

# 2'-cADPR

Regioisomer of the second messenger cADPR and of 3'-cADPR

BLG-C406-025 (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

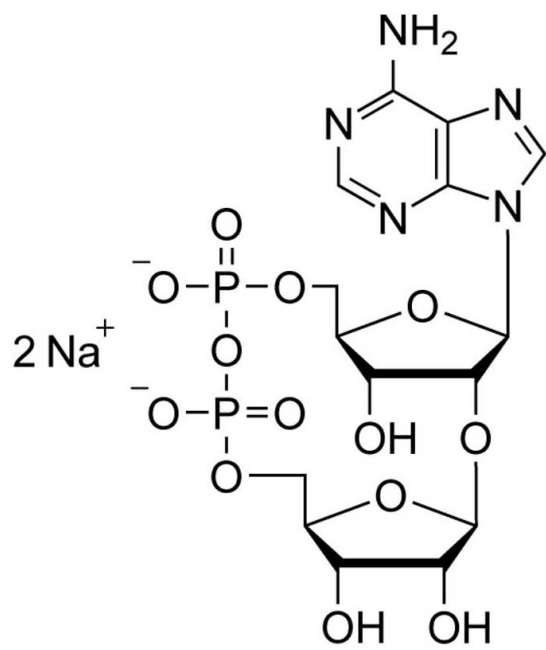
[Order Online »](#)

BLG-C406-005	0.5 $\mu$ mol
BLG-C406-025	5x0.5 $\mu$ mol

## Manuals, SDS & CofA

[View Online »](#)

2'-cADPR is a regioisomer of the canonical second messenger cADPR (Cat. No. BLG-C005) and of 3'-cADPR (Cat. No. BLG-C404). Biosynthesis of 2'-cADPR was identified in Toll/interleukin-1 receptor (TIR)-domain containing proteins in plants and bacteria, and it is hypothesized to have a role in immune signaling.



## Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** (1"-2")-Cyclic adenosine diphosphate ribose . sodium salt, v-cADPR, 1"-2'-gcADPR

**Formula**  $C_{15}H_{21}N_5O_{13}P_2$  (free acid)

**MW** 541.3 (free acid)

**Purity** > 98% HPLC

**Quantity** 0.5 µmol are equivalent to ~ 270 µg



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)