



090PR-03

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

A Geno Technology, Inc. (USA) brand name

dNTPs

(Cat. # 786-460, 786-465, 786-466, 786-467, 786-468)



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

INTRODUCTION

The dNTPs consists of aqueous solutions at pH7.0 of dATP, dCTP, dGTP or dTTP in 100mM concentration, each supplied individually. The dNTP set offers maximum flexibility in the preparation of reaction mixes for different applications, since the nucleotides are supplied separately.

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. #	dNTP	Concentration of each dNTP	Size	
786-460	dNTP Set (dATP, dCTP, dGTP, dTTP)	100mM	4 x 250µl	4 x 25µmol
786-465	dATP	100mM	250µl	25µmol
786-466	dCTP	100mM	250µl	25µmol
786-467	dGTP	100mM	250µl	25µmol
786-468	dTTP	100mM	250µl	25µmol

STORAGE CONDITION

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

TECHNICAL DETAILS

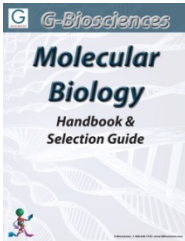
- -dATP: $C_{10}H_{13}N_5O_{12}P_3Na_3$; MW=557.2; $\lambda_{max}=259nm$; $\epsilon=15.2 \times 10^3 M^{-1}cm^{-1}$ at pH 7.0
- -dCTP: $C_9H_{13}N_3O_{13}P_3Na_3$; MW=533.1; $\lambda_{max}=271nm$; $\epsilon=9.3 \times 10^3 M^{-1}cm^{-1}$ at pH 7.0
- -dGTP: $C_{10}H_{13}N_5O_{13}P_3Na_3$; MW=573.2; $\lambda_{max}=253nm$;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

PREPARATION OF dNTP MIXTURE

dNTP Mixture to be prepared	Volumes of dNTP Set(μ l)				Water, nuclease-free(μ l)	Total Volume of dNTP Mixture(μ l)
	100mM dATP	100mM dGTP	100mM dCTP	100mM dTTP		
2mM of each dNTP	10	10	10	10	460	500
	100	100	100	100	4,600	5,000
	250	250	250	250	11,500	12,500
10mM of each dNTP	10	10	10	10	60	100
	100	100	100	100	600	1,000
	250	250	250	250	1,500	2,500
25mM of each dNTP	10	10	10	10	-	40
	100	100	100	100	-	400
	250	250	250	250	-	1,000

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