

HighYield T7 Atto488 RNA Labeling Kit, S pack

Preparation of randomly Atto488-modified RNA probes by in vitro transcription with UTP-ATTO-488

HighYield T7 Atto488 RNA Labeling Kit is designed to produce randomly Atto488-modified RNA probes via in vitro transcription. Such probes are ideally suited for in situ hybridization and Northern Blot experiments. UTP-ATTO-488 is efficiently incorporated into RNA as substitute for its natural counterpart UTP using an optimized reaction buffer and T7 RNA Labeling Polymerase Mix. ATTO488 possesses, in contrast to Fluorescein, a excellent photostability. 35 % UTP-ATTO-488 substitution typically results in an optimal balance between reaction and labeling efficiency. Individual optimization of UTP-ATTO-488/UTP ratio however, can easily be achieved with the single nucleotide format. The resulting Atto488-modified RNA probe can subsequently be detected by fluorescence spectroscopy. The kit contains sufficient reagents for 20 labeling reactions of 20 μ l each (35 % UTP-ATTO-488 substitution, 2.5 mM ATP, GTP, CTP, 0.2 mM UTP, 0.1 mM UTP-ATTO-488).

Ordering Information

[Order Online »](#)

JBS-RNT-101-488-S	20Reactions
-------------------	-------------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Technical Info / Product For the Original Manufacturer's data sheet please [click here](#).

Notes



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)