



092PR-04

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ technical@GBiosciences.com

A Geno Technology, Inc. (USA) brand name

dNTP Mix

10 & 25mM dNTP mixes of dATP, dCTP, dGTP & dTTP

(Cat. # 786-442, 786-443, 786-457, 786-458, 786-459)



think proteins! think G-Biosciences www.GBiosciences.com

INTRODUCTION

The dNTP Mix is a ready-to-use aqueous solution at pH7.0 containing dATP, dCTP, dGTP and dTTP, each at a final concentration of 10 or 25mM. It has been functionally tested in PCR with Taq and Pfu DNA polymerases.

The mix is Endo-, exo-ribonuclease, ribonuclease and phosphatase free.

APPLICATIONS

Not for clinical use. For research use only in PCR, long PCR, RT-PCR, cDNA synthesis, primer extension, DNA sequencing and DNA labeling.

ITEMS INCLUDED

Cat. #	Description	μmoles supplied	Size
786-443	dNTP Mix [10mM]	5	0.5ml
786-442	dNTP Mix [10mM]	25	5 x 0.5ml
786-457	dNTP Mix [25mM]	25	1ml
786-458	dNTP Mix [25mM]	125	5ml
786-459	dNTP Mix [25mM]	250	10ml

STORAGE CONDITION

Shipped at ambient temperature. Upon arrival, store at -20°C.

TECHNICAL DETAILS

- -dATP: $C_{10}H_{13}N_5O_{12}P_3Na_3$; MW=557.2; $\lambda_{max}=259nm$; $\epsilon=15.2 \times 10^3 M^{-1}cm^{-1}$ at pH 7.0
- -dCTP: $C_9H_{13}N_3O_{13}P_3Na_3$; MW=533.1; $\lambda_{max}=271nm$; $\epsilon=9.3 \times 10^3 M^{-1}cm^{-1}$ at pH 7.0
- -dGTP: $C_{10}H_{13}N_5O_{13}P_3Na_3$; MW=573.2; $\lambda_{max}=253nm$; $\epsilon=13.7 \times 10^3 M^{-1}cm^{-1}$ at pH 7.0
- -dTTP: $C_{10}H_{14}N_2O_{14}P_3Na_3$; MW=548.1; $\lambda_{max}=267nm$; $\epsilon=9.6 \times 10^3 M^{-1}cm^{-1}$ at pH 7.0

VOLUMES AND CONCENTRATIONS

20µl Final Reaction Volume

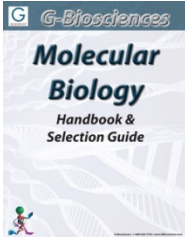
Final dNTP Concentration	10mM dNTP Volume	25mM dNTP Volume
0.2mM	0.4µl	0.16µl
0.5mM	1µl	0.4µl
1.0mM	2µl	0.8µl
1.5mM	3µl	1.2µl

25µl Final Reaction Volume

Final dNTP Concentration	10mM dNTP Volume	25mM dNTP Volume
0.2mM	0.5µl	0.22µl
0.5mM	1.25µl	0.55µl
1.0mM	2.5µl	1µl
1.5mM	3.75µl	1.5µl

RELATED PRODUCTS

Download our Molecular Biology Handbook.



<http://info.gbiosciences.com/complete-molecular-biology-handbook>

For other related products, visit our website at www.GBiosciences.com or contact us.

Last saved: 4/14/2014 CMH



www.GBiosciences.com