

AO Succinimidyl Ester

The amine reactive succinimidyl ester of AO (acridine orange) can be conjugated to peptides, proteins, drugs, polymeric materials and biomolecules with primary amine groups. The conjugates with green fluorescence are able to complex with nucleic acids, resulting in green fluorescence nucleic acid conjugate adduct.



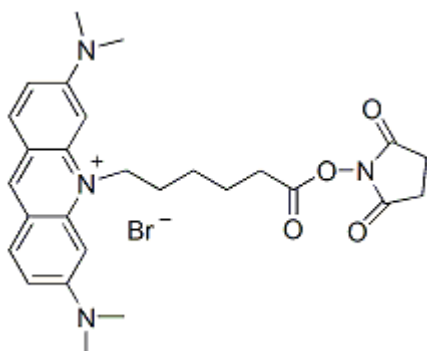
Product attributes

Excitation/Emission	500/526 nm (DNA), 460/650 nm (RNA)
---------------------	------------------------------------

Product Description

The amine reactive succinimidyl ester of green fluorescent AO (acridine orange) can be conjugated to peptides, proteins, drugs, polymeric materials and biomolecules with primary amine groups. The conjugates then are able to complex with nucleic acids, resulting in green fluorescence nucleic acid conjugate adduct, making them potentially useful for studies of nucleic acid binding to various biomolecules, such as DNA-binding proteins. It is also possible that conjugates of other biomolecules might be capable of monitoring their transport into the nucleus. AO dye conjugates of solid or semisolid matrices, such as microspheres, magnetic particles or various resins, might be useful for the detection or affinity isolation of nucleic acids.

- $\lambda_{Ex}/\lambda_{Em}$ (with DNA): 500/526 nm
- Orange solid soluble in DMF or DMSO
- Store at -20°C , desiccated and protected from light
- $\text{C}_{27}\text{H}_{33}\text{BrN}_4\text{O}_4$
- MW: 557.49



See our other [reactive DNA/RNA binding dyes](#).