

Dear Valued Customer,

Thank you for trusting and investing in Molecular Devices for your single-cell dispensing needs in the past. The intent of this letter is to serve as an Advanced Notice of Discontinuation of products purchased through Molecular Devices and to help you with your end-of-year and early 2023 instrument purchase and support planning. The planned discontinued products encompass the Cytena brand Single-Cell Printer (SCP) instruments and disposables listed below.

Molecular Devices continuously evaluates monoclonal verification technologies to ensure that a product's quality, features, and pricing meet our customers' expectations. As a result, we will no longer be offering the Cytena brand CLONESELECT®SINGLE-CELL PRINTER™ product line after December 30th, 2022. In the coming weeks, we will be announcing a new Single-Cell Dispensing technology with options for Trade-in credits toward a new SCP technology platform. As you may know, we currently offer ClonePix 2 with monoclonality as a complete solution with reagents and media for your clone screening, picking, and monoclonality-assurance needs. Please contact your local Molecular Devices sales representative for more information on this or any of our products, including new single-cell dispensing solutions.

## Products Discontinued

### Cytena CLONESELECT®SINGLE-CELL PRINTER™ Systems

- CloneSelect™SCP (PN:5060226)
- CloneSelect Single-Cell Printer® c.sight ( PN: 5069959)
- CloneSelect Single-Cell Printer® f.sight ( PN:5069961)
- CloneSelect Single-Cell Printer® Cartridges (PN: 5060227; PN 5089025)

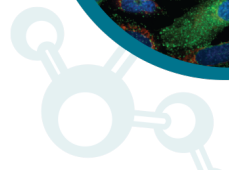
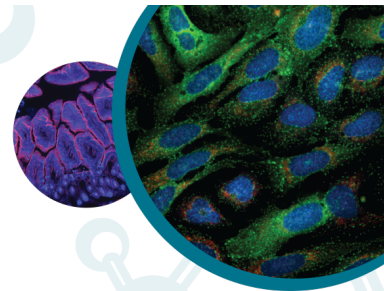
### Important dates for CLONESELECT®SINGLE-CELL PRINTER™ Systems support:

- Effective Immediately (11/23/2022): CLONESELECT®SINGLE-CELL PRINTER™ System service plan is no longer available for purchase. We will continue to support our in-warranty and extended service to customers until the end of their warranty period or service plan. Once your coverage with Molecular Devices expires, you can reach out directly to Cytena by emailing: [support@cytena.com](mailto:support@cytena.com) (United States) / [APAC@cytena.com](mailto:APAC@cytena.com) (APAC) or connect with your local Molecular Devices sales/service representatives for help coordinating support from Cytena.

3860 N First Street  
San Jose, CA 95134  
United States

toll-free 1.800.635.5577  
office 408.747.1700

[moleculardevices.com](https://moleculardevices.com)



- Effective 12/29/2022 CLONESELECT®SINGLE-CELL PRINTER™ Cartridges (PN:5089025) will no longer be available for shipping from Molecular devices. If you'd like to ensure you have enough on hand for you to last through the end of the year, please contact your local Molecular Devices sales representative by 12/15/2022. After this date, you can email [purchaseorders@cytena.com](mailto:purchaseorders@cytena.com) for a new cartridge quotation for purchase or connect with your local sales representative for help coordinating support from Cytena.
- Effective 05/31/2023 CLONESELECT®SINGLE-CELL PRINTER™ Systems will no longer be available for purchase from Molecular Devices. If you are interested in any end of the year (12/31/2022) purchases, we are selling the remaining inventory at deep discounts. Please contact your local Molecular Devices sales representative if you'd like to take advantage of that or if you need any additional information.

**Please direct your questions to any of the following methods:**

Molecular Devices North America Customer Support:

1.800.635.5577 or email [support@moldev.com](mailto:support@moldev.com)

Cytena Support:

Please email [info@cytena.com](mailto:info@cytena.com)

As always, we remain committed to supporting your research and providing you with the highest quality of products and services. Thank you for the opportunity to continue to support and help advance your research and discovery.

Please let us know if we can answer any questions.

Thank you!

3860 N First Street

toll-free 1.800.635.5577

San Jose, CA 95134

office 408.747.1700

United States

[moleculardevices.com](http://moleculardevices.com)