

# Texas Red® cadaverine \*Single Isomer\*

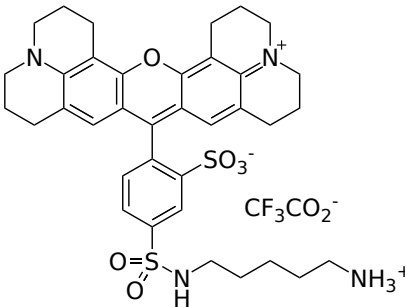
Catalog Number: 482

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	804.89
Soluble In	DMSO
Chemical Structure	

## Spectral Properties

Excitation Wavelength	586 nm
Emission Wavelength	603 nm

## Applications

Texas Red® cadaverine 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.