

---

## Product Information Sheet

### **Ordering Information**

Product Number: 2613  
Product Name: PE-Cy5.5 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: 700

### **Application Notes**

PE-Cy5.5 is a popular color used in flow cytometry. Its primary absorption peak is at 565 nm with emission peak at ~700 nm. The filter sets of 682/33 nm and 695/40 nm are recommended for this tandem color. AAT Bioquest also offers a unique preactivated PE-Cy5.5 to facilitate the PE-Cy5.5 tandem conjugations to antibodies and other proteins such as streptavidin and other secondary reagents. Our preactivated PE-Cy5.5 tandem is ready to conjugate, giving much higher yield than the conventionally tedious SMCC-based conjugation chemistry. In addition, our preactivated PE-Cy5.5 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.