

## Protonex™ Red 670 AM \*Cell-Permeable\*

Catalog Number: 21182

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 | 772.85 |
| Soluble In       | DMSO   |

### Spectral Properties

---

|                       |     |
|-----------------------|-----|
| Excitation Wavelength | N/A |
| Emission Wavelength   | N/A |

### Applications

---

Protonex™ Red 670 AM is the cell-permeable version of Protonex™ Red 670. Protonex™ Red 670 dye works by changing its fluorescence intensity depending on the pH of the environment. Protonex™ Red 670 is minimally fluorescent at a basic pH and maximally fluorescent at an acidic pH. When Protonex™ Red 670 is bound to an acidic intracellular target, it becomes highly fluorescent and emits red light when excited by a red laser such as a 632 nm He-Ne or 647 nm red laser. By measuring the fluorescence intensity of Protonex™ Red 670, one can label or monitor the acidic intracellular targets in live cells.