

7 Dye qPCR Calibration Plate *Optimized for ABI7500 Fast 96-Well*

Catalog number: 67020
Unit size: 1 Set

Component	Storage	Amount (Cat No. 67020)
Component A: ROX Dye qPCR Calibration Plate	Freeze (< -15 °C), Minimize light exposure	1 Plate
Component B: SYBR Dye qPCR Calibration Plate	Freeze (< -15 °C), Minimize light exposure	1 Plate
Component C: TAMRA Dye qPCR Calibration Plate	Freeze (< -15 °C), Minimize light exposure	1 Plate
Component D: NED Dye qPCR Calibration Plate	Freeze (< -15 °C), Minimize light exposure	1 Plate
Component E: FAM Dye qPCR Calibration Plate	Freeze (< -15 °C), Minimize light exposure	1 Plate
Component F: VIC Dye qPCR Calibration Plate	Freeze (< -15 °C), Minimize light exposure	1 Plate
Component G: JOE Dye qPCR Calibration Plate	Freeze (< -15 °C), Minimize light exposure	1 Plate

OVERVIEW

7 Dye qPCR Calibration Plate contains seven spectral calibration plates with seven separate dye standards (TAMRA, SYBR, FAM, JOE, NED, ROX, and VIC). It can be used to maintain your 7500 Real-Time PCR system with Fast 96-well block. For most of qPCR instruments, the necessary calibrations should be run at least every six months. This calibration plate is ready to use without any additional preparation steps required. The qPCR calibration plate might significantly improve qPCR results with multiplexing by more accurately representing fluorescent spectra used in your real-time experiments. Please refer to your instrument's guide for the detailed calibration operation.

Figure 1. ABI 7500 FAST system spectra for seven different dye standards (TAMRA, VIC, ROX, FAM, SYBR, NED, and JOE).

DISCLAIMER

AAT Bioquest provides high-quality reagents and materials for research use only. For proper handling of potentially hazardous chemicals, please consult the Safety Data Sheet (SDS) provided for the product. Chemical analysis and/or reverse engineering of any kit or its components is strictly prohibited without written permission from AAT Bioquest. Please call 408-733-1055 or email info@aatbio.com if you have any questions.

SAMPLE EXPERIMENTAL PROTOCOL

1. Obtain the calibration plate from the spectral calibration kit in the freezer.
2. Allow the calibration plate to warm to room temperature (approximately 5 min).
Note: Do not remove calibration plate from its packaging until you are ready to run it. The fluorescent dye in the wells of the plate is photosensitive. Prolonged exposure to light can diminish the fluorescence from the plate.
3. Remove the calibration plate from its packaging. Leave the optical film on the plate.
4. Centrifuge the plate for 2 min at less than 1500 rpm.
5. Verify that the liquid in each well of the calibration plate is at the bottom of the well. If not, centrifuge the plate again at a higher rpm and for a longer period of time.
6. Follow the instructions of the instrument to run the calibration plate.

EXAMPLE DATA ANALYSIS AND FIGURES

