

6-JOE, SE [6-Carboxy-4',5'-dichloro-2',7'-dimethoxyfluorescein, succinimidyl ester]***CAS#: 113394-23-3***

Catalog number: 203

Unit size: 5 mg

Product Details

Storage Conditions	Freeze (<-15 °C), Minimize light exposure
Expiration Date	12 months upon receiving

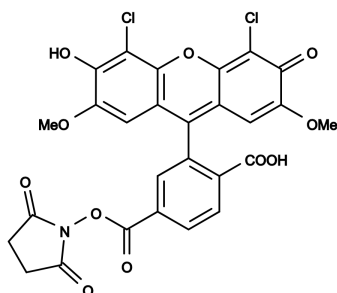
Chemical Properties

Appearance Orange solid

Molecular Weight 602.34

Soluble In DMSO

Chemical Structure

**Spectral Properties**

Excitation Wavelength 520 nm

Emission Wavelength 545 nm

Applications

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.