

96-Well 0.1-mL Spectral Calibration Plates

Catalog Numbers A26336, A26337, A26340

Pub. No. 100025903 **Rev.** A

Cat. no.	Description	Quantity	Storage conditions
A26336	96-Well 0.1-mL Spectral Calibration Plate 1 (containing FAM™, VIC®, ROX™, and SYBR® dyes)	1 plate	Store at -15 to -25°C.
A26337	96-Well 0.1-mL Spectral Calibration Plate 2 (containing ABY®, JUN®, and MUSTANG PURPLE® dyes)	1 plate	
A26340	96-Well 0.1-mL Spectral Calibration Plate 3 (containing TAMRA™, NED™, and Cy®5 dyes)	1 plate	

Note: For Safety and biohazard guidelines, refer to the "Safety" appendix in the appropriate Real-Time PCR Systems User Guide for your instrument. Read the Safety Data Sheets (SDSs) and follow the handling instructions. Wear appropriate protective eyewear, clothing, and gloves.

Description

The 96-Well 0.1-mL Spectral Calibration Plates are used for spectral normalization of Real-Time PCR Systems. Perform calibration as part of routine system maintenance as outlined in the appropriate Real-Time PCR Systems User Guide.

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. Cy is a registered trademark of GE Healthcare.

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

lifetechnologies.com

17 March 2015