaning Log

GloMax® Discover System (Cat.# GM3000)  
 GloMax® Explorer System (Cat.# GM3500, GM3510)

Serial Number: \_\_\_\_\_

Clean the instrument as described in the *GloMax® Discover System Operating Manual #TM397* or the *GloMax® Explorer System Operating Manual #TM433* following any spill or at regular intervals as appropriate to instrument usage. Record activities below.

**Today's Date** (MM/DD/YYYY): \_\_\_\_\_ **Time:** \_\_\_\_\_

**User Name:** \_\_\_\_\_

### Instrument Maintenance

#### Instrument Cleaning:

- Remove Aperture from Aperture Plate.
- Clean the Microplate Sample Tray.
- Clean the Aperture Plate.
- Clean any residue left from various reagents.
- Remove dust using compressed air.
- Wipe clean the exterior of instrument.

#### Tablet PC Cleaning:

- To clean the Tablet PC, turn off the power.
- To avoid scratches, use a soft, lint-free cloth to gently wipe the screen. You can dampen the cloth with water or an eye-glass cleaner, but don't apply liquids directly to the Tablet PC. Don't use window cleaner or other chemical cleaners on the Tablet PC.

### Injector System Maintenance (if present)

#### After Every Use:

- The recommended flush protocol uses a sequence of 1 cycle (1ml) of deionized water, 1 cycle (1ml) of 70% ethanol, 1 cycle (1ml) of deionized water then 1 cycle (1ml) of air to flush the fluid.

#### Monthly Cleaning:

- Place the container of 70% ethanol into the bottle holder. The stainless steel inlet tube should be inserted so that the tube reaches the bottom of the bottle.
- Place an empty Waste Collection Tray on the Microplate Sample Tray.
- From the Tools screen, select *Injectors*. Tap the Prime button to initiate priming.
- Allow the solution to sit in the fluid path for 30 minutes before flushing the injectors.
- Run the Flush protocol nine times with 70% ethanol.
- Empty and clean the Waste Collection Tray, and place it on the Microplate Sample Tray.
- Place a container of deionized water on the bottle holder with the stainless steel inlet tube inserted to reach the bottom of the bottle.
- From the Tools screen, select *Injectors*. Tap the Flush button to initiate flushing.
- Select nine flush cycles.
- When all flush cycles are complete, empty and clean the Waste Collection Tray, and place it on the Microplate Sample Tray.
- Flush the injector three times with air. A small volume of water will remain in the injector after the air purge.

Comments: \_\_\_\_\_

\_\_\_\_\_

\_\_\_\_\_