

# Premium PatchPlate Handling Instructions

---

The Premium PatchPlate™ is available for single-use in the IonWorks Barracuda™ Automated Patch Clamp System for ion channel measurements.

## Storage Precautions

- Keep the Premium PatchPlate frozen until ready to use.
- Keep the Premium PatchPlate in the factory packaging until ready to use.
- Keep the Premium PatchPlate in a vibration-free, and shock-free location.

## Use Precautions

- Do not use the Premium PatchPlate if it was dropped during handling because it may be damaged.
- Do not use the Premium PatchPlate if the low temperature cryopreserving causes plate brittleness and deterioration.
- Dispose of the used Premium PatchPlate according to your company's disposal protocol.

## Premium PatchPlate Handling Instructions

---

### Storage Temperature and Shelf Life

Item	Specifications
Storage Temperature	-70 °C or colder
Shelf Life	3 Months or less from the date on the package.  <b>Note:</b> store in factory packaging in a vibration-free and shock-free location.
Operating Temperature	Room temperature +20 °C to +25 °C  Humidity: 90% RH max.  No condensation

### Thaw Protocol

Take the package out of the freezer and leave it in the unopened packaging for 5 to 10 minutes until it reaches operating temperature.

After the Premium PatchPlate reaches room temperature, remove it from the packaging, place it in the instrument, and immediately start the measurement.

**CAUTION!** Never touch the Premium PatchPlate with bare hands.

---

**For research use only. Not for use in diagnostic procedures.**

The trademarks mentioned herein are the property of Molecular Devices, LLC or their respective owners. These trademarks may not be used in any type of promotion or advertising without the prior written permission of Molecular Devices, LLC.

Patent Pending.

Product manufactured by Molecular Devices, LLC.

1311 Orleans Drive, Sunnyvale, California, United States of America 94089.

Molecular Devices, LLC is ISO 9001 registered.

© 2011 Molecular Devices, LLC.

All rights reserved.