

PE-Texas Red Tandem

Catalog number: 2619

Unit size: 1 mg

Product Details

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

Expiration Date 6 months upon receiving

Chemical Properties

Appearance Purple solid

Molecular Weight ~240000

Soluble In Water

Spectral Properties

Excitation Wavelength 567 nm

Emission Wavelength 615 nm

Applications

PE-Texas Red is a popular color used in flow cytometry. Its primary absorption peak is at 565 nm with emission peak at 600 nm. AAT Bioquest also offers a unique preactivated PE-Texas Red to facilitate the PE-Texas Red tandem conjugations to antibodies and other proteins such as streptavidin and other secondary reagents. Our preactivated PE-Texas Red tandem is ready to conjugate, giving much higher yield than the conventionally tedious SMCC-based conjugation chemistry. In addition, our preactivated PE-Texas Red 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.