

dNTP Master Mix

Technical Data Sheet
 Mix of each dATP, dCTP, dGTP, dTTP, 5mM solutions

References: NU-0010-10
 NU-0010-50
 NU-0010-100

Products and procedures described in this protocol are intended for Research Purposes Only.

Details

NU-0010-10 20 µmol dNTPs
 NU-0010-50 5 x 20 µmol dNTPs
 NU-0010-100 10 x 20 µmol dNTPs

Description

The 20mM dNTP Master Mix contains each dNTP (dA, dC, dG, dT) at a concentration of 5 mM. Each vial contains 20 µmol (1 x 1000µl) dNTP Mix i.e. 5 µmol of each. Ready for immediate use.

Package contents

Reagent	Volume	Description
dNTP Master Mix 20mM total	1 x 1000 µl	NU-0010-10 20 µmol dNTPs
5mM each dNTP	5 x 1000 µl	NU-0010-50 5 x 20 µmol dNTPs
	10 x 1000 µl	NU-0010-100 10x20 µmol dNTPs

Shipping conditions

Dry ice

Storage conditions and stability

Store the Master Mix at -20 °C or -70 °C, in a constant temperature freezer. Avoid multiple freezing/ thawing. For long-term usage, aliquoting is recommended.

Typical analysis

Lithium salts, >99% deoxynucleoside triphosphates (HPLC, area %), <1% deoxynucleoside monophosphates and deoxynucleoside diphosphates.

dNTP characteristics

Deoxynucleotide	Molecular Weight	Molar Extinction
dATP	535.2	1µmol = 15,3 (A ₂₆₀)
dGTP	533.2	1µmol = 13,7 (A ₂₅₂)
dCTP	511.1	1µmol = 9,6 (A ₂₇₂)
dTTP	505.9	1µmol = 8,4 (A ₂₆₀)

Reaction conditions

For direct use in DNA synthesis *in vitro*. Add the Master Mix directly into the reaction mixture. For primer extension technology, add 4µl of dNTP Mix in a total 100µl reaction volume to give a final concentration of 200µM of each dNTP.

Number of reactions: 250

This is a guide only, for long-range application, adjust accordingly.

Related Products:

Reagent	Pack. size	Reference
dNTP set	4 x 5 µmoles	NU-0020-10
dUTP/dNTP Mix	5mM/1.25ml	NU-0040-01
GoldStar DNA polymerase	500U	ME-0064-05
GoldStar Mix	200Rxn	PK0064-02

For any further information required please contact our Customer Help Desk:

For Europe:

E-mail: info@eurogentec.com

Tel: +32 4 372 76 65

For USA:

E-mail: info.usa@eurogentec.com

Tel: +1 858 793 26 61



Eurogentec Headquarters

LIEGE SCIENCE PARK • 4102 SERAING • Belgium • Tel: +32 4 372 76 65
 Fax: +32 4 372 75 00 • info@eurogentec.com • www.eurogentec.com

Eurogentec North America

3347 Industrial Ct • Suite A • San Diego, CA 92121 • USA • Tel: +1 858 793 26 61
 Fax: +1 858 793 26 66 • info.usa@eurogentec.com • www.eurogentec.com