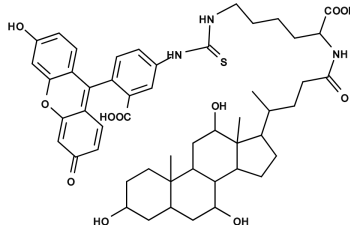


**CLF [Cholyl-Lys-Fluorescein]**Catalog number: 36701  
Unit size: 1 mg**Product Details**

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

**Chemical Properties**

Appearance	Yellow solid
Molecular Weight	926.12
Soluble In	DMSO
Chemical Structure	

**Spectral Properties**

Excitation Wavelength	498 nm
Emission Wavelength	517 nm

**Applications**

CLF is a fluorescein-labeled bile acid analog that closely parallels the cellular binding and uptake properties of cholic acid derivatives. It is a divalent 'unipolar' anionic fluorescent monohydroxy bile salt analog with physical, biological and cholestatic properties that are similar to those of lithocholate, glycolithocholate and their derivatives and thus offers a potentially useful probe for studying mechanisms of monohydroxy bile salt-induced cholestasis at the hepatocellular level. CLF is currently used as a tool for the visualization of bile acid transport within liver tissue.