

## Product Information Sheet

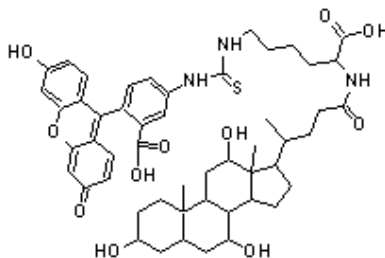
### Ordering Information

Product Number: 36701  
Product Name: CLF [Cholyl-Lys-Fluorescein]  
Unit Size: 1 mg  
Storage Conditions: <-15 °C and kept from light and moisture  
Expiration Date: 12 months upon receiving

### Chemical and Spectral Properties

Molecular Weight: 926.12

Chemical Structure:



Soluble in: DMSO  
Excitation Wavelength: 492 nm  
Emission Wavelength: 516 nm

### Application Notes

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.