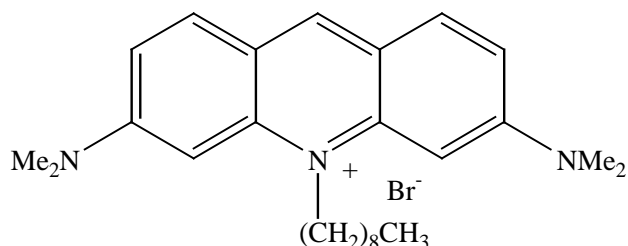


## PRODUCT AND SAFETY DATA SHEET

**PRODUCT NAME:** Nonyl Acridine Orange (NAO)**CATALOG NUMBER:** 70012**MOLECULAR INFORMATION:** C<sub>26</sub>H<sub>38</sub>BrN<sub>3</sub>  
Mwt: 473**PROPERTIES:**

<b>Color &amp; Form</b>	Orange solid
<b>Purity</b>	≥ 95% by HPLC
<b>Solubility</b>	Soluble in DMSO
$\lambda_{ex} \backslash \lambda_{em}$	495nm \ 519nm (MeOH)
<b>Extinction Coefficient (<math>\epsilon</math>)</b>	84,000

**STORAGE AND HANDLING:**

Store at 4 °C and protect from light, especially when in solution.

**APPLICATION:**

NAO is a mitochondrial dye whose staining is not dependent on mitochondrial membrane potential,<sup>1</sup> unlike JC-1 and rhodamine 123 (See **70010** and **70011**). NAO has been used to study multidrug resistance<sup>2</sup> and to measure changes in mitochondrial mass during apoptosis.<sup>3</sup>

Ref.: 1) *Cytometry* **9**, 206(1988); 2) *Cancer Res.* **51**, 4665(1991); 3) *Exp. Cell Res.* **214**, 323(1994).**TOXICITY:**

Unknown

**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.