

A(2',5')pA

Potential metabolite of 2' → 5'-linked oligoadenylates (2-5As)

A(2',5')pA is a potential metabolite of 2' → 5'-linked oligoadenylates (2-5As). A(2',5')pA does not activate RNase L.

BLG-A352-05 (5 x 1 μmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

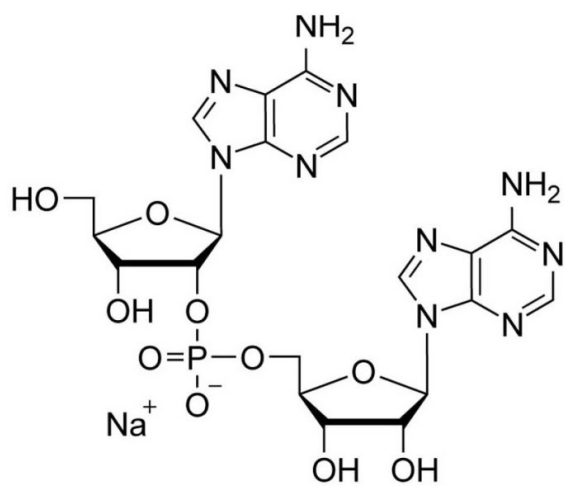
Ordering Information

[Order Online »](#)

BLG-A352-01	1μmol
BLG-A352-05	5x1μmol

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 Adenosine-5'-O-(α , β -methylene)diphosphoribose . sodium salt

Appearance Solid

CAS 2273-76-9

Couple Type Modified nucleotides

Formula $C_{20}H_{25}N_{10}O_{10}P$ (free acid)

MW 596.5

Purity > 95% HPLC

Quantity 1 μ mol is equivalent to ~0.6 mg



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)