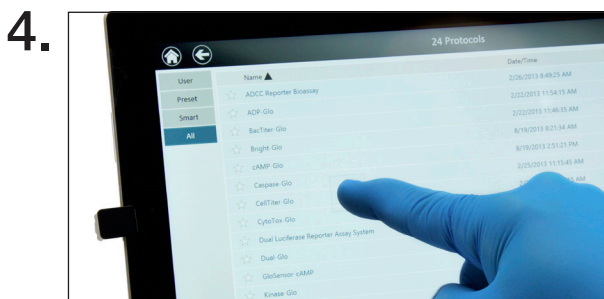
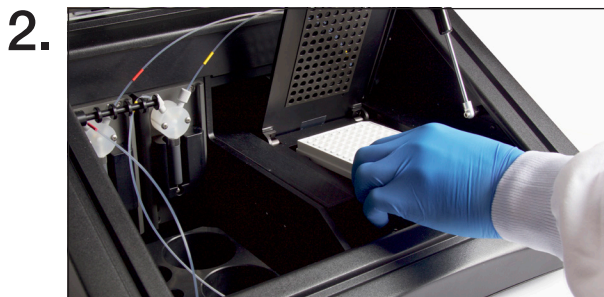
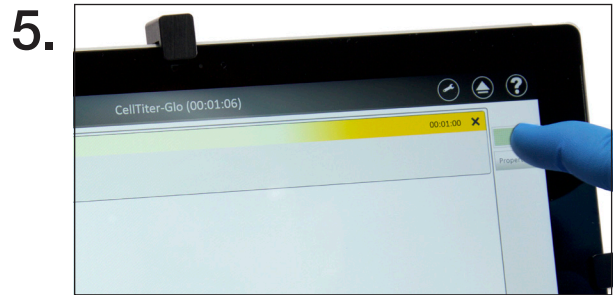
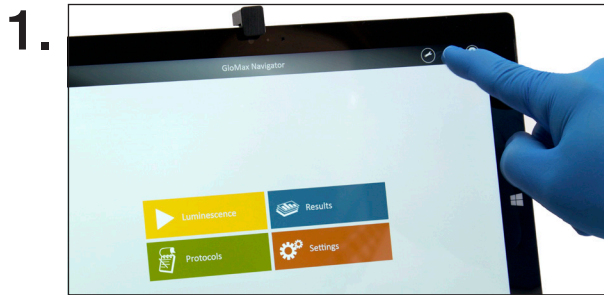


## GloMax® Navigator Detection System

### Quick Start Guide



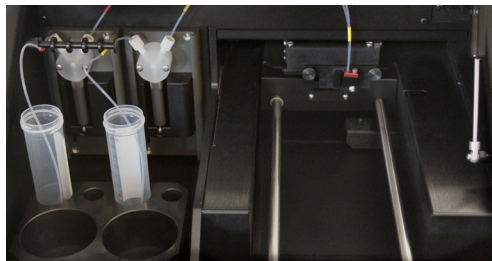
Continued on the back →



Please see the GloMax® Navigator System (TM470) Operating Manual for detailed instructions.

## GloMax® Navigator Detection System

### Cleaning Guide



1. Power OFF the instrument. Push the plate tray to the back of the instrument. Push the read head to the left to create access to the aperture.



2. Clean the Microplate Sample Tray and Aperture using a soft clean cloth dampened with 70% ethanol. Do **NOT** use other solvents or abrasive cleaners.



3. Use a soft clean cloth dampened with deionized water to periodically clean the exterior of the instrument.



4. To clean the Tablet PC, turn off the power. Use a soft clean cloth dampened with water to gently and carefully clean the touch screen.