Glowing products for science

## 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

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

## **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
- λ<sub>Ex</sub>/λ<sub>Em</sub>: 553/577 nm

This datasheet was generated on November 12, 2025 at 12:40:52 AM. Visit product page to check for updated information before use. Product link: <a href="https://biotium.com/product/5-tetramethylrhodamine-dutp-1-mm-in-ph-7-5-tris-hcl-buffer/">https://biotium.com/product/5-tetramethylrhodamine-dutp-1-mm-in-ph-7-5-tris-hcl-buffer/</a>