

CYTO-16 *equivalent to SYTO 16*

Catalog Number: 17568

Unit Size: 0.5 mL

Product Details

Storage Conditions Freeze (< -15 °C), Minimize light exposure

Expiration Date 24 months upon receiving

Chemical Properties

Appearance Liquid

Molecular Weight 565.52

Soluble In DMSO

Spectral Properties

Excitation Wavelength 489 nm

Emission Wavelength 520 nm

Applications

CYTO 16 is manufactured by AAT Bioquest to replace SYTO 16 Green Nucleic Acid Stain. They have virtually the same chemical and biological properties. CYTO 16 is a cell-permeant nucleic acid stain that exhibits bright, green fluorescence upon binding to nucleic acids. The CYTO 16 dye can be used to stain RNA (Ex/Em ~ 494/525 nm) and DNA (Ex/Em ~ 488/518 nm) in both live and dead eukaryotic cells, as well as in Gram-positive and Gram-negative bacteria. Eukaryotic cells incubated with SYTO 16 dye generally show cytoplasmic or mitochondrial staining, as well as nuclear staining.