



Jan 21, 2021

**Subject:** Last time buy opportunities for consumables and discontinuation of repair and maintenance services for IonWorks and PatchXpress instruments

Dear Valued Customer,

We are continually streamlining our existing product portfolio. Our legacy products often rely on outdated and/unavailable third party parts, posing a challenge to providing exceptional services and support to our customers. After careful consideration, we have decided to **cease** all **repair** and **maintenance** services for our **IonWorks® Barracuda**, **IonWorks® Quattro** and **PatchXpress® 7000A** and offer last time buy opportunities for IonWorks consumables (refer to the list of consumables below).

We will continue to provide repair and maintenance services through **December 31, 2021** as long as we are able to support these activities with our current spare parts inventory. As of January 1<sup>st</sup>, 2022 all support and service for these three instruments will cease.

At this time, we are offering Last Time Buy opportunities for IonWorks consumables through **June 30, 2021**. Please note that all sales are final, we are unable to process returns or refunds. List of consumables:

Part Number	Description
5008679	Box, 48 Barracuda Patch Plate, SH
5008680	Box, 48 Barracuda Patch Plate, PPC
9000-0688	Single Hole (SH) PatchPlate for IonWorks Quattro or IonWorks HT
9000-0902	Population PatchPlate (PPC) for IonWorks Quattro or IonWorks HT

We apologize for any inconvenience this may cause. Please contact your Molecular Devices sales representative for quotes. Technical Support can be reached at 1800-635-5577 or by email [support@moldev.com](mailto:support@moldev.com).

As always, we appreciate your loyalty and remain committed to providing you with the highest quality products and services.

Thanks!

**Laxmi Chokhani**

Sustainment Product Manager