

c[A(2',5')pA(2',5')p]

BLG-C188-025 (5 x 0.5 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.

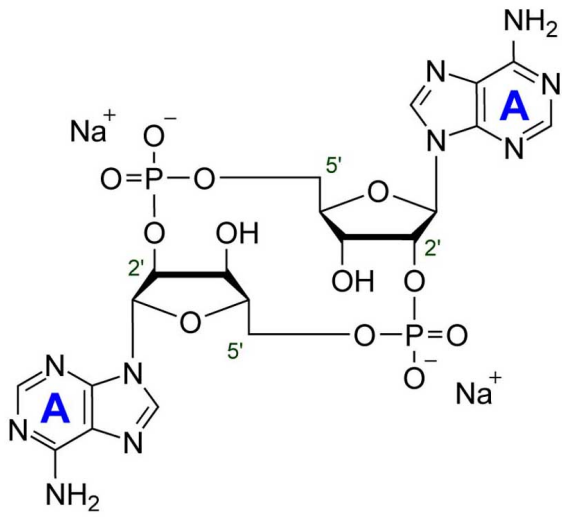
Ordering Information

[Order Online »](#)

BLG-C188-005	0.5 μ mol
BLG-C188-025	5x0.5 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')-monophosphate-adenosine-(2'→5')-monophosphate) . sodium salt, 2'2'-c-diAMP, 2',5'-2',5'-c-diAMP

CAS 1427449-35-1

Formula $C_{20}H_{24}N_{10}O_{12}P_2$ (free acid)

MW 658.4

Purity purity > 95% HPLC

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