

## Protonex™ NIR 770 NHS ester

Catalog Number: 21221

Unit Size: 1 mg

### Product Details

---

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

### Chemical Properties

---

Appearance	Solid
Molecular Weight	535.64
Soluble In	DMSO

### Spectral Properties

---

Excitation Wavelength	748 nm
Emission Wavelength	769 nm

### Applications

---

Protonex™ NIR 770 NHS ester is the only cell-permeable acidotropic pH dye commercially available for developing NIR pH probes. It is an excellent building block that can be readily used for developing site-specific pH sensors for monitoring pH fluctuation in a spatial environment. It can be readily conjugated to amino-containing biomolecules (such as antibodies, peptides and amino-modified oligos). It easily penetrates live cell membranes, thus can be used for monitoring pH changes in live cells. It is complimentary to CypHer7E that is cell-impermeable, and often used to study most classes of cell surface receptors which internalize into the endocytic pathway when activated by a ligand. Protonex™ NIR 770 NHS ester is a flexible, generic fluorophore that can be coupled to a receptor or epitope tag antibody and monitors neutral to acidic intracellular events. It is one of the very few acidotropic pH probes that exhibit NIR fluorescence.