

**PRO-TAO-3 [equivalent to TO-PRO-3]**

Catalog number: 17565

Unit size: 1 mL

**Product Details**

---

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

**Chemical Properties**

---

Appearance	Liquid
Molecular Weight	671.42
Soluble In	Water

**Spectral Properties**

---

Excitation Wavelength	642 nm
Emission Wavelength	657 nm

**Applications**

---

PRO-TAO-3 is the same molecule to TO-PRO®-3 stain. TO-PRO®-3 and the other monomeric cyanine stains have wide applicability due to their low background and bright fluorescence. It is a carbocyanine monomer nucleic acid stain that can be well excited by a red laser at 633 nm or 647 nm with far-red fluorescence similar to Cy®5 dyes. It is useful as a nuclear counterstain and dead cell indicator. It is selective for detecting dsDNA (double-stranded nucleic acids). The cell-impermeable PRO-TTO-3 is widely used for flow cytometry and fluorescence microscopy as a long-wavelength fluorescence well separated from green and red fluorophores in multicolor applications. TO-PRO® is the trademark of ThermoFisher Scientific.