

---

## Product Information Sheet

### Ordering Information

Product Number: 2616  
Product Name: PE-Cy7 Tandem  
Unit Size: 1 mg  
Storage Conditions: Refrigerated (2-4 °C), Desiccated, Avoid Light  
Expiration Date: 6 months upon receiving

### Chemical and Spectral Properties

Appearance: Solid  
Molecular Weight: N/A  
Soluble In: Water  
Excitation Wavelength: 565  
Emission Wavelength: 780

### Application Notes

PE-Cy7 is a popular color used in flow cytometry. Its primary absorption peak is at 565 nm with emission peak at 780 nm. AAT Bioquest also offers a unique preactivated PE-Cy7 to facilitate the PE-Cy7 tandem conjugations to antibodies and other proteins such as streptavidin and other secondary reagents. Our preactivated PE-Cy7 tandem is ready to conjugate, giving much higher yield than the conventionally tedious SMCC-based conjugation chemistry. In addition, our preactivated PE-Cy7 tandem is conjugated to a protein via its amino group that is abundant in proteins while SMCC chemistry targets the thiol group that has to be regenerated by the reduction of antibodies.