

# Biosensor Mount Cleaning Tray for Octet HTX, RED384, and QK384 Systems

*Product Insert (Part No. 18-5133 – Pack of 12)*

## INTENDED USE

Biosensor Mount Cleaning Trays (“Cleaning Trays”) from Pall ForteBio are intended for periodic, regular, automated cleaning of biosensor mounts on Octet HTX, RED384, and QK384 instruments.

With normal instrument use, plastic residue from BLI biosensor hubs can accumulate on the tips of the metal biosensor mounts inside the instrument. This accumulation can potentially lead to incorrect loading of biosensors and inconsistencies in data sensing. It is very important to periodically remove the plastic residue in order to ensure continued optimal performance of the Octet system.

## CLEANING PROCEDURE

- 1 Remove the Cleaning Tray from plastic bag.
- 2 Spray the top of the Cleaning Tray sponge three times with 70% ethanol in a spray bottle. Do not over moisten!
- 3 Open Data Acquisition software and wait for instrument to initialize.
- 4 In the Instrument menu, select “Clean Biosensor Mounts” option. The instrument stage will present automatically.
- 5 Place the moistened Cleaning Tray in the Tray position on the instrument stage and click the “OK” button in the prompt window.
- 6 Follow the simple prompts in the software dialogue to begin the cleaning protocol. The full automated cleaning cycle will take approximately 8 minutes to complete.
- 7 Once the cleaning cycle is complete, the instrument stage will present again. Remove the used Cleaning Tray from the stage and discard.

## HOW OFTEN DO I CLEAN?

For instruments under moderate use (i.e. a few to several assays per week), running the above cleaning procedure once per month is recommended. For heavily utilized instruments (i.e. daily use for screening applications), cleaning may need to be performed as often as once per week.

Minimally, the cleaning procedure should be performed after every 100 biosensor pickups. The Data Acquisition software will display a reminder message in the lower right corner of the acquisition window to run the cleaning procedure after it has recorded 100 biosensor pickups.

## IMPORTANT NOTES

- The biosensor mount cleaning procedure is meant to be a preventative measure and will not remove excessive accumulations of plastic residue.
- If biosensor mounts are particularly dirty (indicated by visible thick white film around the tip of the mounts), OR if the instrument has been used for an extended period of time without cleaning, we recommend first scheduling a Preventative Maintenance visit with one of our service engineers before initiating regular cleanings with Cleaning Trays.
- Cleaning Trays are to be utilized for automated cleaning only with the designated method in Data Acquisition Software Version 9.0 and up.
- Cleaning Trays are recommended for single use only.

Please visit our website at <http://www.fortebio.com/support> for more information and technical support.