

**ReadiUse™ Preactivated PerCP-Cy7
Maleimide**Catalog number: 2725
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

Applications

PerCP-Cy7 is a popular tandem color used in flow cytometry. Its primary absorption peak is at ~480 nm with emission peak at ~780 nm. AAT Bioquest offers this preactivated PerCP-Cy7 to facilitate the PerCP-Cy7 tandem conjugations to reduced antibodies and other biomolecules that contain a thiol group. Our preactivated PerCP-Cy7 maleimide is prepared from the commonly used crosslinker SMCC, and ready to conjugate. 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. Its tandem conjugates (such as PerCP-Cy7) 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- Cyanine 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.