

# Spectral Calibration Plates with ABY<sup>®</sup> or JUN<sup>®</sup> Dye for Multiplexing

Pub. Part no. 100022948 Rev. A Rev. Date January 2014

## 96-well plates

Part Number	Part	Storage Conditions
A24737	96-Well Spectral Calibration Plate with $JUN^{\textcircled{B}}$ Dye for Multiplex qPCR	Store at -15°C to -25°C
A24738	96-Well Spectral Calibration Plate with $ABY^{I\!\!R}$ Dye for Multiplex qPCR	

## Fast 96-well plates

Part Number	Part	Storage Conditions
A24735	Fast 96-Well Spectral Calibration Plate with $JUN^{\textcircled{B}}$ Dye for Multiplex qPCR	Store at -15°C to -25°C
A24734	Fast 96-Well Spectral Calibration Plate with $ABY^{\textcircled{B}}$ Dye for Multiplex qPCR	

## 384-well plates

Part Number	Part	Storage Conditions
A24733	384-Well Spectral Calibration Plate with $JUN^{\textcircled{R}}$ Dye for Multiplex qPCR	Store at -15°C to -25°C
A24736	384-Well Spectral Calibration Plate with $ABY^{\mathbb{R}}$ Dye for Multiplex qPCR	

**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 Spectral Calibration Plates with ABY<sup>®</sup> or JUN<sup>®</sup> Dye for Multiplexing are used for spectral calibration of Applied Biosystems<sup>®</sup> 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: http://www.lifetechnologies.com

## For Research Use Only. Not for use in diagnostic procedures.

## Important Licensing Information

This product may be covered by one or more Limited Use Label Licenses. By use of this product, you accept the terms and conditions of all applicable Limited Use Label Licenses.

## 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.

© 2014 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.



