

OG514 succinimidyl ester [equivalent to Oregon Green® 514 succinimidyl ester]

Catalog Number: 717

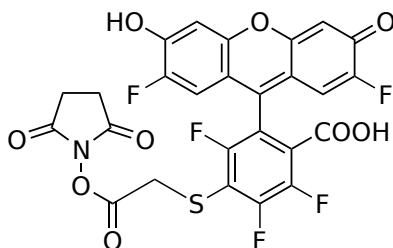
Unit Size: 5 mg

Product Details

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

Chemical Properties

Appearance	Solid orange
Molecular Weight	609.43
Soluble In	DMSO
Chemical Structure	



Spectral Properties

Excitation Wavelength	513 nm
Emission Wavelength	533 nm

Applications

OG514 is the same fluorophore as Oregon Green 514 (Oregon Green® is the trademark of ThermoFisher). OG514 Succinimidyl Ester is an amine-reactive compound used for creating green, fluorescent bioconjugates. It has excitation/emission maxima approximately 489/526 nm. It is an excellent replacement for rhodamine 6G that has poor solubility in most common solvents, thus practically hard to use. Although it is a fluorescein derivative, it has significantly improved photostability and a lower pKa, making its fluorescence essentially pH insensitive in the physiological pH range. OG514 Succinimidyl Ester can be conveniently used to label antibodies, peptides, amino-modified oligos.