

Fluorescein dicaprylate WS1

Catalog number: 22018

Unit size: 5 mg

Product Details

Storage Conditions Freeze (<-15 °C), Minimize light exposure

Expiration Date 12 months upon receiving

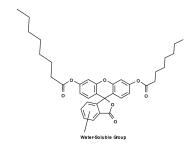
Chemical Properties

Appearance Off- white solid

Molecular Weight 828.97

Soluble In DMSO

Chemical Structure



Spectral Properties

Excitation Wavelength 498 nm

Emission Wavelength 517 nm

Applications

Fluorescein dicaprylate WS1 is a non-fluorescent water-soluble fluorescein derivative. Esterases hydrolyze the dicaprylate group producing the highly fluorescent product fluorescein. This water-soluble fluorescein dicaprylate may stain Salmonella in green fluorescence. It is a fluorogenic substrate for C8 esterase that generates green fluorescence upon esterase cleavage. Compared to the fluorescein dicaprylate esterase substrate, this fluorescein dicaprylate WS1 is much more water-soluble while fluorescein dicaprylate precipitates in aqueous solutions, making it hard to be used in continuous assays.