

8-NO₂-cGMP

Second messenger of NO

The nitrated derivative of cGMP, 8-NO₂-cGMP (8-Nitro-cGMP), is considered to play an important role in NO-mediated signal transduction. Formation of protein Cys-cGMP adducts by 8-NO₂-cGMP was identified as a new post-translational modification of proteins, which is called S-guanylation.

BLG-N011-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

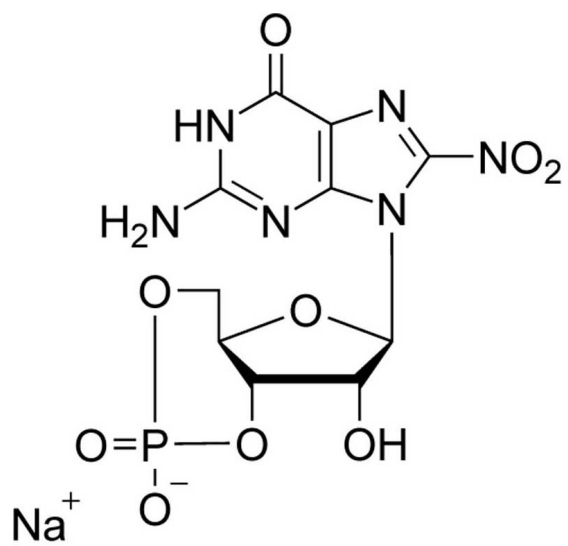
Ordering Information

[Order Online »](#)

BLG-N011-05	5µmol
BLG-N011-25	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 8-Nitroguanosine-3',5'-cyclic monophosphate . sodium salt

CAS 913645-39-3

Couple Type Modified nucleotides

Formula $C_{10}H_{10}N_6O_9P \cdot Na$

MW 412.19

Purity > 98% HPLC

Quantity 5µmol are equivalent to ~2.1 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