

5-Ethynyl-cytidine (5-EC)

Ethynyl-labeled cytidine (5-EC) can be used as a replacement for BrU (5-Bromo-uridine) to measure de novo RNA synthesis in proliferating cells. 5-EC is cell permeable and incorporates into nascent RNA instead of its natural analog cytidine. The resulting ethynyl-functionalized RNA can subsequently be detected via Cu(I)-catalyzed click chemistry that offers the choice to introduce a Biotin group (via Azides of Biotin) for subsequent purification tasks or a fluorescent group (via Azides of fluorescent dyes) for subsequent microscopic imaging.

Ordering Information

[Order Online »](#)

JBS-CLK-087	10mg
-------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

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 65223-78-1

Formula $C_{11}H_{13}N_3O_5$

MW 267.24

Purity $\geq 99\%$ (HPLC)

Solubility 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

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

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

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

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