

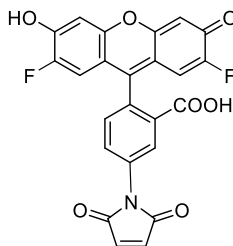
## Product Information Sheet

### Ordering Information

Product Number:	715
Product Name:	OG488 maleimide [equivalent to Oregon Green® 488 maleimide] *Single isomer*
Unit Size:	5 mg
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving

### Chemical and Spectral Properties

Molecular Weight:	463.35
-------------------	--------



Chemical Structure:

Soluble in:	DMSO
Excitation Wavelength:	496 nm
Emission Wavelength:	524 nm

### Application Notes

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.