

**ReadiUse™ PE [R-Phycoerythrin] \*Ammonium Sulfate-Free\***

Catalog Number: 2500, 2501

Unit Size: 1 mg, 10 mg

**Product Details**

---

Storage Conditions	Refrigerated (2-8 °C), Minimize light exposure
--------------------	--

Expiration Date	12 months upon receiving
-----------------	--------------------------

**Chemical Properties**

---

Appearance	Powder red
------------	------------

Molecular Weight	~240000
------------------	---------

Soluble In	Water (Add 100 µL of H <sub>2</sub> O to Cat# 2500 or 1 mL to Cat# 2501 to reconstitute to 10 mg/mL in PBS)
------------	---

**Spectral Properties**

---

Excitation Wavelength	565 nm
-----------------------	--------

Emission Wavelength	574 nm
---------------------	--------

**Applications**

---

R-Phycoerythrin (PE) is isolated from red algae. Its primary absorption peak is at 565 nm with secondary peaks at 496 and 545 nm. All the commercial PE materials are sold in concentrated ammonium sulfate buffers. The commercial PE materials from other vendors require the tedious dialysis or other purifications performed before it can be used for labeling purposes. AAT Bioquest offers this ReadiUse™ PE that can be readily used for any labelings without any purifications required. Our highly purified ReadiUse™ PE facilitates the rapid PE conjugations to antibodies and other proteins such as streptavidin and other secondary reagents.