

PE-iFluor™ 780 TandemCatalog number: 2617
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

Applications

PE-iFluor™ 780 Tandem is a new NIR 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. It is optimized to provide a new unique color for spectral flow cytometry. AAT Bioquest offers the largest number of colors for conventional and spectral flow cytometry applications, including iFluor™, mFluor™ small organic dyes and a variety of their tandems.