

## CypHer5E Maleimide

Catalog Number: 21913

Unit Size: 1 mg

### Product Details

---

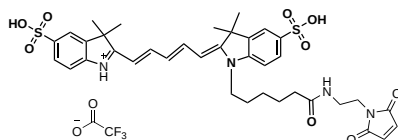
Storage Conditions	Freeze (< -15 °C), Minimize light exposure
Expiration Date	24 months upon receiving

### Chemical Properties

---

Appearance	Solid blue
Molecular Weight	864.90
Soluble In	DMSO

Chemical Structure



### Spectral Properties

---

Excitation Wavelength	643 nm
Emission Wavelength	660 nm

### Applications

---

CypHer5E maleimide is an excellent building block for developing site-specific pH probes. It can be readily conjugated to thiol-containing biomolecules (such as antibodies, peptides and thiol-modified oligos). CypHer5E is a red laser-excitable, pH-sensitive cyanine dye derivative that can be used to study most classes of cell surface receptors which internalize into the endocytic pathway when activated by a ligand. CypHer5E maleimide is a flexible, generic fluorophore that can be coupled to a receptor or epitope tag antibody and monitors neutral to acidic intracellular events. CypHer5E has an excitation peak that perfectly matches the red laser at 633 nm of He-Ne laser and 647 nm of red laser and an emission peak at ~670 nm.