# **DATA SHEET**





# HighFidelity ATTO488 PCR Labeling Kit

Preparation of ATTO488-labeled DNA probes by PCR

Cat. No.	Amount
APP-101-488-S	10 reactions x 20 μl
APP-101-488-L	50 reactions x 20 μl



Structural formula of HighFidelity ATTO488 PCR Labeling Kit



excitation and emission spectrum of ATTO 488

#### 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:  $\lambda_{exc}$  500 nm,  $\lambda_{em}$  520 nm,  $\epsilon$  90.0 L mmol<sup>-1</sup> cm<sup>-1</sup> (Tris-HCl pH 7.5)

#### **Description:**

HighFidelity ATTO488 PCR Labeling Kit is designed to produce randomly ATTO488-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-XX-ATTO-488 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-XX-ATTO-488 substitution typically results in an optimal balance between reaction and labeling efficiency. Individual optimization of dUTP-XX-ATTO-488/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  $\mu$ l each (50% dUTP-XX-ATTO-488 substitution, 100  $\mu$ M dATP/dGTP/dCTP, 50  $\mu$ M dTTP, 50  $\mu$ M dUTP-XX-ATTO-488).

## Content:

#### **High Fidelity Polymerase**

in storage buffer with 50% glycerol (v/v) #APP-101-488-S: 1x 40 μl (100 units, 2.5 units/μl) #APP-101-488-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)



Jena Bioscience GmbH Löbstedter Str. 71 | 07749 Jena, Germany | Tel.:+49-3641-6285 000 https://www.jenabioscience.com







# HighFidelity ATTO488 PCR Labeling Kit

Preparation of ATTO488-labeled DNA probes by PCR

**dTTP - Solution** 1x 20 μl (100 mM)

### dUTP-XX-ATTO-488

#APP-101-488-S: 1x 10 μl (1 mM) #APP-101-488-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-XX-ATTO-488

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 concenctra- tion	
PCR-grade water	ΧμΙ		
High Fidelity La- beling Buffer (10x)	2 µl	1x	
1 mM dATP/dCTP/ dGTP working so- lution (s. 1.1)	2 µl	100 µM	
1 mM dTTP working solution (s. 1.2)	1 µl	50 μM	
1 mM dUTP-XX- ATTO-488	1 µl	50 µM	
forward primer (10 µM)	Χ μί	0.1 - 1 μM (e.g. 0.3 μM 500 bp forward primer)	
reverse primer (10 µM)	Χ μί	0.1 - 1 μM (e.g. 0.3 μM 500 bp reverse primer)	
template DNA	Χ μΙ	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



# **DATA SHEET**





# HighFidelity ATTO488 PCR Labeling Kit

Preparation of ATTO488-labeled DNA probes by PCR

amplified. A time of 2 min/kbp is recommended. Elongation at 72°C works as well.

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-XX-ATTO-488, #NU-803-XX-488

