

8-NO₂-Guo

Reference for analysis of hydrolysates of RNA stressed by nitric oxide/superoxide and peroxyxynitrite, respectively. Corresponding nucleobase available (Prod. No. [BLG-N003](#)). Instable, can contain 8-nitroguanine.

BLG-N004-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

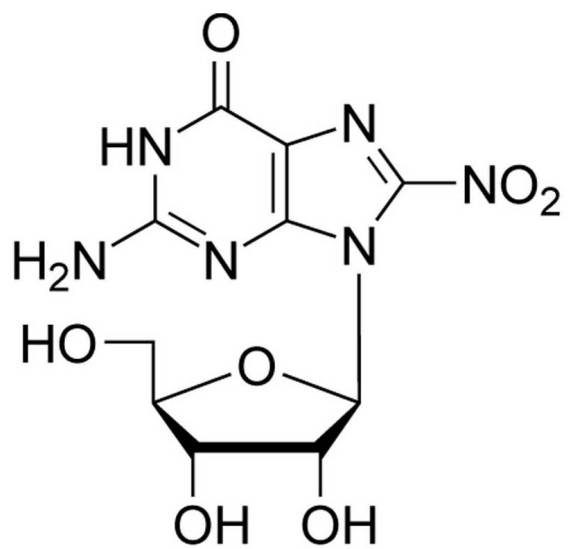
Ordering Information

[Order Online »](#)

BLG-N004-05	5µmol
BLG-N004-25	5x5µmol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	8-Nitroguanosine
CAS	337536-53-5
Couple Type	Modified nucleotides
Formula	$C_{10}H_{12}N_6O_7$
MW	328.3
Purity	> 95% HPLC
Quantity	5µmol≈1.6mg

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