

8-Br-N¹-cIDPR

Analogue of cADPR with higher permeability and stability

8-Br-N¹-cIDPR is a membrane-permeant and more stable analogue of the second messenger cADPR

BLG-B071-025 (5 x 0.5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

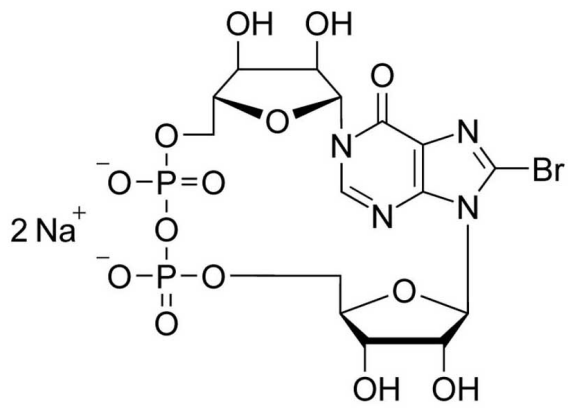
Ordering Information

[Order Online »](#)

BLG-B071-005	0.5µmol
BLG-B071-025	5x0.5µmol

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-Bromo-cyclic inosine diphosphate ribose . sodium salt

CAS 638195-70-7

Couple Type Modified nucleotides

Formula $C_{15}H_{19}BrN_4O_{14}P_2$ (free acid)

MW 621.2 (free acid)

Purity > 97% HPLC

Quantity 0.5µmol≈~311µg

Solubility Excellent solubility in water and aqueous buffers.

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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com