

PE-iFluor™ 770 Tandem

Catalog Number: 2608

Unit Size: 1 mg

Product Details

Storage Conditions	Refrigerated (2-8 °C), Minimize light exposure,
Expiration Date	6 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	N/A
Soluble In	Water

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

PE-iFluor™ 770 Tandem is a new color used in flow cytometry. Its primary absorption peak is at 565 nm with emission peak at ~800 nm. It has been validated with a spectral flow cytometer. PE-iFluor™ 770 Tandem gives a significantly improved staining index than the other spectrally similar PE-IR dye tandems. It is optimized to provide a new unique color for spectral flow cytometry. AAT Bioquest offer the largest number of colors for conventional and spectral flow cytometry applications, including iFluor™, mFluor™ small organic dyes and a variety of their tandems.