

mFluor Violet 480 SE

Catalog Number: 1158

Unit Size: 1 mg

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure,
Expiration Date	12 months upon receiving

Chemical Properties

Appearance	Solid
Molecular Weight	N/A
Soluble In	DMSO

Spectral Properties

Excitation Wavelength	N/A
Emission Wavelength	N/A

Applications

mFluor™ Violet 480 SE is reasonably stable and shows good reactivity and selectivity with protein amino groups. mFluor™ Violet 480 SE provides a convenient tool to label monoclonal, polyclonal antibodies or other proteins (>10 kDa) for flow cytometric applications with the violet laser excitation. These dyes have fluorescence excitation and emission maxima of ~405 nm and ~480 nm respectively. These spectral characteristics make them a unique color for flow cytometry application. mFluor™ dyes are developed for multicolor flow cytometry-focused applications. These dyes have large Stokes Shifts, and can be well excited by the laser lines of flow cytometers (e.g., 350 nm, 405 nm, 488 nm and 633 nm). mFluor™ Violet dyes are optimized to be excited with a violet laser at 405 nm. AAT Bioquest offers the largest collection of fluorescent dyes that are excited by violet laser.