Acridine Orange, 10 mg/mL in Water (High Purity) (AO)

Acridine orange (AO) stains dsDNA green (525 nm) and RNA or single stranded DNA red (650 nm). The dye is membrane-permeant and its nucleic acid binding property has been used for cell-cycle studies.



Product attributes

Probe cellular localization	Nucleus & cytoplasm
For live or fixed cells	For fixed cells, For live/intact cells
Assay type/options	No-wash staining
Cell permeability	Membrane permeant
Colors	Green/Red
Excitation/Emission	500/526 pm (DNA) 460/650 pm (RNA)

Call us: 800-304-5357 Email: btinfo@biotium.com

Product Description

Acridine orange (AO) stains dsDNA green (525 nm) and RNA or single stranded DNA red (650 nm). The dye is membrane-permeant and its nucleic acid binding property has been used for cell-cycle studies.^{2,3} Acridine orange has also been used for the detection of microorganisms in cerebrospinal fluid and other clinical specimens. We offer a highly purified form of acridine orange while most of the other commercially available grades of AO are either in zinc chloride complex form or of low purity.

- $\lambda_{Ex}/\lambda_{Em} = 500/526 \text{ nm (DNA)}, 460/650 \text{ nm (RNA)}$
- Yellow solution
- Store at 4°C and protect from light
- C₁₇H₂₀CIN₃
- MW: 301.82

References

- 1. Cytometry 12(4), 330 (1991).
- 2. Methods Cell Biol 33, 285 (1990).
- 3. J Cell Physiol 143(2), 279 (1990).
- 4. J. Clin. Microbiol 14(2), 201 (1981).

This datasheet was generated on November 17, 2025 at 08:07:10 AM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/acridine-orange-10-mgml-solution-in-water-high-purity-ao/