

XFD647 AlkyneCatalog number: 1897
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	1010.09
Soluble In	DMSO

Spectral Properties

Excitation Wavelength	650 nm
Emission Wavelength	671 nm

Applications

XFD647 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 647 (Alexa Fluor® is the trademark of ThermoFisher). It is a bright deep red fluorescent dye. XFD647 dye is water soluble and pH-insensitive from pH 4 to pH 10. XFD647 alkyne is used for conjugation with a biological molecule that contains an azide via the click chemistry. It provides a convenient way to install the fluorophore of Alexa Fluor® 647 into a biological molecule that lacks alkyne and azide groups.