## Cyanine Dye dUTP

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



## **Product attributes**

Call us: 800-304-5357 Email: btinfo@biotium.com

## **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 November 15, 2025 at 05:26:51 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/cyanine-dye-dutp/">https://biotium.com/product/cyanine-dye-dutp/</a>