

HighYield T7 P&L RNA NMR Kit (2F-ATP), L pack

Synthesis of 2-Fluoro-modified RNA

HighYield T7 P&L RNA NMR Kit (2F-ATP) is designed to produce large amounts of 2-Fluoro-modified RNA via in vitro transcription with T7 RNA polymerase. The resulting internally modified (m)RNA can subsequently be used for structure determination via nuclear magnetic resonance (NMR) spectroscopy. 2F-ATP is incorporated instead of its natural counterpart in a non-perturbing way (intact base pairing properties). 2-Fluoro-modification simultaneously functions as a sensitive NMR reporter group (larger chemical shift dispersion than 1H). The modified T7 RNA polymerase (T7 P&L RNA polymerase) with proline 266 replaced by leucine (P266L) has been associated with increased 5' homogeneity of transcripts especially when synthesized from A-initiating phi2.5 promoter. The kit contains sufficient reagents for 15 reactions (S-Pack) or 50 reactions (L-Pack) á 20 µl (7.5 mM GTP, 7.5 mM 2-Fluoro-ATP, 7.5 mM UTP, 7.5 mM CTP). An individual optimization of 2-Fluoro-ATP concentration can easily be achieved with the single nucleotide format. A 20 µl reaction yields about 80 µg RNA after 30 min incubation (1 µg T7 control template, 1.4 kb RNA transcript). Yields may however vary depending on the template (promotor design, sequence length, secondary structure formation).

Ordering Information

[Order Online »](#)

JBS-RNT-203-L	50Reactions
---------------	-------------

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)