

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

Catalog Number: 67017

Unit Size: 1 Plate

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Liquid
Molecular Weight	N/A
Soluble In	DMSO

Spectral Properties

Excitation Wavelength	559 nm
Emission Wavelength	570 nm

Applications

The ABY Dye qPCR Calibration Plate is designed for use with the 7500 Real-Time PCR system featuring a Fast 96-well block. It facilitates the calibration and maintenance of your qPCR instruments. To ensure optimal performance, it is recommended to conduct these calibrations biannually. This ready-to-use calibration plate requires no further preparation, streamlining the calibration process. Utilizing this plate may significantly enhance the accuracy of fluorescent spectra representation in multiplex qPCR experiments, leading to improved results. For detailed calibration procedures, please consult the user manual of your specific instrument.