

**Protonex™ Red 780 AM \*Cell-Permeable\***

Catalog Number: 21186

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	798.89
Soluble In	DMSO

**Spectral Properties**

---

Excitation Wavelength	748 nm
Emission Wavelength	769 nm

**Applications**

---

Protonex™ Red 780 AM is the cell-permeable version of Protonex™ Red 780. Protonex™ Red 780 dye works by changing its fluorescence intensity depending on the pH of the environment. Protonex™ Red 780 is minimally fluorescent at a basic pH and maximally fluorescent at an acidic pH. When Protonex™ Red 780 is bound to an acidic intracellular target, it becomes highly fluorescent and emits near-infrared light when excited by a red laser. By measuring the fluorescence intensity of Protonex™ Red 780, one can label or monitor the acidic intracellular targets in live cells.