

AMPCPR

**Analogue of ADPR with improved stability
well suited for protein crystallization studies
and characterization of receptor proteins**

AMPCPR is an ADPR mimic with improved hydrolytic stability, suitable for the characterization and crystallization of NUDIX hydrolases. NUDIX hydrolases are a superfamily of hydrolytic enzymes capable of cleaving nucleoside diphosphates linked to another moiety X. Within the superfamily, several categories are classified based on their substrate specificities, like the MutT family hydrolases, the ADPR pyrophosphatases, the NADH hydrolases, and the ApnA hydrolases. AMPCPR was reported to interact with the nudix-type motif 9 (NUDT9)-homolog domain of transient receptor potential melastatin 2 (TRPM2), a frequently studied cation channel involved in multiple physiologic and pathophysiologic processes.

*BLG-A290-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

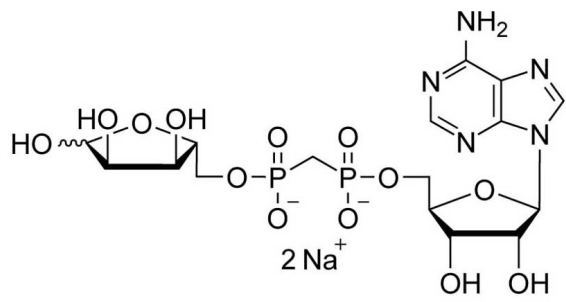
Ordering Information

[Order Online »](#)

BLG-A290-01	1 μ mol
BLG-A290-05	5x1 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 Adenosine-5'-O-(α , β -methylene)diphosphoribose . sodium salt

Appearance Solid

CAS 182950-50-1

Couple Type Modified nucleotides

Formula $C_{16}H_{25}N_5O_{13}P_2$ (free acid)

MW 557.3

Purity > 95% HPLC

Quantity 1 μ mol is equivalent to ~0.6 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