



092PR-04

G-Biosciences ♦ 1-800-628-7730 ♦ 1-314-991-6034 ♦ [technical@GBiosciences.com](mailto: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](http://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*

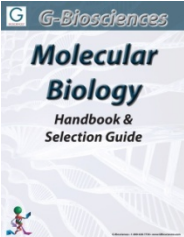
| Final dNTP Concentration | 10mM dNTP Volume | 25mM dNTP Volume |
|--------------------------|------------------|------------------|
| 0.2mM                    | 0.4µl            | 0.16µl           |
| 0.5mM                    | 1µl              | 0.4µl            |
| 1.0mM                    | 2µl              | 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](http://www.GBiosciences.com) or contact us.

Last saved: 4/14/2014 CMH



[www.GBiosciences.com](http://www.GBiosciences.com)