

2',2''-Di-F-c-didGMP

Analogue of the bacterial second messenger
c-diGMP

Analogue of the bacterial second messenger and STING activator c-diGMP (Cat. No. BLG-C057) in which the ribose 2'-hydroxy groups have been replaced by fluorine atoms. 2',2''-Di-F-c-didGMP may be useful in studies on ligand-receptor interactions.

BLG-D210-005 (5 x 0.1 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

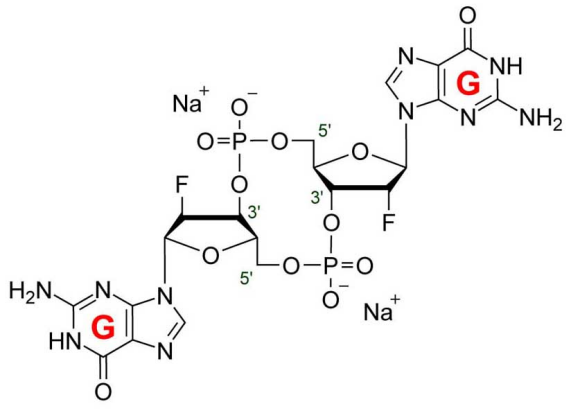
Ordering Information

[Order Online »](#)

BLG-D210-001	0.1µmol
BLG-D210-005	5x0.1µmol

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'-,2''-Dideoxy-2'-,2''-difluoro-cyclic diguanosine monophosphate . sodium salt

Appearance Solid

CAS 1334145-18-4

Formula $C_{20}H_{22}F_2N_{10}O_{12}P_2$ (free acid)

MW 694.4 (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