

Biotin-dCTP-PCR

Non-fluorescently labeled propargylamino-dCTP

DNA labeling and modification

	Cat. No.	Amount	Conc.
	PP-324S-BIO	200 µl	1 mM
	PP-324L-BIO	1 ml	1 mM

For *in vitro* use only
Quality guaranteed for 12 months
Store at -20°C, avoid frequent thawing and freezing

Biotin-dCTP-PCR

5-propargylamino-2'-deoxy-cytidine-5'-triphosphate
labeled with Biotin, triethylammonium salt, pH 7.5

Purity
>95%

Description

Biotin-dCTP-PCR is recommended for direct enzymatic labeling of DNA. The labeled-dCTP is specially optimized for incorporation into DNA by a wide variety of DNA polymerases. Recommended labeling methods include PCR, DOP-PCR, Nick Translation, 3'-end labeling and RT-PCR.

In PCR labeling, repeated cycles of denaturation, annealing and extension allow the amplification of a specific DNA fragment. Extension of the annealed primers with Taq polymerase results in a duplication of the DNA fragment in each cycle. When dCTP is partially substituted by Biotin-dCTP a Biotin labeled double-stranded DNA is generated.

The resultant DNA is suited for a variety of hybridization experiments, including Southern and Northern blots, colony hybridizations and *in situ* hybridizations.

Recommended PCR assay

20 µl PCR labeling assay			
component	stock conc.	amount	final conc.
High Yield Buffer complete (Cat.-No. PCR-201)	10x	2 µl	1x
dATP	1 mM	2 µl	100 µM
dCTP	1 mM	1 µl ¹⁾	50 µM ¹⁾
dGTP	1 mM	2 µl	100 µM
dTTP	1 mM	2 µl	100 µM
Biotin-dCTP-PCR	1 mM	1 µl ¹⁾	50 µM ¹⁾
forward Primer	10 µM	1 µl	500 nM
reverse Primer	10 µM	1 µl	500 nM
Template DNA		0.1-10 ng	5-500 pg/µl
Taq Pol (Cat.-No. PCR-201)	5 units/µl	0.2 µl	1 unit/assay
PCR grade H ₂ O		fill up to 20 µl	

1) The optimal final concentration of the labeled nucleotide may vary depending on the application.

Biotin-dCTP-PCR

Non-fluorescently labeled propagylamino-dCTP

DNA labeling and modification

Recommended cycling conditions

Initial denaturation	94°C	2 min	1x
Denaturation	94°C	30 sec	25-30x
Annealing ¹⁾	50-60°C	30 sec	
Elongation ²⁾	72°C	1 min	
Final elongation	72°C	5 min	1x

- 1) The annealing temperature depends on the melting temperature of the primers used.
- 2) The elongation time depends on the length of the fragments to be amplified. A time of 2 min/kb is recommended.

For optimal amplification results and high incorporation rates an individual optimization of the recommended PCR assay and cycling conditions may be necessary for each new primer-template pair.

Related products

Kits for DNA labeling
Standard PCR / Thermophilic Polymerases
Deoxynucleotides (dNTPs)
Primers and Oligonucleotides
DNA Ladders

For detailed information please visit
www.jenabioscience.com/pcr