

8-PhEtT-N1-cIDPR

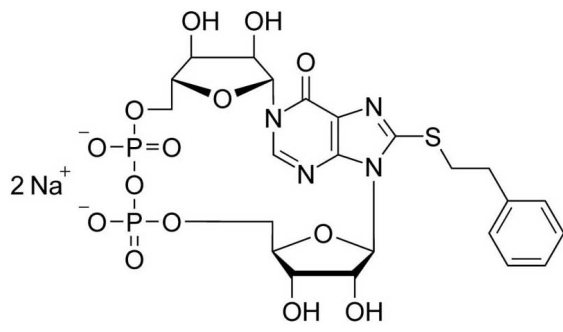
Hydrolysis-resistant putative agonist of the second messenger cyclic ADP ribose (Cat. No. C 005) with significantly increased lipophilicity and membrane permeability. The parent compound N1-cIDPR (Cat. No. C 098) is offered as well.

BLG-P034-025 (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-P034-005	0.5 μ mol
BLG-P034-025	5x0.5 μ 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-Phenylethylthio-cyclic inosine diphosphate ribose . sodium salt

Formula $C_{23}H_{28}N_4O_{14}P_2S$ (free acid)

MW 678.5

Purity > 97% HPLC



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