

# 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)

Expiration Date 12 months upon receiving

## **Chemical Properties**

Appearance Solid

Molecular Weight ~240000

Soluble In Water

## **Spectral Properties**

Excitation Wavelength 566 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.