

# 8-PT-ADPR

Analogue of the second messenger ADP-ribose (ADPR)

8-PT-ADPR is a putative antagonist of the second messenger ADP-ribose (ADPR), inhibiting ADPR-activated cation transport through TRPM2.

*BLG-P079-05 (5 x 1  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

## Ordering Information

[Order Online »](#)

BLG-P079-01	1 $\mu$ mol
BLG-P079-05	5x1 $\mu$ 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** 8- Phenylthioadenosine- 5'- O- diphosphoribose . sodium salt

**Appearance** Solid

**Couple Type** Modified nucleotides

**Formula**  $C_{21}H_{27}N_5O_{14}P_2S$  (free acid)

**MW** 667.5 (free acid)

**Purity** > 95% HPLC

**Quantity** 1µmol is equivalent to ~0.7 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)