

TO Succinimidyl Ester

The amine reactive succinimidyl ester of TO (thiazole orange) can be conjugated to peptides, proteins, drugs, polymeric materials and biomolecules with primary amine groups. The conjugates are expected to be essentially nonfluorescent until they are able to complex with nucleic acids, resulting in green fluorescence.



Product attributes

Excitation/Emission 512/533 nm (with DNA)

Call us: 800-304-5357

Product Description

The amine reactive succinimidyl ester of TO (thiazole orange) can be conjugated to peptides, proteins, drugs, polymeric materials and biomolecules with primary amine groups. The conjugates are expected to be essentially nonfluorescent until they are able to complex with nucleic acids, resulting in green fluorescence. 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. TO 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.

- λ_{Ex}/λ_{Em} (with DNA): 512/533 nm
- Orange solid soluble in DMF or DMSO
- C₂₈H₂₈N₃O₄S
- Store at -20°C, desiccated and protected from light
- MW: 503

See our other reactive DNA/RNA binding dyes.

This datasheet was generated on November 21, 2025 at 02:31:59 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/to-succinimidyl-ester/