

c[A(2',5')pS[Rp]- A(3',5')pS[Rp]]

BLG-C224-005 (5 x 0.1 μ mol pack size) is not sold in the U.S. or Canada.

Please [contact us](#) for available options.

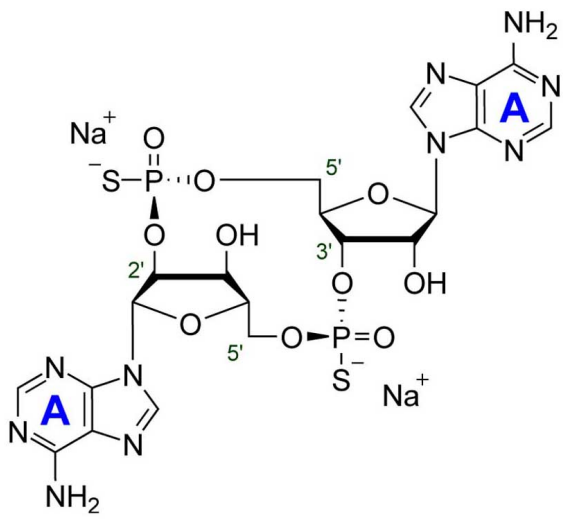
Ordering Information

[Order Online »](#)

BLG-C224-005	5x0.1 μ mol
--------------	-----------------

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 Cyclic (adenosine-(2'→5')-monophosphorothioate[Rp]-adenosine-(3'→5')-monophosphorothioate[Rp]) . sodium salt, Rp,Rp-2'3'-c-diAMPSS

CAS 1638750-95-4

Formula $C_{20}H_{24}N_{10}O_{10}P_2S_2$ (free acid)

MW 690.6

Purity > 95% HPLC

Quantity 0.1umol=~69ug

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