

2'/3'-O-Ac-ADPR / ADP-2'/3'-O-Ac-R

Analogue of ADPR and putative second messenger for characterization of receptor proteins

2'/3'-O-Ac-ADPR is an analogue of ADPR, that is generated from NAD⁺ by sirtuin-mediated protein deacetylation. 2'/3'-O-Ac-ADPR is a putative intracellular second messenger and has been reported to play a role in starfish oocyte maturation, to bind to a number of intracellular proteins, including heterochromatic proteins linked to X-chromosome inactivation and to activate the TRPM2 ion channel.

BLG-A405-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

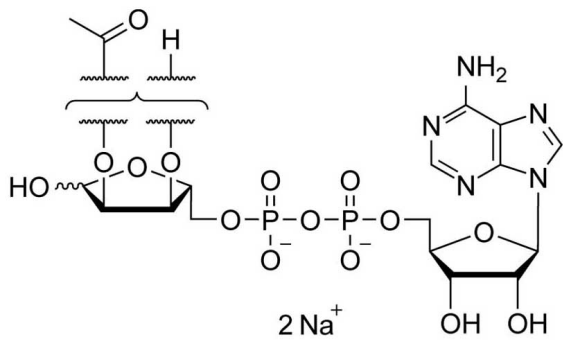
Ordering Information

[Order Online »](#)

BLG-A405-05	5µmol
BLG-A405-25	5x5µmol

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'-/ 3'-O-Acetyladenosine-5'-O-diphosphoribose, Adenosine-5'-O-diphospho-2'-/ 3'-O-acetylribose . sodium salt

Appearance Solid

CAS 642090-67-3

Couple Type Modified nucleotides

Formula $C_{17}H_{25}N_5O_{15}P_2$ (free acid)

MW 601.4 (free acid)

Purity > 95% HPLC

Quantity 5µmol are equivalent to ~3 mg

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