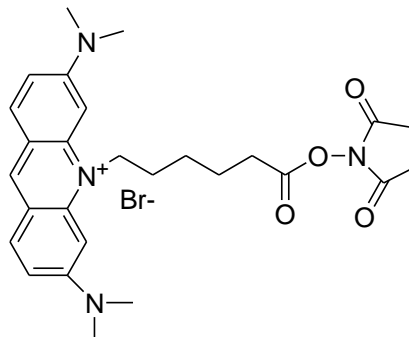


PRODUCT AND SAFETY DATA SHEET

PRODUCT NAME: AO Succinimidyl ester

CATALOG # 40074

MOLECULAR INFORMATION: C₂₇H₃₃BrN₄O₄
MWt: 557.49



PROPERTIES:

Color & Form	Orange solid
Solubility	Soluble in DMSO or DMF
Absorption / Emission	500nm / 526nm (with DNA)
Correction factor 260nm	0.78
Correction factor 280nm	0.43

STORAGE AND HANDLING: Stored desiccated at -20°C and protect from light, especially when in solution.

APPLICATION: 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. The conjugates might be 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.

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.