

96-Well Spectral Calibration Plates

Catalog Numbers 4432327, 4432334, 4432340, 4432346, 4432358, 44432370, 4432352

Pub. No. 4432395 **Rev.** C

Cat. no.	Description	Storage conditions
4432327	96-Well Spectral Calibration Plate with FAM™ dye	Store at –15 to –25°C.
4432334	96-Well Spectral Calibration Plate with VIC® dye	
4432340	96-Well Spectral Calibration Plate with ROX™ dye	
4432346	96-Well Spectral Calibration Plate with SYBR® Green dye	
4432352	96-Well Spectral Calibration Plate with TAMRA™ dye	
4432358	96-Well Spectral Calibration Plate with NED™ dye	
4432370	96-Well Spectral Calibration Plate with FAM™/ROX™ and VIC®/ ROX™ dyes]	

Note: For safety and biohazard guidelines, refer to the “Safety” section in the appropriate user guide for your instrument. For every chemical, read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Description

The 96-Well Spectral Calibration Plates are used for spectral calibration and normalization of Applied Biosystems® Real-Time PCR instruments. Perform calibration during installation and as part of routine instrument maintenance.

For information on instrument setup, the calibration process, and data analysis, refer to the appropriate user guide for your instrument.

For more information, visit our web site: www.lifetechnologies.com

Limited product warranty

Life Technologies Corporation and/or its affiliate(s) warrant their products as set forth in the Life Technologies' General Terms and Conditions of Sale found on Life Technologies' website at www.lifetechnologies.com/termsandconditions. If you have any questions, please contact Life Technologies at www.lifetechnologies.com/support.

The information in this guide is subject to change without notice.

DISCLAIMER

TO THE EXTENT ALLOWED BY LAW, LIFE TECHNOLOGIES AND/OR ITS AFFILIATE(S) WILL NOT BE LIABLE FOR SPECIAL, INCIDENTAL, INDIRECT, PUNITIVE, MULTIPLE, OR CONSEQUENTIAL DAMAGES IN CONNECTION WITH OR ARISING FROM THIS DOCUMENT, INCLUDING YOUR USE OF IT.

Important Licensing Information: These products may be covered by one or more Limited Use Label Licenses. By use of these products, you accept the terms and conditions of all applicable Limited Use Label Licenses.

©2015 Thermo Fisher Scientific Inc. All rights reserved. All trademarks are the property of Thermo Fisher Scientific and its subsidiaries unless otherwise specified.

For support visit lifetechnologies.com/support or email techsupport@lifetech.com
lifetechnologies.com

