

Product Information

dNTP Set, 100 mM each

Catalog Number: 40052

Unit Size: 4 x 250 uL

Kit Contents

Component	Size
40052A: dATP, 100 mM	250 uL
40052C: dCTP, 100 mM	250 uL
40052G: dGTP, 100 mM	250 uL
40052T: dTTP, 100 mM	250 uL

Color and Form: Clear solutions

Storage and Handling

Store at -20°C. Product is stable for at least 1 year from date of receipt when stored as recommended.

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.

Related Products

Catalog number	Product
29050	Cheetah™ HotStart Taq DNA Polymerase, 500 U
31000-T	EvaGreen® Dye, 20X in water, 1 mL
29052	ROX Reference Dye, 25 uM in TE Buffer, 5 x 1 mL
31041-T	Forget-Me-Not™ qPCR Master Mix, 100 reactions
31042-T	Forget-Me-Not™ qPCR Master Mix with ROX, 100 reactions
31003-T	Fast EvaGreen® qPCR Master Mix, 100 reactions
29054-T	HotStart Polymerase Modification Kit, Trial Size
40008	CF™488A-dUTP, 25 nmol
40006	CF™594-dUTP, 25 nmol
40007	CF™640R-dUTP, 25 nmol
40029	Biotin-11-dUTP, 1 mM solution, 50 uL
41003-T	GelRed Nucleic Acid Gel Stain, 10,000X in water, 0.1 mL
41005-T	GelGreen Nucleic Acid Gel Stain, 10,000X in water, 0.1 mL
31030-50	DNA Gel Extraction Kit, 50 columns
31021	1 KB DNA Ladder, 30 ug/300 uL
31031	100 bp DNA Ladder, 30 ug/300 uL
41006	TBE Buffer, 5X, 4 L

Please visit our website at www.biotium.com for information on our life science research products, including environmentally friendly EvaGreen® qPCR master mixes, fluorescent CF™ dye antibody conjugates and reactive dyes, apoptosis reagents, fluorescent probes, and kits for cell biology research.

Materials from Biotium are sold for research use only, and are not intended for food, drug, household, or cosmetic use.