

Texas Red® hydrazide *Single Isomer*

Catalog Number: 481

Unit Size: 5 mg

Product Details

Storage Conditions Freeze (< -15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

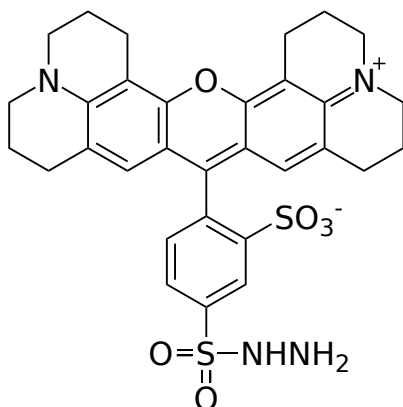
Chemical Properties

Appearance Solid purple

Molecular Weight 620.74

Soluble In DMSO

Chemical Structure

**Spectral Properties**

Excitation Wavelength 586 nm

Emission Wavelength 603 nm

Applications

Texas Red® hydrazide reacts with carbonyl compounds such as aldehyde, ketones and acids to give bright red fluorescent conjugates. It is widely used for labeling glycoproteins and carbohydrates through reductive amination.