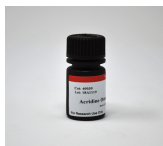


## Acridine Orange, 10 mg/mL in Water (High Purity) (AO)

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.



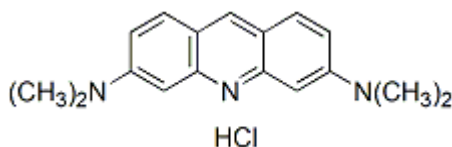
### Product attributes

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 nm (DNA), 460/650 nm (RNA)

## 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.<sup>2,3</sup> 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 nm (DNA), 460/650 nm (RNA)
- Yellow solution
- Store at 4 °C and protect from light
- $C_{17}H_{20}ClN_3$
- 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 January 8, 2026 at 12:39:58 PM. 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/>