

**Protonex™ Lyso-NIR 770**

Catalog Number: 21242

Unit Size: 1 mg

**Product Details**

---

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

**Chemical Properties**

---

Appearance	Solid
Molecular Weight	736.76
Soluble In	DMSO

**Spectral Properties**

---

Excitation Wavelength	748 nm
Emission Wavelength	769 nm

**Applications**

---

Protonex™ Lyso-NIR 770 is a fluorescent derivative prepared from our Protonex™ NIR 770 NHS ester. It is cell-permeable and selectively localized in acidic organelles such as lysosomes. It is the only cell-permeable acidotropic pH dye with NIR fluorescence for monitoring pH changes in acidic organelles. Its fluorescence can be readily captured with the common Cy7 filter set that is equipped with most of the fluorescence instruments. Due to its significantly red-shifted excitation and emission wavelengths, it has much lower background than the existing pH sensors (such as LysoSensors). It might be used in vivo for monitoring pH changes in animals due to the significantly improved skin-penetration capability of NIR excitation.