

Product Information Sheet

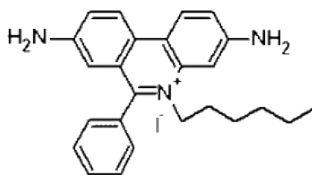
Ordering Information

Product Number:	17562
Product Name:	Hexidium Iodide
Unit Size:	5 mg
Storage Conditions:	<-15 °C and kept from light and moisture
Expiration Date:	12 months upon receiving

Chemical and Spectral Properties

Molecular Weight: 497.41

Chemical Structure:



Soluble in: DMSO

Excitation Wavelength: 518 nm

Emission Wavelength: 600 nm

Application Notes

The nucleic acid stain, hexidium iodide is permeant to mammalian cells and selectively stains almost all gram-positive bacteria in the presence of gram-negative bacteria. Hexidium iodide exhibits excitation/emission maxima of ~518/600 nm upon binding DNA. Generally, both the cytoplasm and nuclei of eukaryotic cells show staining with hexidium iodide. We do not recommend its use as a nuclear stain since it also stains other subcellular compartments.