

xtraFluor™ Violet 420 Amine

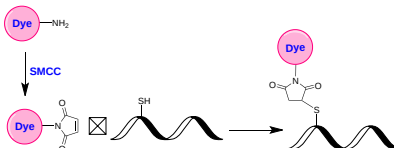
Catalog Number: 70806

Unit Size: 100 ug

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	1 year upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	N/A
Soluble In	DMSO
Chemical Structure	

Spectral Properties

Excitation Wavelength	407 nm
Emission Wavelength	424 nm

Applications

xtraFluor™ Violet 420 Amine is a violet laser-excitable fluorescent dye with excitation and emission maxima of approximately 407 nm and 424 nm. The dye contains a primary amine functionality that supports in situ maleimide generation through reaction with heterobifunctional crosslinkers such as SMCC. This reaction produces thiol-reactive maleimide-functional XV420 intermediates that can subsequently conjugate with sulfhydryl-containing biomolecules including reduced antibodies, peptides, and proteins.

XV420 provides strong fluorescence signal intensity for improved detection of dim or weakly expressed targets in complex samples. The dye is compatible with conventional and spectral flow cytometry workflows using appropriate violet laser detection configurations. Its spectral properties and high brightness support multicolor panel design requiring reduced spillover and improved population discrimination. The dye can be incorporated into custom thiol-reactive conjugation workflows requiring flexible fluorescent labeling strategies under standard bioconjugation conditions.