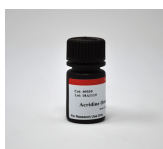


Acridine Orange, 10 mg/mL in Water (High Purity) (AO)

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.



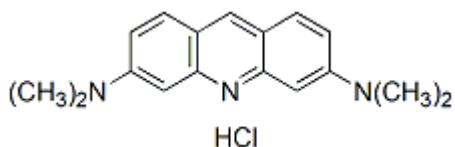
Product attributes

| | |
|-----------------------------|--|
| 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.^{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 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 June 4, 2026 at 10:42:29 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/>