

2',2'',2''',2''''-Tetra-O-Me-c-tetraAMP

Analogue of c-tetraAMP containing O-methyl modifications in all ribose 2'-positions

Cyclic oligoadenylates such as c-tetraAMP (Biolog Cat. No. C 335) were found to be novel bacterial second messengers involved in the Type III CRISPR-Cas-associated detection and degradation of invasive genetic elements in many prokaryotes. 2',2'',2''',2''''-Tetra-O-Me-c-tetraAMP is an analogue of c-tetraAMP containing O-methyl modifications in all ribose 2'-positions, which are expected to make the compound resistant to degradation by nucleases.

BLG-T066-025 (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-T066-005	0.5µmol
BLG-T066-025	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 2'-, 2'', 2''', 2''''-Tetra-O-methyl-cyclic tetraadenosine monophosphate . sodium salt

Couple Type Modified nucleotides

Formula $C_{44}H_{56}N_{20}O_{24}P_4$ (free acid)

MW 1372.94 (free acid)

Purity > 95% HPLC

Quantity 0.5µmol = ~0.7mg

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)