

April 1, 2020

Dear Valued Customer,

In an effort to optimize our portfolio, we are discontinuing our [XP Media](#) and [CloneMedia](#) for Mouse Hybridoma Generation products, effective on July 12th 2020. Parts [K8861](#), [K8862](#), [K8863](#), [K8864](#), [K8865](#), [K8866](#), [K8867](#), and [K8868](#) will no longer be available for purchase after that date.

In an effort to facilitate your transition, we are offering a one-time promotion* (subject to the terms and conditions below), for a 15% discount on the products listed above. Orders must be placed on or before May 14, 2020 with promo code **SpringBPD2020** to guarantee fulfillment. To order, please visit our [online store](#), contact your local sales rep, or call our toll free number at 1-800-635-5577. Last time delivery conditions may vary by product. Molecular Devices' final delivery date(s) may be affected by limited supply conditions.

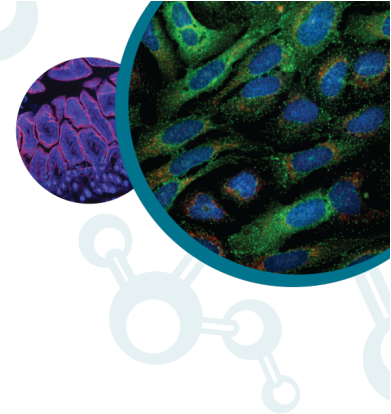
We will continue to take orders from May 15th to a final discontinuation date of July 12, 2020 as long as stock is available but cannot guarantee products will remain available during that time. Products which are out of stock prior to the final discontinuation date will no longer be offered for sale but can be purchased directly from our supplier [StemCell Technologies](#). Please refer to the comparison table provided at the end of this letter to determine equivalent part numbers. Any orders placed after July 12th will not be accepted.

All discontinued product orders subject to this notice carry Molecular Devices' standard warranty as contained in the order acknowledgement form except where noticed below. Molecular Devices will honor all warranty claims for credit only and will not provide replacement parts past the discontinuation date, or until items are out of stock. The applicable warranty period for making any return claims for discontinued products is subject to the standard [Molecular Devices policy](#).

You are a valued customer and want to thank you for your understanding. Our company is dedicated to providing you with the highest quality products and services to meet your needs.

Sincerely,

Steve Wiltgen, Ph.D.
Senior Manager, Science & Technology Strategy



Molecular Devices Product Description	Molecular Devices Part Number	StemCell Technologies Product Description	StemCell Technologies Catalog Number
XP Media and CloneMedia Complete Kit for Mouse Hybridoma Generation	K8861	ClonaCell™-HY Hybridoma Kit	#03800
XP Media Pre-Fusion Myeloma Growth Medium and Hybridoma Expansion Medium (without HT)	K8862	ClonaCell™-HY Medium A	#03801
XP Media Hybridoma Fusion Medium	K8863	ClonaCell™-HY Medium B	#03802
XP Media Hybridoma Fusion Recovery Medium	K8864	ClonaCell™-HY Medium C	#03803
CloneMedia Hybridoma Semi-Solid Selection and Cloning Medium (with HAT)	K8865	ClonaCell™ -HY Medium D	#03804
XP Media Hybridoma Growth Medium (with HT)	K8866	ClonaCell™ -HY Medium E	#03805
CloneMedia Hybridoma Semi-Solid Selection and Cloning Medium (without HAT)	K8867	ClonaCell™ -HY Medium D without HAT	#03810
Hybridoma Polyethylene Glycol (PEG) for Cell Fusion	K8868	ClonaCell™ -HY PEG	#03806

Terms and Conditions: Purchase order must reference promotion code **SpringBPD2020**. The offer is 15% off the list price of all qualifying Molecular Devices Reagents and Supplies products, **valid March 19, 2020 through May 14, 2020**. The offer does not include any other options and accessories. This promotion cannot be combined with any other promotion or discounts. No substitutions or swap outs. The discount is not convertible to cash. All prices exclude delivery charges and local tax which are the responsibility of the customer. Purchase orders must be received **by 5:00 PM, Pacific Standard Time on May 14, 2020**. Promotion is not valid for standing orders. The purchase is subject to Molecular Devices standard terms and conditions of sale. Promotion is available in regions where product can be sold, when buying directly from Molecular Devices. All products are research use only. Void where prohibited.