

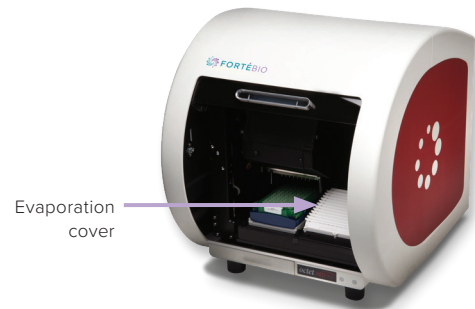
# Microplate evaporation cover

Product insert (part no. 18-5132, 3 per box)

## General information

The evaporation cover minimizes sample evaporation by only exposing one column on the sample plate at a time. Biosensors will dip into the open sample well column, and once the read is complete, the cover will return to its original position (Figure 1). The use of the evaporation cover can extend the experimental run time to up to 12 hours, with minimal sample evaporation, allowing for maximum sample recovery after concluding an assay.

- For use only on the Octet® RED96e system.
- Compatible with Greiner 96-well regular microplates (Part No. 655209).
- Extends the length of total experiment time up to 12 hours with less than 20% evaporation at 25°C.
- Single-use only and should not be cleaned or re-used.
- Can be used over a 15–40°C assay temperature range.
- Solvent resistant but should not be subjected to 100% DMSO.



- Sensitive to greater than 0.1% Triton X-100 surfactant in assay buffer. For assay buffers containing Triton X-100, reduce shake speed from 1000 RPM to 900 RPM for kinetic assays.
- Store unused evaporation covers at room temperature in a dry location in the original packaging.



Figure 1: One column flap of the evaporation cover is opened to allow for the biosensors to dip into the sample wells..

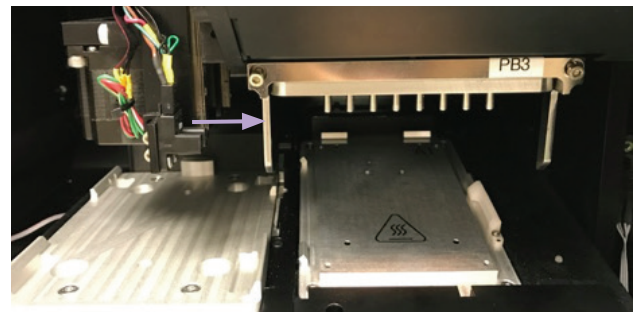
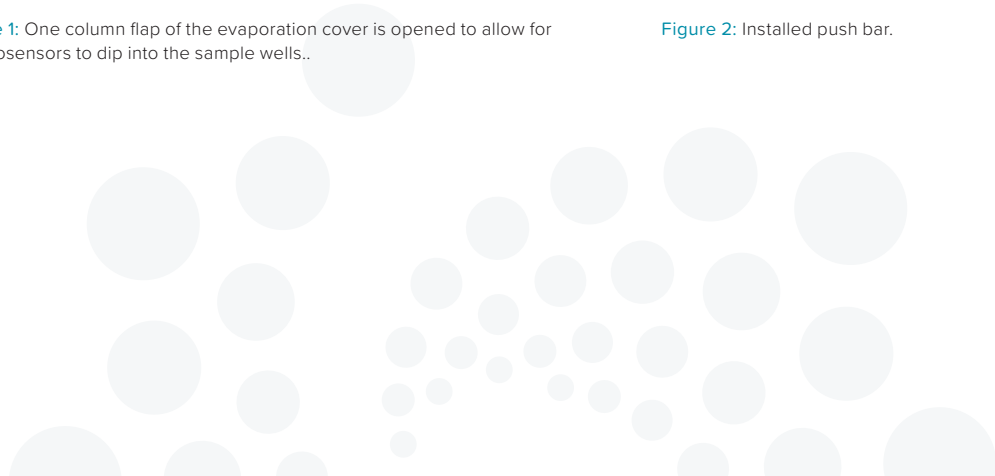


Figure 2: Installed push bar.



## Proper use of evaporation covers

- 1 Before using the evaporation cover, ensure that the push bar is installed near the sensor pickers as shown in Figure 1, otherwise the evaporation cover will not function properly.
- 2 For best results, place the 96-well microplate in the Octet RED96e instrument immediately after preparation then seat the evaporation cover on the plate.
- 3 Make sure that all four corners are pressed down onto the plate. The LED light next to the plate will illuminate and turn solid blue if the evaporation cover is installed properly (Figure 2). If the cover is not installed properly, the LED light will blink and you won't be able to start the experiment.

For your reference, the LED status information is located inside the instrument below the home position of the sensor pickers (Figure 3).

Start the experiment after placing the biosensor tray on the tray holder and giving the samples enough time to equilibrate to the desired temperature.



Figure 3: LED lights up when the evaporation cover is seated properly.

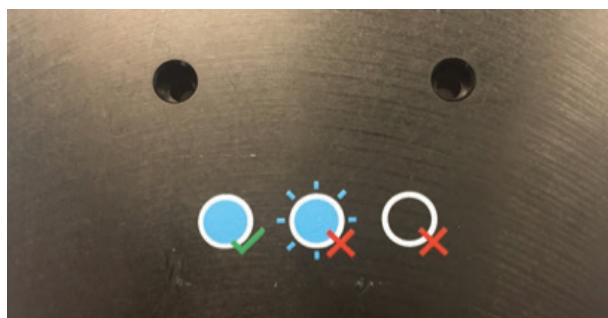


Figure 4: LED status information.