

# 8-MA-N1-cIDPR

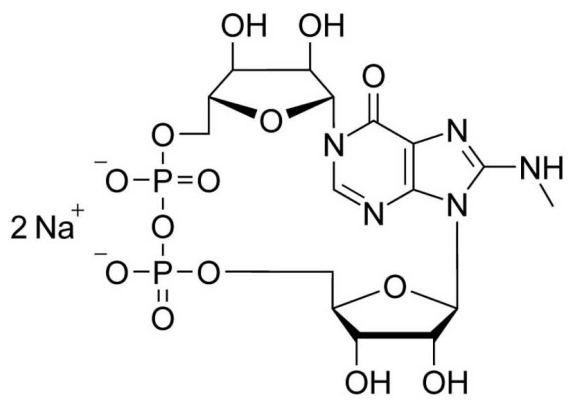
Hydrolysis-resistant putative agonist of the second messenger cyclic ADP ribose (Cat. No. C 005) with increased lipophilicity and membrane permeability. The parent compound N1-cIDPR (Cat. No. C 098) is offered as well.

*BLG-M063-025 (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 »](#)

<b>BLG-M063-005</b>	0.5 $\mu$ mol
<b>BLG-M063-025</b>	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** 8-Methylamino-cyclic inosine diphosphate ribose . sodium salt

**Formula**  $C_{16}H_{23}N_5O_{14}P_2$  (free acid)

**MW** 571.3

**Purity** > 97% HPLC



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)