

www.biotium.com

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

## **PRODUCT AND SAFETY DATA SHEET**

**PRODUCT NAME:** 

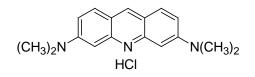
Acridine Orange, 10mg/ml solution in water (high purity)

CATALOG #

**MOLECULAR INFORMATION:** 

 $C_{17}H_{20}ClN_3$ MWt: 301.82 [65-61-2]

40039



**PROPERTIES:** 

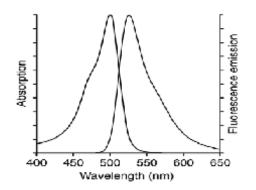
| Color & Form        | Yellow solution  |
|---------------------|------------------|
| Purity              | ≥98% by HPLC     |
| Solubility          | Soluble in water |
| Absorption/Emission | 493/526nm        |
| Extinction          | >35,000          |
| Coefficient         |                  |

STORAGE AND HANDLING:

Stored at 4 °C and protect from light

**APPLICATION:** 

Acridine orange (AO) stains dsDNA green (525 nm) and RNA or single stranded DNA red (650 nm)<sup>1</sup>. 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.<sup>4</sup> 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.



Absorption and emission of acridine orange (AO) bound to DNA.

Ref.: 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).

**TOXICITY:** 

Unknown

## Biotium

46117 Landing Parkway, Fremont, CA 94538 U.S.A. Tel: 1-510-265-1027; Fax: 1-510-265-1352

www.biotium.com

| FIRST AID: | Potentially harmful. Avoid prolonged or repeated exposure. Avoid getting in eyes, on skin, or on  |
|------------|---|
|            | clothing. Wash thoroughly after handling. If eye or skin contact occurs, wash affected areas with |
|            | plenty of water for 15 minutes and seek medical advice. In case of inhaling or swallowing,        |
|            | move individual to fresh air and seek medical advice immediately.                                 |
|            |   |

**Disclaimer:** *Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use. Biotium is not liable for any damage resulting from handling or contact with this product.*