

# 8-Br-7-CH-cADPR

Membrane-permeant and hydrolysis-resistant cADPR antagonist

8-Br-7-CH-cADPR is a membrane-permeant and hydrolysis-resistant antagonist of the second messenger cADPR.

*BLG-B100-025 (5 x 0.5  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

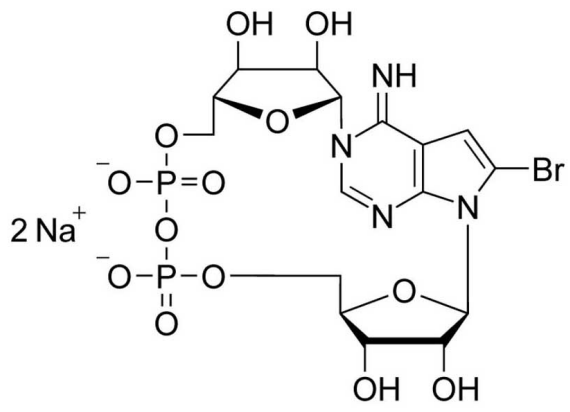
## Ordering Information

[Order Online »](#)

BLG-B100-005	0.5 $\mu$ mol
BLG-B100-025	5x0.5 $\mu$ mol

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 8-Bromo-7-deaza-cyclic adenosine diphosphate ribose . sodium salt, 7-Deaza-8-bromo-cyclic ADP-ribose

**CAS** 189876-06-0

**Couple Type** Modified nucleotides

**Formula**  $C_{16}H_{21}BrN_4O_{13}P_2$  (free acid)

**MW** 619.2 (free acid)

**Purity** > 97% HPLC

**Quantity** 0.5µmol≈310µg

**Solubility** Soluble in water (>50mM).

**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)