

## IonWorks HT™ Technical Tips

### Tip #1

#### **Increased seal resistances by placing PatchPlates™ in a dessicator**

A marginal increase in seal resistance performance of approximately 50-150 MΩ was observed by placing PatchPlates in a dessicator. The following observations were made at the Molecular Devices Corporation Sunnyvale, California facilities. Customers should conduct their own validation studies to verify if any appreciable benefit can be derived from the treatment.

1. Place unopened PatchPlate sleeves in a clean, empty dessicator set at 4-8% relative humidity.
2. Incubate overnight or up to 48 hours.

#### Notes:

Incubation times longer than 48 hours were detrimental, resulting in a greater number of “No seal” drop-outs.

Treatment seemed most effective during the wet season—treatment during the dry season was less effective in increasing seal resistances. Therefore, customer sites in areas with higher humidity may benefit more from desiccation treatment.