

## PerCP-iFluor® 710

Catalog Number: 2652

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 (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® 710 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® 710 conjugates potentially useful fluorochromes for multicolor analysis with FITC, PE and other fluorochromes. PerCP tandem structure may make it more photostable than PerCP alone, which generally photobleaches rapidly with more powerful water-cooled gas lasers.