

3'-cADPR

Regioisomer of the second messenger
cADPR and of 2'-cADPR

*BLG-C404-025 (5 x 0.5 µmol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

Ordering Information

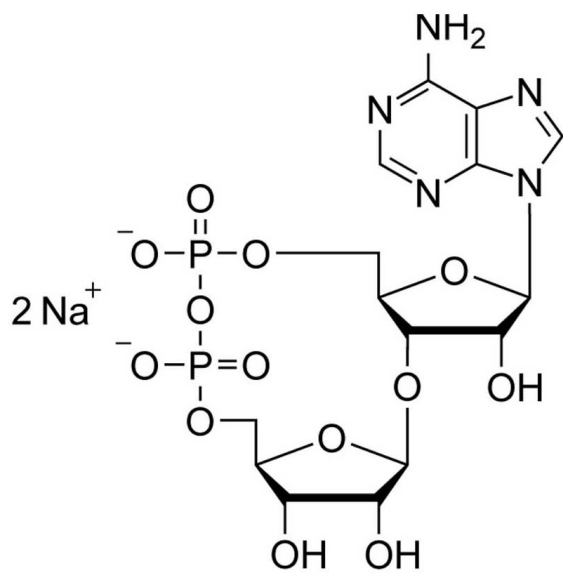
[Order Online »](#)

BLG-C404-005	0.5µmol
BLG-C404-025	5x0.5µmol

Manuals, SDS & CofA

[View Online »](#)

3'-cADPR is a regioisomer of the canonical second messenger cADPR (Cat. No. BLG-C005) and of 2'-cADPR (Cat. No. BLG-C406). 3'-cADPR was identified as novel signaling molecule involved in the innate immunity in plants and bacteria.



Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (1"-3")-Cyclic adenosine diphosphate ribose . sodium salt, v2-cADPR, 1"-3'-gcADPR

Formula $C_{15}H_{21}N_5O_{13}P_2$ (free acid)

MW 541.3 (free acid)

Purity > 98% HPLC

Quantity 0.5 µmol are equivalent to ~ 270 µg



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