

# His-ADPR / Histidine-ADPR

## Analogue of ADPR identified as new second messenger in type II Thois defence system in bacteria

His-ADPR is a unique signaling molecule produced by the Toll/interleukin-1 receptor (TIR) domain protein of the type II Thois defense system in bacteria. It is formed in response to phage infection and activates the corresponding Thois effector by binding to a macro domain located at the C-terminus of the effector protein. His-ADPR is the first described member of a new class of TIR-derived immune signaling molecules that combine nucleotide and amino acid subunits.

*BLG-H042-025 (5 x 0.5  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

### Ordering Information

[Order Online »](#)

BLG-H042-005	0.5 $\mu$ mol
BLG-H042-025	5x0.5 $\mu$ 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** Adenosine-5'-O-diphospho-1"-β-(NE1(τ)-L-histidine)-ribose . sodium salt

**Appearance** Solid

**Couple Type** Modified nucleotides

**Formula**  $C_{21}H_{30}N_8O_{15}P_2$  (free acid)

**MW** 696.5 (free acid)

**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)