

PerCP-iFluor® 720

Catalog Number: 2653

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	747 nm

Applications

PerCP-iFluor® 720 is a tandem fluorochrome specifically engineered for high-parameter flow cytometry applications. It consists of a peridinin chlorophyll protein (PerCP) donor covalently linked to an iFluor® 720 acceptor, enabling efficient energy transfer via Förster resonance energy transfer (FRET) upon excitation with a 488 nm blue laser. This results in a strong and well-defined emission peak near 747 nm, making it an ideal far-red reporter in multicolor panels.

The spectral characteristics of PerCP-iFluor® 720 allow for minimal spillover into adjacent channels, such as PE-Cy7 and APC, facilitating cleaner resolution in complex staining panels. Compared to traditional PerCP tandems, PerCP-iFluor® 720 exhibits enhanced photostability and higher signal-to-noise ratios, improving sensitivity for the detection of low-abundance targets. Its consistent performance across fixation and permeabilization conditions makes it suitable for both surface and intracellular staining protocols. Conjugation to antibodies or streptavidin further expands its utility in immunophenotyping, cell subset discrimination, and high-dimensional cytometry.