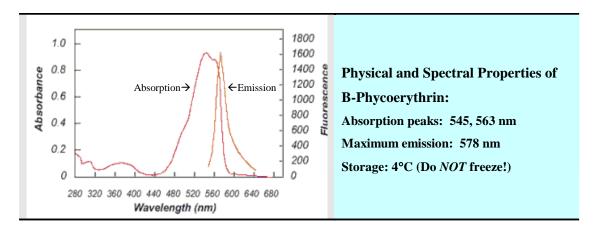
## **SMCC Activated B-Phycoerythrin**



Catalog Number: AS- 72109 Concentration: 1mg/mL Size: 1 mg

**Description:** B-PE (B-Phycoerythrin), <sup>1</sup> a fluorescent protein, belongs to the phycobiliprotein family of highly soluble and fluorescent proteins derived from cyanobacteria and eukaryotic algae. B-PE is made of  $\alpha$ ,  $\beta$  and  $\gamma$  subunits and is present as  $(\alpha\beta)_6\gamma$ . The absorption bands have peaks at 545 nm ( $\epsilon_M$  =2.41 X 10<sup>6</sup> M<sup>-1</sup>cm<sup>-1</sup>) and 563 ( $\epsilon_M$  =2.33 X 10<sup>6</sup> M<sup>-1</sup>cm<sup>-1</sup>). <sup>1,2</sup> B-PE and the closely related R-PE are the most intensely fluorescent phycobiliproteins. They are significantly brighter and more photostable than conventional organic fluorophores. <sup>3</sup> B-PE labeled streptavidins, primary and secondary antibodies have been widely used in applications such as flow cytometry and multi-color immunofluorescent staining.

SMCC Activated B-PE, with a maleimide group introduced, allows it to be conveniently conjugated to the thiol groups of proteins, without the need for additional activation.

SMCC Activated B-PE is supplied in MES buffer, pH 6.0 with EDTA and preservative. Store at 4°C and keep from light.

## References

- 1. Glazer, AN. and CS. Hixson J. Biol. Chem. 252, 32 (1977).
- 2. Lundell, DJ. et al. J. Biol. Chem. 259, 5472 (1984).
- 3. Oi, VT. et al. J. Cell Biol. 93, 981 (1982).