

APC-iFluor™ 720 Tandem

Catalog Number: 2633

Unit Size: 1 mg

۲	r	0	a	u	C	τ	υ	e	τ	a	I	IS	>

Storage Conditions Refrigerated (2-8 °C), Minimize light exposure,

Expiration Date 6 months upon recieving

Chemical Properties

Appearance Solid

Molecular Weight N/A

Soluble In Water

Spectral Properties

Excitation Wavelength N/A

Emission Wavelength N/A

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™ 720 is a new unique tandem color for multicolor applications, in particular, for designing spectral flow cytometric panels. It has been validated with Cytek's Aurora flow cytometer. Its primary absorption peak is at 651 nm with emission peak at ~740 nm. It is one of the brightest NIR fluorophores that can be used with flow cytometry and fluorescence imaging applications.