



Glowing products for science

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

Cyanine Dye dUTP

Fluorescent dUTP can be used to synthesize labeled DNA probes for in-situ hybridization, microarray or blotting techniques.



Product attributes

Product Description

Cyanine 555-dUTP and Cyanine 647-dUTP can be used to synthesize labeled DNA probes for in-situ hybridization, microarray or blotting techniques.

Also see our [CF® Dye dUTP Conjugates](#), available with a wide selection of bright and photostable CF® Dyes.

This datasheet was generated on June 13, 2026 at 09:08:03 AM. Visit product page to check for updated information before use.
Product link: <https://biotium.com/product/cyanine-dye-dutp/>