

## Nonyl Acridine Orange (NAO)

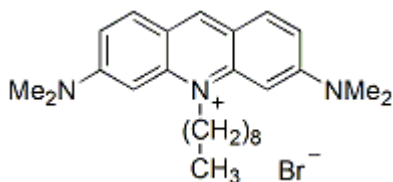
NAO is a green fluorescent mitochondrial dye whose staining is not dependent on mitochondrial membrane potential.



## Product Description

NAO is a green fluorescent mitochondrial dye whose staining is not dependent on mitochondrial membrane potential (1), unlike JC-1 ([70011](#)) or Rhodamine 123 ([70010](#)). NAO has been used to study multidrug resistance (2) and to measure changes in mitochondrial mass during apoptosis (3). The dye is reported to bind cardiolipin, a phosphoprotein in the inner mitochondrial membrane (4). Also see our [MitoView™ Dyes and other mitochondrial dyes](#).

- $\lambda_{Ex}/\lambda_{Em}$  (MeOH) = 495/522 nm
- $\epsilon$  (MeOH) = 63,000
- Orange solid soluble in DMSO or DMF
- Store at 4 °C and protect from light, especially in solution
- $C_{26}H_{38}BrN_3$
- MW: 473
- [75168-11-5]



## References

1. Cytometry 9, 206 (1988).
2. Cancer Res 51, 4665 (1991).
3. Exp Cell Res 214, 323 (1994).
4. FEBS Letters 260, 236 (1990).

This datasheet was generated on January 4, 2026 at 08:21:21 PM. Visit product page to check for updated information before use.  
Product link: <https://biotium.com/product/nonyl-acridine-orange-nao/>

## Product attributes

CAS number	75168-11-5
Probe cellular localization	Mitochondria
For live or fixed cells	For live/intact cells
Assay type/options	No-wash staining, Real-time imaging
Potential dependence	Mitochondrial potential-independent
Colors	Red
Excitation/Emission	495/522 nm