

PerCP-iFluor® 780

Catalog Number: 2655

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	480 nm
Emission Wavelength	806 nm

Applications

PerCP (Peridinin-chlorophyll-protein complex) is isolated from Dinophyceae sp. It has an extremely high extinction coefficient, a high quantum efficiency and a large Stokes shift. It is well excited with the Argon laser at 488 nm with its maximum emission peak at 677 nm. PerCP protein is commonly used for fluorescent immunolabeling, particularly in applications involving fluorescent-activated cell sorting (FACS). PerCP-iFluor® 780 tandem conjugates can be excited with a standard 488 nm laser and emits in the far red at a longer wavelength for multicolor flow cytometric analysis of cells. These multiple emission wavelengths make PerCP-iFluor® 780 conjugates potentially useful fluorochromes for multicolor analysis with FITC, PE and other fluorochromes. This PerCP tandem might have the longest emission among all the PerCP tandems.