

Cy3.5 Dye qPCR Calibration Plate
Optimized for ABI7500 Fast 96-WellCatalog number: 67014
Unit size: 1 Plate**Product Details**

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

Chemical Properties

Appearance	Liquid
Soluble In	DMSO

Applications

Cy3.5 Dye qPCR Calibration Plate can be used to maintain your 7500 Real-Time PCR System with a Fast 96-well block. For most qPCR instruments, it is necessary to calibrate the device at least every six months. This calibration plate is ready to use without any additional preparation steps. 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.