

isoGDP

Analogue of GDP

Analogue of adenosine-5'-O-diphosphate which can be useful in receptor mapping studies. The corresponding triphosphate form is offered as well (isoGTP, Cat. No. BLG-I039).

BLG-I038-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

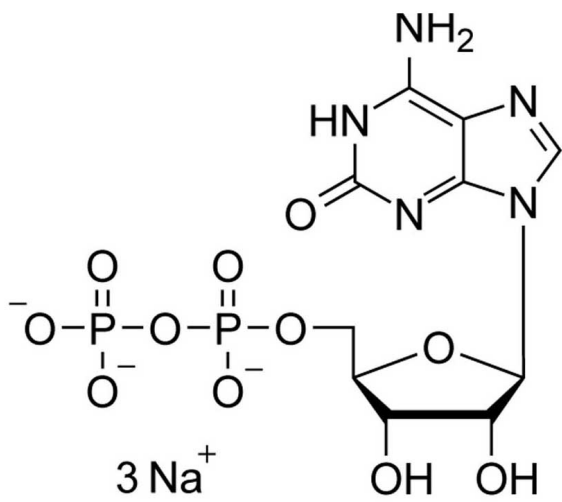
Ordering Information

[Order Online »](#)

BLG-I038-05	5x1 μ mol
-------------	---------------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Isoguanosine-5'-O-diphosphate . sodium salt

CAS 58083-94-6

Formula $C_{10}H_{15}N_5O_{11}P_2$ (free acid)

Formulation 10 mM aqueous solution

MW 443.2 (free acid)

Purity > 95% HPLC



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