

pA-3'-p / pAp

Used to study the kinetic properties and structure of sulfotransferases

In pA-3'-p phosphate groups are attached to both the ribose 3' position and the ribose 5' position of adenosine. This ubiquitous nucleotide is present in almost all organisms. It is a product of 3'-phosphoadenosine 5'-phosphosulfate (PAPS), a cofactor of sulfotransferases (ST / SULTs), and can be used to study the kinetic properties and structure of sulfotransferases. pA-3'-p can also act as an inhibitor of exoribonucleases, such as Xrn1, Rat1p and DXO, and of poly(ADP-ribose) polymerase 1 (PARP-1). Furthermore, it inhibits ADP-induced platelet shape change and aggregation but does not affect ADP-induced inhibition of adenylyl cyclase.

BLG-A444-50 (5 x 10 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

Ordering Information

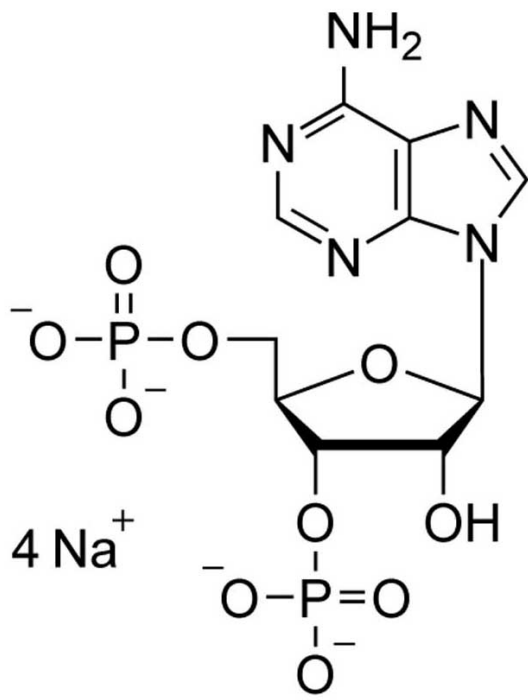
[Order Online »](#)

BLG-A444-50

5x10µmol

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-3', 5'-O-bisphosphate . sodium salt

CAS 75431-54-8

Couple Type Modified nucleotides

Formula $C_{10}H_{15}N_5O_{10}P_2$ (free acid)

MW 427.2 (free acid)

Purity > 95% HPLC

Quantity 10µmol = ~4.3mg

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