

dNTP Set, 100 mM Each

Highly purified dNTPs (>99% by HPLC) in purified water. Each deoxynucleotide solution (dATP, dCTP, dGTP, dTTP) is provided in a separate vial at 100 mM.



Product attributes

Product Description

Highly purified dNTPs (>99% by HPLC) in purified water. Each deoxynucleotide solution (dATP, dCTP, dGTP, dTTP) is provided in a separate vial at 100 mM. The dNTPs are suitable for use in PCR, cDNA synthesis, sequencing, and other molecular biology applications.

This datasheet was generated on November 3, 2025 at 07:06:56 PM. Visit product page to check for updated information before use. Product link: https://biotium.com/product/dntp-set-100-mm-each/