

OG514 acid [equivalent to Oregon Green® 514 carboxylic acid]

Catalog Number: 716

Unit Size: 25 mg

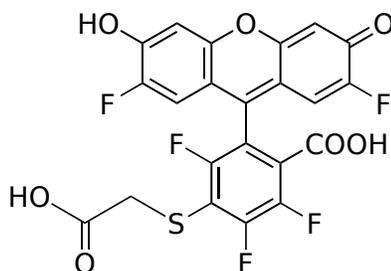
Product Details

| | |
|--------------------|--------------------------------------------|
| Storage Conditions | Freeze (< -15 °C), Minimize light exposure |
| Expiration Date | 24 months upon receiving |

Chemical Properties

| | |
|------------------|--------------|
| Appearance | Solid yellow |
| Molecular Weight | 512.36 |
| 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 acid 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.