

APC-iFluor® 780 Tandem

Catalog Number: 2639

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	787 nm
Emission Wavelength	812 nm

Applications

Tandem dyes are a unique class of fluorescent molecules that consist of two different covalently linked fluorophores, a donor (e.g. PE or APC) and a longer-wavelength emitting fluorescence acceptor (e.g. Texas Red, Cy5, Cy7, iFluor® 594 or iFluor® 750). APC-iFluor® 780 is a new unique tandem color for multicolor applications, in particular, for designing spectral flow cytometric panels. It has been validated with Cytex's Aurora flow cytometer. Its primary absorption peak is at 651 nm with emission peak at ~804 nm. It is one of the brightest NIR fluorophores that can be used with flow cytometry and fluorescence imaging applications.