

# ADP-HPD

**Analogue of ADPR and highly potent, noncompetitive and specific inhibitor of PARG**

ADP-HPD is a highly potent, noncompetitive and specific inhibitor of poly-(ADP-ribose) glycohydrolase (PARG; IC<sub>50</sub> = 120 nM), displaying a 1000-fold lower IC<sub>50</sub> compared to native hydrolase product ADPR (IC<sub>50</sub> = 120 μM). The NAD glycohydrolase from Bungarus fuscus venom is less sensitive to inhibition by ADP-HPD, exhibiting an IC<sub>50</sub> = 260 μM. ADP-HPD does not inhibit either poly(ADP-ribose) polymerase (PARP) or NAD:arginine mono(ADP-ribosyl)-transferase A at inhibitor concentrations up to 1 mM. ADP-HPD has been widely used in PAR research, besides numerous other applications, for example in crystallization experiments with PARG, to investigate the structure and function of PARG, and to measure PARP activity in permeabilized cells.

*BLG-A471-025 (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-A471-005	0.5μmol
BLG-A471-025	5x0.5μmol

## 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-diphosphate-(hydroxymethyl)-pyrrolidinediol . sodium salt

**Appearance** Solid

**CAS** 148252-73-7

**Couple Type** Modified nucleotides

**Formula**  $C_{15}H_{24}N_6O_{12}P_2$

**MW** 542.3

**Purity** > 98 % HPLC

**Quantity** 0.5 µmol are equivalent to ~0.3 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](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)