

5-Ethynyl-uridine (5-EU)

Ordering Information

[Order Online »](#)

JBS-CLK-N002-10	10mg
-----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Handling Protect from moisture. Keep under inert gas.

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Appearance white to off-white solid

Application Notes RNA synthesis monitoring

CAS 69075-42-9

Couple Type Modified nucleotides

Formula $C_{11}H_{12}N_2O_6$

MW 268.22

Purity ≥ 99 % (HPLC)

Solubility Up to 200 mM in 1 x PBS or water by heating the obtained suspension for 1 minute by 70 °C, less soluble in DMSO

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



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)