

Cy5 Dye qPCR Calibration Plate *Optimized for ABI7500 Fast 96-Well*

Catalog Number: 67016

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	N/A

Spectral Properties

Excitation Wavelength	651 nm
Emission Wavelength	670 nm

Applications

A Cy5 dye qPCR calibration plate is a type of plate used in quantitative PCR (polymerase chain reaction) experiments. It is designed to calibrate the real-time PCR instrument, which measures the fluorescence signal emitted by the Cy5 dye during the PCR reaction. The Cy5 dye qPCR calibration plate contains a set of pre-determined concentrations of a DNA template that serves as a reference for the quantification of unknown samples. The plate is loaded into the PCR instrument along with the unknown samples, and the instrument measures the fluorescence signal from each well during the PCR amplification process. By comparing the fluorescence signal of the unknown samples to the known concentrations in the calibration plate, researchers can determine the amount of DNA present in the unknown samples. This information can be used to quantify gene expression levels, detect the presence of pathogens, or perform other types of molecular analysis.