



## HighFidelity Cy5 PCR Labeling Kit

Preparation of Cy5-labeled DNA probes by PCR

Cat. No.	Amount
APP-101-CY5-S	10 reactions x 20 µl
APP-101-CY5-L	50 reactions x 20 µl

**For general laboratory use.**

**Shipping:** shipped on gel packs

**Storage Conditions:** store at -20 °C

**Additional Storage Conditions:** avoid freeze/thaw cycles, store dark

**Shelf Life:** 12 months

**Spectroscopic Properties:**  $\lambda_{\text{exc}}$  649 nm,  $\lambda_{\text{em}}$  670 nm,  
 $\epsilon$  250.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

**Description:**

HighFidelity Cy5 PCR Labeling Kit is designed to produce randomly Cy5-modified DNA probes by PCR. Such probes are ideally suited for Fluorescence *in situ* hybridization (FISH) and Northern Blot experiments. PCR-based labeling is superior to random-primed labeling with Klenow fragment if template amounts are limited or amplification of a specific DNA fragments is required. Amplification of probes up to 4kbp is feasible.

dUTP-Cy5 is efficiently incorporated into DNA as substitute for its natural counterpart dTTP using an optimized reaction buffer and a High Fidelity Polymerase blend consisting of *Taq* polymerase and a proofreading enzyme. 50 % dUTP-Cy5 substitution typically results in an optimal balance between reaction and labeling efficiency. Individual optimization of dUTP-Cy5/dTTP ratio however, can easily be achieved with the single nucleotide format.

The kit contains sufficient reagents for 10 labeling reactions (S-Pack) or 50 labeling reactions (L-Pack) of 20 µl each (50% dUTP-Cy5 substitution, 100 µM dATP/dGTP/dCTP, 50 µM dTTP, 50 µM dUTP-Cy5).

**Content:**

**High Fidelity Polymerase**

in storage buffer with 50% glycerol (v/v)

#APP-101-Cy5-S: 1x 40 µl (100 units, 2.5 units/µl)

#APP-101-Cy5-L: 2x 40 µl (2x 100 units, 2.5 units/µl)

**High Fidelity Labeling Buffer**

1x 500 µl (10x)

**dATP - Solution**

1x 20 µl (100 mM)

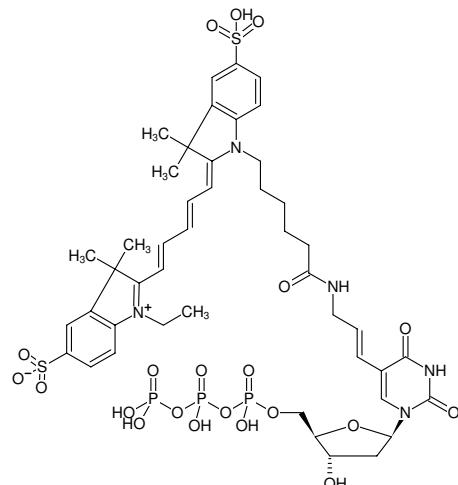
**dGTP - Solution**

1x 20 µl (100 mM)

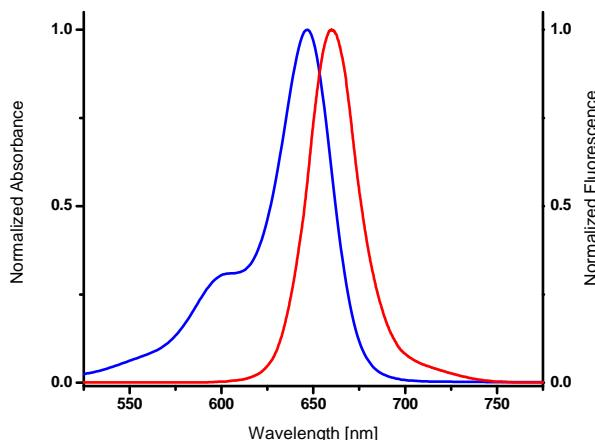
**dCTP - Solution**

1x 20 µl (100 mM)

**dTTP - Solution**



Structural formula of HighFidelity Cy5 PCR Labeling Kit



excitation and emission spectrum of Cy5



## HighFidelity Cy5 PCR Labeling Kit

Preparation of Cy5-labeled DNA probes by PCR

1x 20 µl (100 mM)

### dUTP-Cy5

#APP-101-Cy5-S: 1x 10 µl (1 mM)  
#APP-101-Cy5-L: 5x 10 µl (1 mM)

### Lambda DNA

1x 20 µl (100 ng/µl)

### 500 bp forward primer

1x 20 µl (10 µM)

### 500 bp reverse primer

1x 20 µl (10 µM)

### PCR-grade water

1x 1.2 ml

### To be provided by user

DNA template  
Primer  
DNA purification tools (optional)

## 1. Preparation of working solutions

### 1.1 Preparation of 1 mM dATP/dCTP/dGTP working solution

- Thaw 100 mM dATP, 100 mM dCTP and 100 mM dGTP solutions on ice, voretex and spin-down briefly.
- Prepare a 1:100 dilution with PCR-grade water to achieve a final concentration of 1 mM (e.g. 2 µl 100 mM dATP + 2 µl 100 mM dCTP + 2 µl 100 mM dGTP + 194 µl PCR-grade water).
- 1 mM ATP/CTP/GTP working solution can be stored at -20°C. Prepare aliquots to avoid freeze/thaw cycles.

### 1.2 Preparation of 1 mM dTTP working solution

- Thaw 100 mM dTTP solution on ice, voretex and spin-down briefly.
- Prepare a 1:100 dilution with PCR-grade water to achieve a final concentration of 1 mM (e.g. 2 µl 100 mM dTTP + 198 µl PCR-grade water).
- 1 mM dTTP working solution can be stored at -20 °C. Prepare aliquots to avoid freeze/thaw cycles.

## 3. Standard PCR Labeling protocol

The standard protocol is set-up for labeling of a 500 bp DNA fragment. An optimal balance between reaction and labeling efficiency is typically achieved with 50% dUTP-Cy5 substitution following the standard protocol below however, individual optimization might

improve results for individual applications.

- Assemble the PCR on ice in the order stated below (DNase-free reaction tube).
- Voretex and spin-down briefly.
- Perform assay set-up and reaction under low-light conditions.

Component	Volume	Final concentration
PCR-grade water	X µl	
High Fidelity Labeling Buffer (10x)	2 µl	1x
1 mM dATP/dCTP/dGTP working solution (s. 1.1)	2 µl	100 µM
1 mM dTTP working solution (s. 1.2)	1 µl	50 µM
1 mM dUTP-Cy5	1 µl	50 µM
forward primer (10 µM)	X µl	0.1 - 1 µM (e.g. 0.3 µM 500 bp forward primer)
reverse primer (10 µM)	X µl	0.1 - 1 µM (e.g. 0.3 µM 500 bp reverse primer)
template DNA	X µl	1 - 10 ng genomic DNA (e.g. 1 ng Lambda DNA)
High Fidelity Polymerase (2.5 units/µl)	1 µl	2.5 units
Total volume	20 µl	

## Recommended cycling conditions

Cycle step	Temperature	Time	Cycles
Initial denaturation	95°C	2 min	1x
Denaturation Annealing <sup>1)</sup> Elongation <sup>2)</sup>	95°C 58°C 68°C	20 sec 30 sec 60 sec	30x
Final Elongation	68°C	2 min	1x

<sup>1)</sup>The annealing temperature depends on the melting temperature of primers used.

<sup>2)</sup>The elongation time depends on the length of fragments to be amplified. A time of 2 min/kbp is recommended. Elongation at 72°C works as well.



## HighFidelity Cy5 PCR Labeling Kit

Preparation of Cy5-labeled DNA probes by PCR

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.

#### 4. Probe purification:

Probe purification is not required for most hybridization experiments. If a downstream application requires purification (e.g. concentration determination by absorbance measurement) we recommend silica-membrane or gel filtration-based purification.

#### Related Products:

Aminoallyl-dUTP-Cy5, #NU-803-Cy5