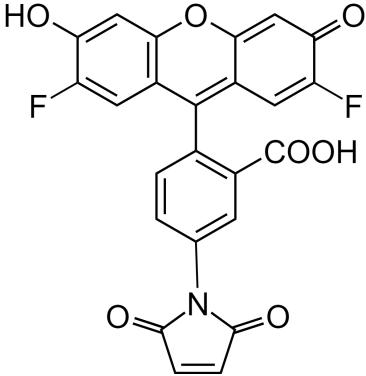
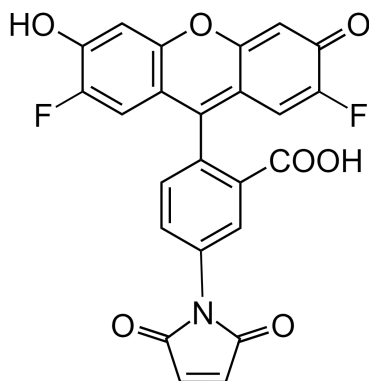


OG488 maleimide [equivalent to Oregon Green® 488 maleimide]Catalog number: 715
Unit size: 5 mg**Product Details**

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

Chemical Properties

Appearance	Red solid
Molecular Weight	463.35
Soluble In	DMSO
Chemical Structure	

**Spectral Properties**

Excitation Wavelength	498 nm
Emission Wavelength	526 nm

Applications

OG488 maleimide is the same molecule to Oregon Green® 488 maleimide (Invitrogen #O6034, Oregon Green® is the Trademark of Invitrogen). This fluorinated analog of fluorescein has higher photostability and a lower pKa ~ 4.7 while fluorescein maleimide has pKa ~6.4. The decreased pKa of OG488 maleimide makes its fluorescence essentially pH insensitive in the physiological pH range. This thiol-reactive OG488 maleimide can be used to make bright green fluorescent bioconjugates with excitation/emission maxima ~496/524 nm. It is an excellent replacement to fluorescein maleimide that has severe pH dependence and low photostability.