

Product Information

Acridine orange

Catalog Number	Unit Size
C014	10 mg/mL, 1 mL

Storage upon receipt:

- -20°C
- Protect from light

Ex/Em: 500/526 nm, bound to DNA

460/650 nm, bound to RNA

Product Description

Acridine orange is a cell-permeant nucleic acid binding dye that emits green fluorescence when bound to dsDNA and red fluorescence when bound to ssDNA or RNA. This unique characteristic makes acridine orange useful for cell-cycle studies. Acridine orange has also been used as a lysosomal dye.

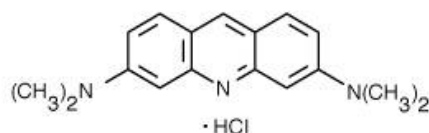
Specifications

Chemical Name: acridine orange

Molecular Formula: C₁₇H₂₀ClN₃

Molecular Weight: 301.82

CAS Number: 65-61-2



Application

Nuclear counterstaining.