

## PerCP-Cy7

Catalog Number: 2654

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	780 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-Cy7 tandem conjugate is a common color used in flow cytometry. It 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-Cy7 conjugates an extremely useful fluorochrome for multicolor analysis with FITC, PE and other fluorochromes.