

**DiPO™-1 [equivalent to POPO®-1] \*1 mM  
DMSO Solution\***Catalog number: 17582  
Unit size: 100 ul**Product Details**

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

**Chemical Properties**

Appearance	Liquid
Molecular Weight	1170.54
Soluble In	Water

**Spectral Properties**

Excitation Wavelength	433 nm
Emission Wavelength	457 nm

**Applications**

DiPO™-1 is the same compound to POPO®-1 (POPO® is the trademark of ThermoFisher). DiPO™-1 is a high-affinity nucleic acid stain based upon a symmetric cyanine dye dimer that is essentially nonfluorescent in the absence of nucleic acids that exhibits excitation/emission maxima ~434/456 nm when bound to nucleic acids. POPO-1 is a Blue emitting dye that can be excited by the 405 laser and emission captured by the 450/50 filter. It is commonly used as a nucleic acid stain in flow cytometry and fluorescence microscopy.