

2'-CMP

Metabolite formed during enzymatic hydrolysis of cytidine-2',3'-cyclic monophosphate by 2',3'-cyclic nucleotide 3'-phosphodiesterase

BLG-C093-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

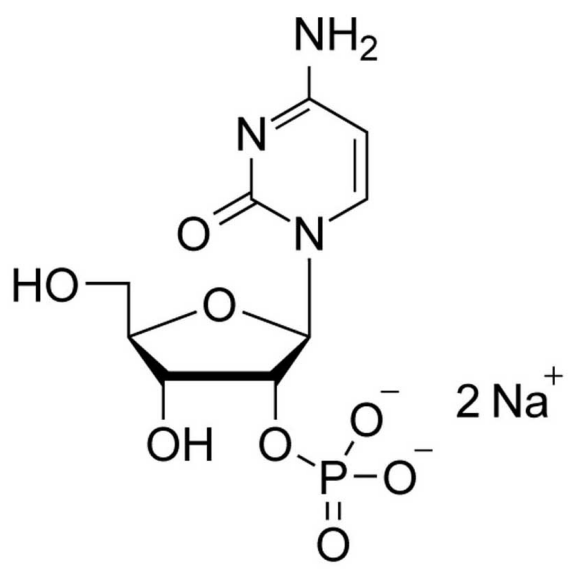
Ordering Information

[Order Online »](#)

BLG-C093-05	5µmol
BLG-C093-25	5x5µmol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Cytidine-2'-O-monophosphate . sodium salt

CAS 85-94-9

Couple Type Modified nucleotides

Formula $C_9H_{14}N_3O_8P$ (free acid)

MW 323.2

Purity > 95% HPLC

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