

2'-pRib-AMP

Isomeric analogue of the natural signal molecule ADPR

BLG-P128-025 (5 x 0.5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

Ordering Information

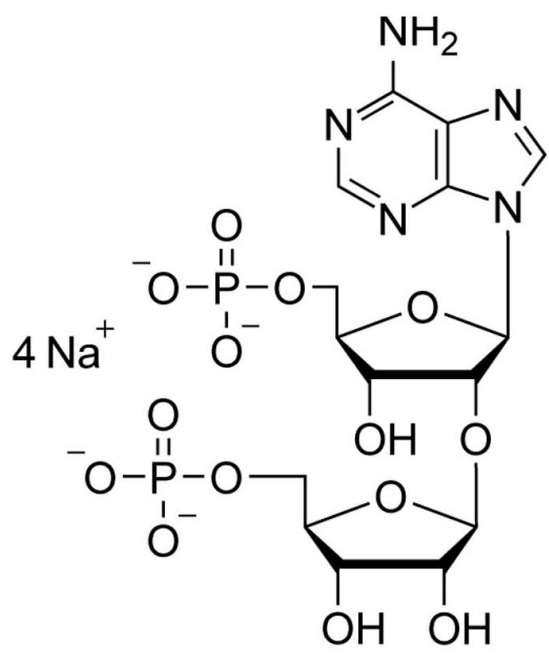
[Order Online »](#)

BLG-P128-005	0.5µmol
BLG-P128-025	5x0.5µmol

Manuals, SDS & CofA

[View Online »](#)

2'-pRib-AMP is the monomeric metabolite of poly(ADP-ribose) (PAR) and can be used to detect and analyse PAR-specific binding properties of poly(ADP-ribose) polymerases. Recent studies have shown that plant nucleotide-binding leucine-rich repeat (NLR) receptors with an N-terminal Toll/interleukin-1 receptor (TIR) domain recognise pathogen effectors. The TIR-enabled nicotinamideadenine dinucleotide hydrolase (NADase) activity is used to mediate the immune response. It has been shown that TIR-containing proteins can biocatalyse the production of 2'-pRib-AMP from NAD⁺ in plants, and it has been postulated that 2'-pRib-AMP may have a second messenger function in TIR signalling of plant immunity.



Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 2'-(5"-Phosphoribosyl)-adenosine-5'-O-monophosphate . sodium salt, 2'-pRib-5'-AMP, iso-ADPR

CAS 1672716-76-5

Formula $C_{15}H_{23}N_5O_{14}P_2$ (free acid)

MW 559.3 (free acid)

Purity > 95% HPLC

Quantity 0.5 µmol = ~280 µg

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