



Dear Valued Customer,

You are receiving this email because our records indicate that you have a FLIPR® Tetra High-Throughput Cellular Screening System. After careful consideration, we have decided to cease manufacturing the ICCD camera for this system. This letter serves as an Advanced Notice of product discontinuation.

Please note: This discontinuation is specific to the ICCD camera for FLIPR Tetra systems only. It does not in any way affect FLIPR systems with EMCCD and HS-EMCCD cameras.

Last day to purchase a service plan for the FLIPR Tetra system with ICCD camera: **December 31, 2021**

All service plan coverage will end by December 31, 2022, after which our ability to perform repairs and provide support will be on a reasonable effort basis.

The discontinuance of the ICCD camera for FLIPR Tetra systems is in keeping with our practice of continuous improvement and providing our users with new products that have improved design and functionality. As such, our high speed, high sensitivity EMCCD (HS-EMCCD) camera replaces the ICCD camera.

We recommend that ICCD camera customers upgrade to the HS-EMCCD camera. Visit our [website](#) for additional information and part number.

For FLIPR- ICCD Refurb camera, trade-in, or upgrading to the HS-EMCCD camera, please contact Technical Support:

- **North America:** 1.800.635.5577 or email [support@moldev.com](mailto:support@moldev.com)
- **Europe:** 00800.665.32860 or email [infoboxeu@moldev.com](mailto:infoboxeu@moldev.com)
- **China:** 400.820.3586 or email [support.china@moldev.com](mailto:support.china@moldev.com)
- For all other regions, [click here](#).

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

Thank you,

**Laxmi Chokhani**

Sustainment Product Manager