

8-Br-cADPR

Antagonist of the second messenger cyclic
ADP ribose

Analogue and antagonist of the second messenger cyclic adenosine diphosphate ribose (cADPR, [Prod. No. BLG-C005](#)) in which the hydrogen in position 8 of the heterocyclic nucleobase is replaced by bromine.

BLG-B065-025 (5 x 0.5 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

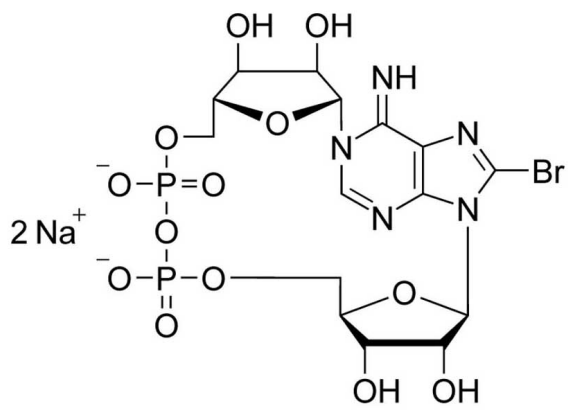
Ordering Information

[Order Online »](#)

BLG-B065-025	5x0.5 μ 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-cyclic adenosine diphosphate ribose . sodium salt

CAS 151898-26-9

Couple Type Modified nucleotides

Formula $C_{15}H_{20}BrN_5O_{13}P_2$ (for free acid)

MW 620.2

Purity > 95% HPLC

Quantity 0.5µmol≈~310µg

Solubility Soluble 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