

xtraFluor™ Violet 420 Labeling Dye

Catalog Number: 70800 70801,

Unit Size: 100 ug 1 mg,

Product Details

Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	24 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

xtraFluor™ Violet 420 (XV420) is an extremely bright violet laser-excitable fluorescent labeling dye. With a maximum excitation well matched the 405 nm violet laser and an emission peak at ~420 nm, XV420 can be conveniently used to replace BD Horizon Brilliant™ Violet 421 (BV421). It can be used on flow cytometers equipped with a violet laser and appropriate filters for large multicolor flow cytometry panels. It is a brighter alternative to Pacific Blue (ThermoFisher) and BD Horizon™ V450 (BD Biosciences). XV420 enables improved resolution for dim population identification. It is particularly useful in multicolor applications where it can be used to better resolve dim populations. XV420 dye offers new options for multicolor panel design through the use of a violet laser. Additionally, the dye's spillover properties, stability under light and in standard buffers, and compatibility with blood collection tubes offer ease-of-use for a range of applications. XV420 can be readily conjugated to antibodies through the well-known Buccutite™ chemistry that was developed by the scientists of AAT Bioquest. Compared to other conjugation chemistry, Buccutite™ chemistry can be run under extremely mild conditions (e.g., room temperature at neutral pH) with high conjugation yield.