



Glowing products for science

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

dNTP Mix, 25 mM Each

Highly purified mix of dNTPs (>99% by HPLC) in 0.6 mM Tris-HCl (pH 7.5); each dNTP is at a concentration of 25 mM in the solution provided.



Product attributes

Product Description

Highly purified mix of dNTPs (>99% by HPLC) in 0.6 mM Tris-HCl (pH 7.5); each dNTP is at a concentration of 25 mM in the solution provided. The dNTPs are suitable for use in PCR, cDNA synthesis, sequencing, and other molecular biology applications.

This datasheet was generated on January 14, 2026 at 07:54:13 AM. Visit product page to check for updated information before use.

Product link: <https://biotium.com/product/dntp-mix-25-mm/>