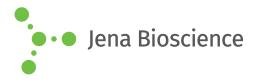
DATA SHEET

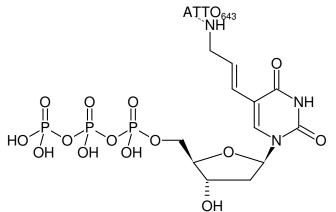




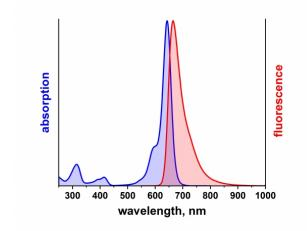
HighFidelity ATTO643 PCR Labeling Kit

Preparation of ATTO643-labeled DNA probes by PCR

| Cat. No. | Amount |
|---------------|----------------------|
| APP-101-643-S | 10 reactions x 20 μl |
| APP-101-643-L | 50 reactions x 20 μl |
| APP-101-643-L | 50 reactions x 20 μl |



Structural formula of HighFidelity ATTO643 PCR Labeling Kit



excitation and emission spectrum of ATTO 643

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: λ_{exc} 643 nm, λ_{em} 665 nm, ϵ 150.0 L mmol⁻¹ cm⁻¹ (Tris-HCl pH 7.5)

Description:

HighFidelity ATTO643 PCR Labeling Kit is designed to produce randomly ATTO643-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-ATTO-643 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-ATTO-643 substitution typically results in an optimal balance between reaction and labeling efficiency. Individual optimization of dUTP-ATTO-643/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 μ l each (50% dUTP-ATTO-643 substitution, 100 μ M dATP/dGTP/dCTP, 50 μ M dTTP, 50 μ M dUTP-ATTO-643).

Content:

High Fidelity Polymerase

in storage buffer with 50% glycerol (v/v) #APP-101-643-S: 1x 40 μl (100 units, 2.5 units/μl) #APP-101-643-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)

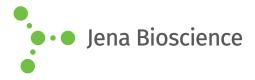
dGTP - Solution 1x 20 μl (100 mM)

dCTP - Solution 1x 20 μl (100 mM)



FIA AG Certified QMS and EMS according to DIN EN ISO 9001 and DIN EN ISO 14001 Ren –No. ICV03597 034 and ICV03597 534 lena Bioscience GmbH Löbstedter Str. 71 | 07749 Jena, Germany | Tel.:+49-3641-6285 000 https://www.jenabioscience.com







HighFidelity ATTO643 PCR Labeling Kit

Preparation of ATTO643-labeled DNA probes by PCR

dTTP - Solution 1x 20 μl (100 mM)

dUTP-ATTO-643

#APP-101-643-S: 1x 10 μl (1 mM) #APP-101-643-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-ATTO-643 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-ATTO- 643 | 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 μί | 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

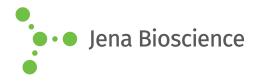
| Cycle step | Temperature | Time | Cycles |
|---|----------------------|----------------------------|--------|
| Initial denaturation | 95°C | 2 min | 1x |
| Denaturation Annealing ¹⁾ Elongation ²⁾ | 95°C 58°C 68°C | 20 sec 30 sec 60 sec | 30x |
| Final Elongation | 68°C | 2 min | 1x |

¹⁾The annealing temperature depends on the melting temperature of primers used.

 $^{(2)}$ The elongation time depends on the length of fragments to be



DATA SHEET





HighFidelity ATTO643 PCR Labeling Kit

Preparation of ATTO643-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-ATTO-643, #NU-803-643

