

7-cADPR / N7-cADPR

Analogue of canonical cADPR identified as new second messenger in type IV Thoeis defence system in bacteria

For other salt forms please inquire. 7-cADPR, a recently identified structural isomer of the canonical second messenger cADPR (Cat. No. C 005), is one of several new, biologically relevant endogenous cyclic nucleotides, including 2'-cADPR (Cat. No. C 406) and 3'-cADPR (Cat. No. C 404). 7-cADPR has been identified as a signaling molecule involved in the innate immunity defense system of bacteria. Following recognition of phage invasion, bacterial Toll/interleukin-1 receptor (TIR)-domain proteins produce the previously unknown 7-cADPR. 7-cADPR activates the bacterial caspase-like protease, which indiscriminately degrades cellular proteins, halting phage replication. Rousett et al. denote this TIR-caspase defense system as type IV Thoeis, which is abundant in bacteria and efficiently protects against phage propagation.

BLG-C411-005 (5 x 0.1 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

Ordering Information

[Order Online »](#)

BLG-C411-001	0.1 µmol
BLG-C411-005	5x0.1 µmol

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 (1''-7)-Cyclic adenosine diphosphate ribose . sodium salt

Appearance Solid

Couple Type Modified nucleotides

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

MW 541.3 (free acid)

Purity > 90 % HPLC

Quantity 0.1 µmol are equivalent to ~54 µg

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