

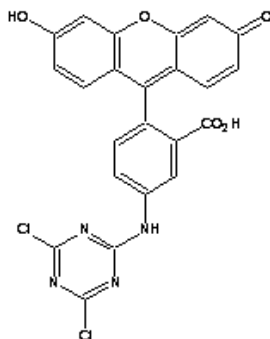
Product Information Sheet

Ordering Information

Product Number:	200
Product Name:	5-DTAF [5-(4,6-Dichlorotriazinyl)aminofluorescein]
Unit Size:	25 mg
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving

Chemical, Physical and Spectral Properties

Molecular Weight:	495.28
Appearance:	Yellow solid
Chemical Structure:	



Soluble in:	DMSO
Excitation Wavelength:	492 nm
Emission Wavelength:	517 nm

Application Notes

The 5-isomer of fluorescein dichlorotriazine (5-DTAF) is highly reactive with proteins. Unlike other reactive fluoresceins, 5-DTAF not only reacts with amino groups, but also reacts with thiol groups and even directly reacts with hydroxy groups such as polysaccharides and other alcohols in aqueous solution at pH above 9. Due to its high reactivity, 5-DTAF is also used to label carbohydrates.