

4-N₃-PET-8-Biotin-11-cGMP

Biotinylated cGMP analogue for photoaffinity labelling of cGMP-responsive receptors.

BLG-A186-005 (5 x 0.1 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

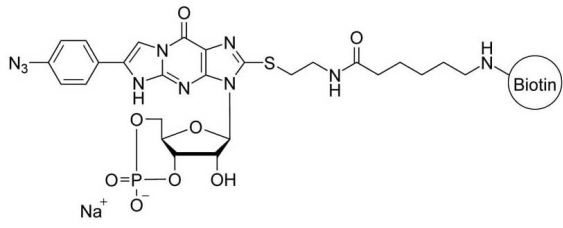
Ordering Information

[Order Online »](#)

BLG-A186-001	0.1µmol
BLG-A186-005	5x0.1µmol

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	β -(4-Azidophenyl)-1, N2-etheno-8-(2-(6-[biotinyl]aminohexanoyl)aminoethylthio)guanosine-3',5'-cyclic monophosphate . sodium salt
Formula	$C_{36}H_{44}N_{12}O_{10}PS_2 \cdot Na$
MW	922.9
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