

XFD700 NHS Ester *Same Structure to Alexa Fluor™ 700 NHS Ester*Catalog number: 1836, 71911
Unit size: 1 mg, 5 mg**Product Details**

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

Chemical Properties

Appearance	Dark blue solid
Molecular Weight	~1400
Soluble In	DMSO

Spectral Properties

Excitation Wavelength	696 nm
Emission Wavelength	719 nm

Applications

XFD700 is manufactured by AAT Bioquest, and it has the same chemical structure of Alexa Fluor® 700 (Alexa Fluor® is the trademark of ThermoFisher). It is a bright red fluorescent dye. XFD700 dye is water soluble and pH-insensitive from pH 4 to pH 10. The NHS ester (or succinimidyl ester) of XFD700 is the most convenient amine-reactive form for conjugating this dye to a protein or antibody.