



Glowing products for science

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

Cyanine NS Dye Succinimidyl Ester

Cyanine NS succinimidyl esters are non-sulfonated cyanine dyes that can be used to label proteins, oligonucleotides and nucleic acids.



Product attributes

Product Description

Cyanine NS succinimidyl esters are non-sulfonated cyanine dyes that can be used to label proteins, oligonucleotides and nucleic acids. Available in 3 colors, Cyanine 488NS, Cyanine 555NS, and Cyanine 647NS are non-sulfonated versions of Cyanine 488, Cyanine 555, and Cyanine 647, respectively. They are designed to match by charge and molecular weight for 2D DIGE platform. Biotium offers these dyes as high quality amine-reactive succinimidyl esters.

Also see sulfonated [Cyanine Dye Succinimidyl Esters](#) and our [CF@ Dye Succinimidyl Esters](#), available with more than 20 bright and photostable CF® Dyes.

This datasheet was generated on June 5, 2026 at 05:45:48 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cyanine-ns-dye-succinimidyl-ester/>