

QPix[™] Chroma Filter

Thin Film Optical Filter for
Blue/White Colony Selection

The QPix™ Chroma Filter is a thin film translucent optical filter that provides a robust method to confidently select and pick rare expressing bacterial colonies based on color intensity in white light. The optical properties of the filter have been selected and extensively tested for the ability to distinguish between blue and white colonies, even when the blue colonies are newly formed and powder blue.

Description	Part number
QPix™ Chroma Filter Thin film optical filter for Blue/White colony selection, 1 filter	X1101
QPix™ Chroma Filters Thin film optical filters for Blue/White colony selection, 25 filters	X1102
QPix™ Chroma Colorimetric Colony Selection Software Kit. Blue/White Colony Selection Software Kit for QPix 400 Series. Includes software license and 2 QPix™ Chroma Filters (X1101)	X1103

Table 1-1. Product configurations

Contents

	Page
Chapter 1: About the QPix™ Chroma Filter	4
Product Specifications	4
Blue/White Colony Detection Software Module	4
Storage and Handling	4
Chapter 2: Ordering Information	5
Chapter 3: Obtaining Support	6



This document is provided to customers who have purchased Molecular Devices equipment, software, reagents, and consumables to use in the operation of such Molecular Devices equipment, software, reagents, and consumables. This document is copyright protected and any reproduction of this document, in whole or any part, is strictly prohibited, except as Molecular Devices may authorize in writing.

Software that may be described in this document is furnished under a non-transferrable license. It is against the law to copy, modify, or distribute the software on any medium, except as specifically allowed in the license agreement. Furthermore, the license agreement may prohibit the software from being disassembled, reverse engineered, or decompiled for any purpose.

Portions of this document may make reference to other manufacturers and/or their products, which may contain parts whose names are registered as trademarks and/or function as trademarks of their respective owners. Any such usage is intended only to designate those manufacturers' products as supplied by Molecular Devices for incorporation into its equipment and does not imply any right and/or license to use or permit others to use such manufacturers' and/or their product names as trademarks.

Each product is shipped with documentation stating specifications and other technical information. Molecular Devices products are warranted to meet the stated specifications. Molecular Devices makes no other warranties or representations express or implied, including but not limited to, the fitness of this product for any particular purpose and assumes no responsibility or contingent liability, including indirect or consequential damages, for any use to which the purchaser may put the equipment described herein, or for any adverse circumstances arising therefrom. The sole obligation of Molecular Devices and the customer's sole remedy are limited to repair or replacement of the product in the event that the product fails to do as warranted.

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.

Patents: <http://www.moleculardevices.com/productpatents>

Chapter 1: About the QPix™ Chroma Filter

The QPix™ Chroma Filter is a thin film translucent optical filter that provides a robust method to confidently select and pick rare expressing bacterial colonies based on color intensity in white light. The optical properties of the filter have been selected and extensively tested for the ability to distinguish between blue or white colonies, even when the blue colonies are newly formed and powder blue.

The QPix™ Chroma Filter design has the same perimeter as either a source plate Qtray, or any of the adjustable holders containing source plate receptacles. The filter easily places on or removes from the imaging bed of the instrument to fit between the light source and specimen tray. The optical properties of the filter are selected and extensively tested for the ability to distinguish between blue and white colonies.

The QPix™ Chroma Filter only works in conjunction with the QPix Software version 2.0 or later, and QPix™ Chroma Colorimetric Colony Selection Software License: Blue/White Colony Selection Software License for QPix 400 Series.

Product Specifications

Dimensions: 23 cm x 23 cm

Thickness: 1 mm

Assembly: A heavy, black, bond, paper border frame around the filter acts as a handling stiffener and prevents light scattering.

Note: Due to inherent biological variations, some picked white colonies lose their insertions and might later be viewed to be blue. These effects are inherent to the biology and are not due to the picking or selection processes. When selecting the insertion methodology, take care to make sure that the desired results are achieved.

Blue/White Colony Detection Software Module

The QPix Software Version 2.0 offers both regular and regional picking for blue or white bacterial colonies only with the installation of the optional QPix™ Chroma Colorimetric Colony Selection Software License: Blue/White Colony Selection Software License for QPix 400 Series.

When stepping through the Blue/White Picking process, the Test Image screen is where you visualize and select either blue or white colonies. The detection controls enable the manual or automatic selection of blue or white colonies. See the *QPix Colony Picking Software User Guide* for your instrument for more details.

Storage and Handling

The QPix™ Chroma Filter ships at room temperature in a rigid envelope. The rigid envelope is also intended for filter storage when not in use. Carefully handle the filter, touching only the black frame of the filter.

To prevent formation of slight creases and dents in the film, which can cause distortions in the optical path, store the filters in the rigid envelopes when not in use.

The filter is validated for up to ten use cycles with proper handling. To log the number of filter uses, place a mark in a usage tracking check box provided on the frame edge of the filter.

Using a soft, lint-free cloth, gently clean the filter with ethyl or isopropyl alcohol, as needed. It can also be sterilized under the QPix UV sterilization cycle.

Chapter 2: Ordering Information

The following software and consumable are recommended to operate the QPix instrument with Software Version 2.0 or later.

Description	Part number
Software	
QPix™ Chroma Colorimetric Colony Selection Software License. Blue/White Colony Selection Software License for QPix 420, 450 or 460 Colony Picking Systems	SL9400-A06
Zone of Inhibition Detection Software License for QPix 420, 450 or 460 Colony Picking Systems	SL9400-A07
QPix 420 Software Version 2.0 Upgrade	SL9400-A08
QPix 450/460 Software Version 2.0 Upgrade	SL9400-A09
Consumables	
QPix™ Chroma Filter, thin film optical filter for Blue/White colony selection, qty 1	X1101
QPix™ Chroma Filter, thin film optical filter for Blue/White colony selection, qty 25	X1102
QPix™ Chroma Colorimetric Colony Selection Software Kit. Blue/White Colony Selection Software Kit for QPix 420, 450 or 460 Colony Picking Systems. Includes software license and 2 QPix™ Chroma Filters (X1101).	X1103
1 Way Adjustable Petri Dish Holder Plate diameter compatibility: 138 mm to 141 mm	X9401
2 Way Adjustable Microplate Holder Plate diameter compatibility: Any ANSI/SLAS 1-2004 microplates	X9402
4 Way Adjustable Petri Dish Holder Plate diameter compatibility: 97 mm to 100.5 mm	X9403
5 Way Adjustable Petri Dish Holder Plate diameter compatibility: 87 mm to 90.5 mm	X9404

Table 2-1. Available products



For more information about QPix™ instruments and accessories visit:
www.moleculardevices.com/qpix-sw2.0

Chapter 3: Obtaining Support

Molecular Devices is a leading worldwide manufacturer and distributor of analytical instrumentation, software and reagents. We are committed to the quality of our products and to fully supporting our customers with the highest possible level of technical service.

Our support web site, www.moleculardevices.com/support, has a link to the Knowledge Base with technical notes, software upgrades, safety data sheets, and other resources. If you do not find the answers you are seeking, follow the links to the Technical Support Service Request Form to send an email message to a pool of technical support representatives.

You can contact your local representative or contact Molecular Devices Technical Support by telephone at 800-635-5577 (U.S. only) or +1 408-747-1700. In Europe call +44 (0) 118 944 8000.

Please have the product name, part number, and lot number available when you call.

Contact Us

Phone: [+1.800.635.5577](tel:+18006355577)
 Web: www.moleculardevices.com
 Email: info@moldev.com

Check our website for a current listing of worldwide distributors.

Regional Offices

USA and Canada	+1.800.635.5577	China (Beijing)	+86.10.6410.8669	Japan	+81.3.6362.9109
United Kingdom	+44.118.944.8000	China (Shanghai)	+86.21.3372.1088	South Korea	+82.2.3471.9531
Europe*	00800.665.32860	Hong Kong	+852.3971.3530		

**Austria, Belgium, Denmark, Finland, France, Germany, Ireland, Netherlands, Spain, Sweden and Switzerland*