



Glowing products for science

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

## 5-TAMRA-dUTP, 1 mM Solution

5-TAMRA-dUTP (5-Tetramethylrhodamine-dUTP) can be enzymatically incorporated into DNA. Note: 5-TAMRA-dUTP may not be compatible with PCR labeling of DNA probes.



### Product attributes

Excitation/Emission	553/577 nm
---------------------	------------

## Product Description

5-TAMRA-dUTP (5-Tetramethylrhodamine-dUTP) can be enzymatically incorporated into DNA. Note: 5-TAMRA-dUTP may not be compatible with PCR labeling of DNA probes.

- 1 mM solution in Tris pH 7.5 buffer
- MW: 990.5
- $\lambda_{\text{Ex}}/\lambda_{\text{Em}}$ : 553/577 nm

This datasheet was generated on January 9, 2026 at 07:11:08 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/5-tetramethylrhodamine-dutp-1-mm-in-ph-7-5-tris-hcl-buffer/>