

2'-[mPEG(11)]-AHC-c-diGMP

c-diGMP analogue

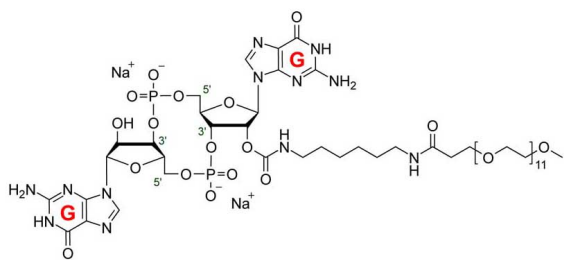
Analogue of the bacterial second messenger c-diGMP (Cat. No. BLG-C057), where one of the two ribose 2'-hydroxy groups are connected to a polyethylene glycol chain. Its increased lipophilicity could be advantageous for passing cellular membranes.

BLG-M073-005 (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-M073-001	0.1 µmol
BLG-M073-005	5x0.1 µmol

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 2'-O-(6-[mPEG(11)]amino)hexylcarbamoyl)-cyclic diguanosine monophosphate . sodium salt

Appearance Solid

Formula $C_{53}H_{88}N_{12}O_{28}P_2$ (free acid)

MW 1403.3 (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