

## OG488 Taxol Conjugate

Catalog Number: 23175

Unit Size: 5x50 ug

### Product Details

---

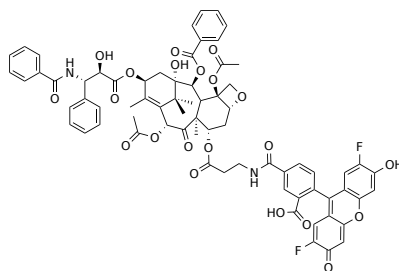
Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	24 months upon receiving

### Chemical Properties

---

Appearance	Solid orange
Molecular Weight	1319.28
Soluble In	DMSO

Chemical Structure



### Spectral Properties

---

Excitation Wavelength	498 nm
Emission Wavelength	526 nm

### Applications

---

OG488 Taxol Conjugate is the same compound as Oregon Green® 488 paclitaxel (also known as Flutax-2). Taxol, also known as Paclitaxel, is an anticancer agent. Taxol is known to promote tubulin assembly into stable aggregated structures. The green fluorescent OG488 Taxol Conjugate is an effective fluorescent probe for studying microtubules in cells and in solutions. The fluorescent label OG488 is attached to paclitaxel by derivatization of the 7-β-hydroxy group of native paclitaxel, minimizing the interference of OG488 fluorophore on the binding of the probe to microtubules. OG488 Taxol Conjugate has been successfully used for live-cell tubulin labeling. The binding of OG488 Taxol Conjugate to microtubules can be conveniently analyzed either by FACS or fluorescence imaging.