

100 mM dNTP Set, PCR Grade

Cat. nos.	Size	Store at -30°C to -10°C
10297-018	4 × 250 μL	
10297-117	8 × 1.25 mL	
Pub. Part no. 10297.pps	MAN0000810	Rev. Date 7 Sep 2011

Description

100 mM dNTP (2'-deoxynucleoside 5'-triphosphate) Set consists of all four deoxynucleotides (dATP, dCTP, dGTP, dTTP), each at a concentration of 100 mM. The deoxynucleotides are suitable for use in polymerase chain reaction (PCR), sequencing, fill-in, nick translation, cDNA synthesis, and TdT tailing reactions. The product is supplied as ready-to-use solutions.

Cat. no. 10297-018 includes 100 mM each of dATP, dCTP, dGTP, and dTTP. Each vial contains 250 μL 25 μmol dNTP in purified water (pH 7.5).

Cat. no. 10297-117 includes 100 mM each of dATP, dCTP, dGTP, and dTTP. Each dNTP is provided in 2 × 1.25-mL vials containing 250 μmol dNTP in purified water (pH 7.5).

Volumes and Concentrations

20 μL Final Reaction Volume

Final dNTP Conc	dNTP Volume	Rxns per kit	
		Cat. no. 10297-018	Cat. no. 10297-117
0.2 mM	0.04 μL	6,250	62,500
0.5 mM	0.1 μL	2,500	25,000
1.0 mM	0.2 μL	1,250	12,500
1.5 mM	0.3 μL	833	8,333

25 μL Final Reaction Volume

Final dNTP conc	dNTP Volume	Rxns per kit	
		Cat. no. 10297-018	Cat. no. 10297-117
0.2 mM	0.05 μL	5,000	50,000
0.5 mM	0.125 μL	2,000	20,000
1.0 mM	0.25 μL	1,000	10,000
1.5 mM	0.375 μL	666	6,666

Intended Use: For research use only. Not for human or animal therapeutic or diagnostic use.

Volumes and Concentrations, Continued

50 μ L Final Reaction Volume

Final dNTP conc	dNTP Volume	Rxns per kit	
		Cat. no. 10297-018	Cat. no. 10297-117
0.2 mM	0.1 μ L	2,500	25,000
0.5 mM	0.25 μ L	1,000	10,000
1.0 mM	0.5 μ L	500	5,000
1.5 mM	0.75 μ L	333	3,333

100 μ L Final Reaction Volume

Final dNTP Conc	dNTP Volume	Rxns per kit	
		Cat. no. 10297-018	Cat. no. 10297-117
0.2 mM	0.2 μ L	1,250	12,500
0.5 mM	0.5 μ L	500	5,000
1.0 mM	1.0 μ L	250	2,500
1.5 mM	1.5 μ L	166	1,666

Certificate of Analysis

The Certificate of Analysis (CofA) is available at www.invitrogen.com/cofa, and it is searchable by product lot number, which is printed on each box.

Limited Use Label License No. 358: Research Use Only

The purchase of this product conveys to the purchaser the limited, non-transferable right to use the purchased amount of the product only to perform internal research for the sole benefit of the purchaser. No right to resell this product or any of its components is conveyed expressly, by implication, or by estoppel. This product is for internal research purposes only and is not for use in commercial applications of any kind, including, without limitation, quality control and commercial services such as reporting the results of purchaser's activities for a fee or other form of consideration. For information on obtaining additional rights, please contact outlicensing@lifetech.com or Out Licensing, Life Technologies, 5791 Van Allen Way, Carlsbad, California 92008.

©2011 Life Technologies Corporation. All rights reserved. The trademarks mentioned herein are the property of Life Technologies Corporation or their respective owners.

For support visit www.lifetechnologies.com/support or email techsupport@lifetech.com
www.lifetechnologies.com

