Highly purified mix of dNTPs (>99% by HPLC) in 0.6 mM Tris-HCl (pH 7.5); each dNTP is at a concentration of 10 mM in the solution provided.



Product attributes

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

Product Description

dNTP Mix, 10 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 10 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 November 11, 2025 at 07:33:23 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/dntp-mix-10-mm/