

5-FITC [Fluorescein-5-isothiocyanate] *CAS 3326-32-7*

Catalog number: 120, 121, 122 Unit size: 100 mg, 1 g, 10 g

Product Details	
Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving
Chemical Properties	
Appearance	Yellow solid
Molecular Weight	389.38
Soluble In	DMSO
Chemical Structure	$ \begin{array}{c} H O \\ + \\ + \\ + \\ + \\ + \\ + \\ + \\ + \\ + \\$
Spectral Properties	
Excitation Wavelength	491 nm
Emission Wavelength	516 nm

Applications

5-FITC and 6-FITC have very similar absorption and fluorescence spectra. However, the isomers may differ in their binding and reactivities to proteins, and the conjugates may migrate differently in electrophoretic gels. Thus, we offer highly purified single isomers. Cautions must be exercised for the storage of FITC conjugates.