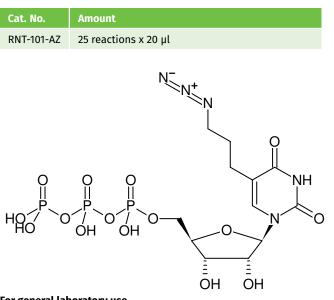




HighYield T7 Azide RNA Labeling Kit (UTP-based)

Preparation of randomly Azide-modified RNA probes by in vitro transcription with 5-Azido-C₃-UTP



For general laboratory use.

Shipping: shipped on gel packs

Storage Conditions: store at -20 °C

Additional Storage Conditions: avoid freeze/thaw cycles

Shelf Life: 12 months

Description:

HighYield T7 Azide RNA Labeling Kit (UTP-based) is designed to produce randomly Azide-modified RNA probes via *in vitro* transcription. Such probes are ideally suited for *in situ* hybridization, Northern Blot or crosslinking experiments.

5-Azido- C_3 -UTP is efficiently incorporated into RNA as substitute for its natural counterpart UTP using an optimized reaction buffer and T7 RNA Labeling Polymerase Mix. 35 % 5-Azido- C_3 -UTP substitution typically results in an optimal balance between reaction and labeling efficiency.

Individual optimization of 5-Azido-C₃-UTP/UTP ratio however, can easily be achieved with the single nucleotide format.

The resulting Azide-modified RNA probe can subsequently be labeled by copper-free Click Chemistry with Dibenzocyclooctyne (DBCO)-functionalized detection reagents (e.g DBCO-functionalized fluorescent dyes or (Desthio)Biotin) thereby forming a stable, covalently labeled RNA probe^[1].

The kit contains sufficient reagents for 25 labeling reactions of 20 μ l each (35 % 5-Azido-C₃-UTP substitution, 1 mM ATP, GTP, CTP, 0.65 mM UTP, 0.35 mM 5-Azido-C₃-UTP).

Content:

HighYield T7 RNA Labeling Polymerase Mix $2x 40 \mu$ l, incl. RNase inhibitor and 50 % glycerol (v/v)

HighYield T7 Reaction Buffer

1x 200 µl (10x), HEPES-based

ATP - Solution 1x 100 μl (100 mM)

GTP - Solution 1x 100 µl (100 mM)

CTP - Solution 1x 100 µl (100 mM)

UTP - Solution

1x 100 µl (100 mM)

5-Azido-C₃-UTP 1x 20 μl (10 mM)

T7 G-initiating control template (1.4 kbp)

1x 10 μl (200 ng/μl), 1.4 kbp PCR fragment plus T7 class III phi6.5 promotor resulting in approx. 1400 nt RNA transcript

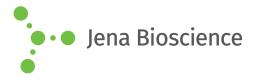
PCR-grade water

1x 1.2 ml



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







HighYield T7 Azide RNA Labeling Kit (UTP-based)

Preparation of randomly Azide-modified RNA probes by in vitro transcription with 5-Azido-C₃-UTP

DTT 1x 150 μl (100 mM)

To be provided by user

T7 Promotor-containing DNA template DBCO-functionalized detection molecule RNA purification tools RNAse-free DNAse I (optional)

1. Important Notes (Read before starting)

1.1 Prevention of RNAse contamination

Although a potent RNase Inhibitor is included, creating a RNAse-free work environment and maintaining RNAse-free solutions is critical for performing successful in vitro transcription reactions. We therefore recommend

- to perform all reactions in sterile, RNAse-free tubes using sterile pipette tips.
- to wear gloves when handling samples containing RNA.
- to keep all components tightly sealed both during storage and reaction procedure.

1.2 Template requirements

• <u>Template type</u>: Linearized plasmid DNA or PCR products containing a double-stranded T7 class II phi2.5 or class III phi6.5 promotor region upstream of the target sequence. Transcription initiation from T7 class III promotor is generally more efficient than initiation from T7 class II promotor. Minimum T7 promotor sequences:

T7 class III phi6.5 promotor 5'-<u>TAATACGACTCACTATAGNN</u>...-3' Bold: First base incorporated into RNA, *NN*: ideally *CG*

or

T7 class II phi2.5 promotor 5'-<u>TAATACGACTCACTATTAGNN</u>...-3' Bold: First base incorporated into RNA, *NN*: ideally *CG*

Template quality: DNA template quality directly influences yield and quality of transcription reaction. Linearized plasmid DNA needs to be fully digested and to be free of contaminating RNase, protein and salts. We recommend selecting restriction enzymes that generate blunt ends or 5'-overhangs and purification by phenol/chloroform extraction. A PCR mixture can be used directly however, better yields will usually be obtained with purified PCR products (e.g. via silica-membrane based purification columns).

2. Preparation of working solutions

2.1 Preparation of 10 mM ATP/CTP/GTP working solution

- Thaw 100 mM ATP, 100 mM CTP and 100 mM GTP solutions on ice, voretex and spin-down briefly.
- Prepare a 1:10 dilution with PCR-grade water to achieve a final concentration of 10 mM (e.g. 5 µl 100 mM ATP + 5 µl 100 mM CTP + 5 µl 100 mM GTP + 35 µl PCR-grade water).
- 10 mM ATP/CTP/GTP working solution can be stored at -20°C. Avoid freeze/thaw cycles.

2.2 Preparation of 10 mM UTP working solution

- Thaw 100 mM UTP solution on ice, voretex and spin-down briefly.
- Prepare a 1:10 dilution with PCR-grade water to achieve a final concentration of 10 mM (e.g. 5 μl 100 mM UTP + 45 μl PCR-grade water).
- 10 mM UTP working solution can be stored at -20 °C. Avoid freeze/thaw cycles.

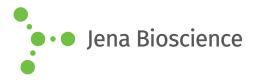
3. In vitro Transcription protocol

The protocol is optimized for 0.5 μ g - 1 μ g DNA template. An optimal balance between reaction and labeling efficiency is typically achieved with 35% 5-Azido-C₃-UTP substitution following the standard protocol below however, individual optimization might improve results for individual applications.

- Place HighYield T7 RNA Labeling Polymerase Mix on ice.
- Thaw all remaining components at room temperature (RT), mix by voretexing and spin down briefly.
- Assemble all components at RT to a nuclease-free microtube (sterile pipette tips) in the following order:
- Mix PCR-grade water, HighYield T7 Reaction Buffer and DTT by voretexing and spin down briefly.
- Add nucleotide solutions and template DNA, vortex and spin down briefly.
- Add HighYield T7 RNA Labeling Polymerase Mix vortex and spin down briefly.
- Incubate for 30 min at 37 °C in the dark (e.g. PCR cycler). Depending on the RNA probe individual optimization may increase product yield (2h 4h at 37 °C).









HighYield T7 Azide RNA Labeling Kit (UTP-based)

Preparation of randomly Azide-modified RNA probes by in vitro transcription with 5-Azido-C₃-UTP

Component	Volume	Final concenctra- tion
PCR-grade water	Xμl	
HighYield T7 Reac- tion Buffer (10x)	2 µl	1x
100 mM DTT	2 µl	
10 mM ATP/CTP/ GTP working solu- tion (s. 2.1)	2 μl	1 mM
10 mM UTP working solution (s. 2.2)	1.3 µl	0.65 mM
10 mM 5-Azido-C ₃ - UTP	0.7 μl	0.35 mM
Template DNA	Χ μl	0.5 - 1 µg
HighYield T7 RNA Labeling Poly- merase Mix	2 μl	
Total volume	20 µl	

Please note: Reagents for the following steps are not provided within this kit.

DNA template removal (optional)

Depending on the down-stream application of labeled RNA probe, removal of template DNA might be required. We recommend a salt-resistant, high efficiency DNAase such as Turbo™DNAse (ThermoFisher). Follow the manufacturer instructions.

General protocol for Copper-free CLICK labeling of Azide-modified RNA

For a detailled protocol, please refer to our background information: https://www.jenabioscience.com/images/741d0cd7d0/Click_Labeling_Azide_RNA.pdf

Related Products:

5-Azido-C3-UTP, #NU-157

Selected References:

[1] Miller et al. (2018) Near-infrared fluorescent northern blot. RNA 24(12):1871.

