



Glowing products for science

Call us : [800-304-5357](tel:800-304-5357)

Cyanine Dye Succinimidyl Ester

Biotium offers high quality amine-reactive succinimidyl ester (SE) dyes used to label proteins, oligonucleotides, and nucleic acids. We offer Cyanine 555, Cyanine 647, Cyanine 680, and Cyanine 750 as high quality succinimidyl esters.



Product attributes

Product Description

Biotium offers high quality amine-reactive succinimidyl ester (SE, or NHS-Ester) dyes used to label proteins, oligonucleotides, and nucleic acids. We offer Cyanine 555, Cyanine 647, Cyanine 680, and Cyanine 750 as high quality succinimidyl esters. Also see our [CF® Dye Succinimidyl Esters](#), available with more than 20 bright and photostable CF® Dyes.

This datasheet was generated on June 10, 2026 at 12:10:15 PM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/cyanine-dye-succinimidyl-ester/>