

## xtraFluor™ Violet 420 Streptavidin Conjugate

Catalog Number: 70805

Unit Size: 100 ug

### Product Details

---

Storage Conditions	Minimize light exposure, Refrigerated (2-8 °C)
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 Streptavidin Conjugate might be one of the brightest fluorescent dye-labeled streptavidin conjugates commercially available. It can be conveniently used for detecting biotin-tagged biomolecules such as biotinylated antibodies, peptides and oligos. XV420 is an extremely bright violet laser-excitable fluorophore. 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.