

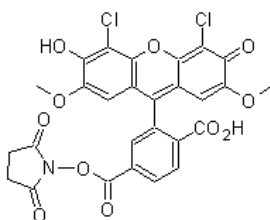
Product Information Sheet

Ordering Information

Product Number:	203
Product Name:	6-JOE, SE [6-Carboxy-4',5'-dichloro-2',7'-dimethoxyfluorescein, succinimidyl ester]
Unit Size:	5 mg
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving

Chemical, Physical and Spectral Properties

Molecular Weight:	602.34
Appearance:	Red solid
Chemical Structure:	



Soluble in:	DMF or DMSO
Excitation Wavelength:	520 nm
Emission Wavelength:	548 nm

Application Notes

6-JOE has similar absorption and emission spectra to those of rhodamine 6G, and its SE is predominantly used in automated DNA sequencing. It is also used to label nucleotides.