

2',3'-cIMP

2',3'-cIMP is structurally related to the naturally occurring 2',3'-cNMPs 2',3'-cAMP (Cat. No. A 307) and 2',3'-cGMP (Cat. No. G 025). Although 2',3'-cIMP was detected in a number of mammalian organs by Jia et al. (2014), its biological role is still elusive.

*BLG-I032-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

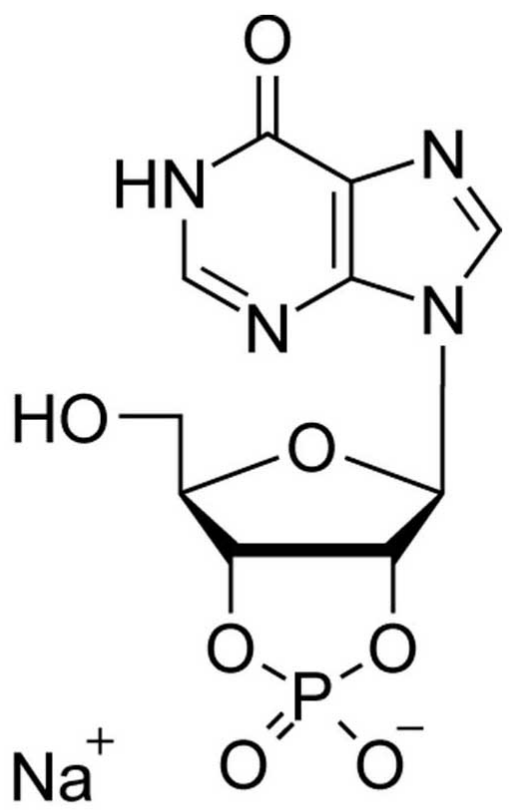
Ordering Information

[Order Online »](#)

BLG-I032-25	5x5 μ mol
-------------	---------------

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 Inosine-2',3'-cyclic monophosphate . sodium salt

CAS 15569-30-9

Couple Type Modified nucleotides

Formula $C_{10}H_{10}N_4O_7P \cdot Na$

MW 352.18

Purity > 97% HPLC

Quantity 5µmol are equivalent to ~1.8 mg

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