

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

Catalog number: 67012 Unit size: 1 Plate

Product Details	
Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving
Chemical Properties Appearance	Liquid

Applications

JOE Dye qPCR Calibration Plate 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.

Note: After calibration, JOE dye qPCR calibration plates can be stored at -20 °C, protected from light. Avoid freeze-thawing plates. It is not recommended to use the plates more than 3 times.