

APC-XFD750 Tandem

Catalog number: 2627

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
Soluble In	Water

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-XFD750 tandem has the structure and properties to the commonly used APC-Alexa Fluor® 750. Its primary absorption peak is at 651 nm with emission peak at ~780 nm.