

8-AET-N¹-cIDPR

1-cIDPR functionalized for conjugation with markers and dyes or for immobilization

8-AET-N¹-cIDPR is a metabolically stable analogue of cADPR which is suitable as a ligand in affinity chromatography or for coupling of various labelling structures including fluorophores. The parent compound N¹-cIDPR is offered as well (Prod. No. BLG-C098).

BLG-A140-025 (5 x 0.5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

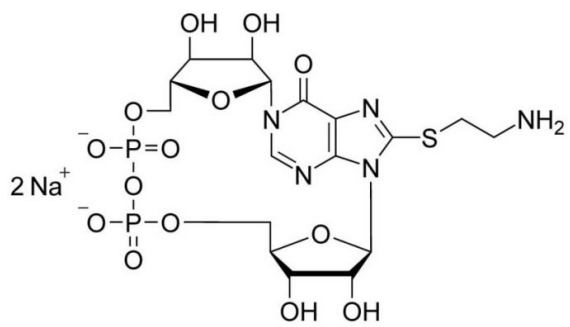
Ordering Information

[Order Online »](#)

BLG-A140-005	0.5µmol
BLG-A140-025	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-(2-Aminoethylthio)-cyclic inosine diphosphate ribose . sodium salt

Couple Type Modified nucleotides

Formula $C_{17}H_{25}N_5O_{14}P_2S$ (free acid)

MW 617.4 (free acid)

Purity > 97% HPLC

Quantity 0.5µmol≈309µg

Solubility Soluble in water or 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](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)