

---

## Product Information Sheet

### **Ordering Information**

Product Number: 2500

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

Unit Size: 1 mg

Storage Conditions: Refrigerated (2-4 °C), Desiccated, Avoid Light

Expiration Date: 12 months upon receiving

### **Chemical and Spectral Properties**

Appearance: Solid

Molecular Weight: N/A

Soluble In: Water

Excitation Wavelength: 565

Emission Wavelength: 575

### **Application Notes**

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.