

PerCP-iFluor® 750

Catalog Number: 2656

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	482 nm
Emission Wavelength	784 nm

Applications

PerCP-iFluor® 750 is a tandem fluorophore engineered for high-parameter flow cytometry, providing strong far-red fluorescence with minimal spectral spillover. It comprises a peridinin chlorophyll protein (PerCP) donor covalently linked to an iFluor® 750 acceptor, allowing efficient energy transfer through Förster resonance energy transfer (FRET). Upon excitation with a 488 nm blue laser, the conjugate emits maximally near 784 nm, yielding a large Stokes shift and enabling sensitive detection in the near-infrared region. Its distinct spectral characteristics make PerCP-iFluor® 750 highly suitable for use in complex multicolor panels, where it helps minimize compensation and improves resolution of closely spaced populations. The dye is ideal for applications such as immunophenotyping, rare cell analysis, and other advanced flow cytometry workflows requiring high signal-to-noise ratios and precise population discrimination. PerCP-iFluor® 750 can be conjugated to proteins and antibodies using SMCC chemistry or through site-specific conjugation with Buccutite™ MTA and Buccutite™ FOL.