

Fluorescein *CAS 2321-07-5*

Catalog Number: 1

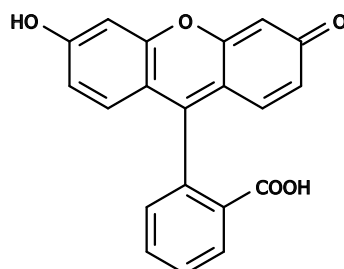
Unit Size: 1 g

Product Details

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

Chemical Properties

Appearance	Solid orange
Molecular Weight	332.31
Soluble In	DMSO
Chemical Structure	

**Spectral Properties**

Excitation Wavelength	498 nm
Emission Wavelength	517 nm

Applications

Fluorescein is a fluorophore widely used in microscopy, forensics and serology to detect latent blood stains. Fluorescein has an absorption maximum at 490 nm and emission maximum of 514 nm (in water). It has a pKa at 6.4 and multiple ionization equilibria. This leads to pH dependent absorption and emission over the range of 5 to 9. It has an isoabsorptive point at 460 nm.