

2'-[DY-547]-AHC-c-diGMP

Fluorescent analogue of c-diGMP

Fluorescent analogue of the bacterial second messenger c-diGMP (Cat. No. BLG-C057);

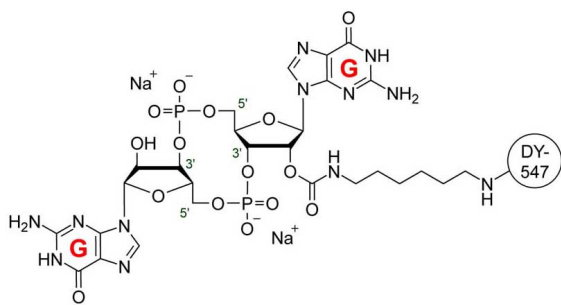
Ordering Information

[Order Online »](#)

BLG-D116-001	0.1µmol
--------------	---------

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	2'-O-(6-[DY-547]-aminohexylcarbamoyl)-cyclic diguanosine monophosphate . sodium salt
Couple Type	Modified nucleotides
Formula	$C_{57}H_{71}N_{14}O_{22}P_2S_2 \cdot Na$
MW	1453.3 (free acid nucleotide)
Purity	> 95% HPLC
Quantity	0.1 μmol ≈ 0.1 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