

c-pentaAMP

BLG-C394-025 (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

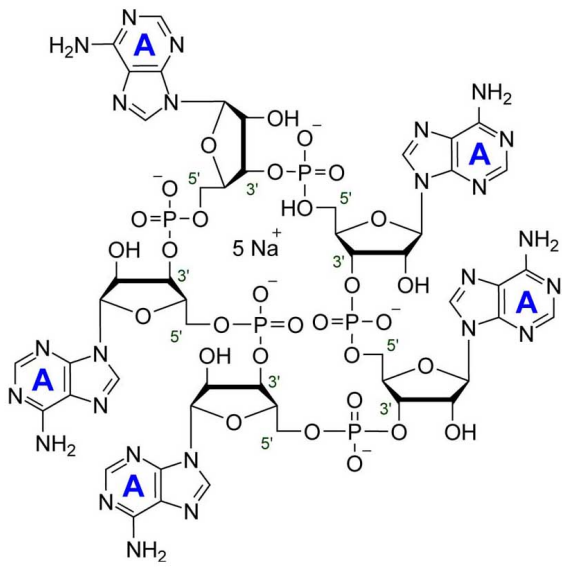
[Order Online »](#)

BLG-C394-005	0.5 μ mol
BLG-C394-025	5x0.5 μ mol

Manuals, SDS & CofA

[View Online »](#)

Cyclic oligoadenylates 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. The *Streptococcus thermophilus* type III-A CRISPR-Cas system produces c-pentaAMP as well as a range of differently sized cyclic oligoadenylates both in vitro and in vivo, although only the least abundant c-hexaAMP (Biolog Cat. No. C 332) activated the Csm6 ribonuclease of this organism.



Handling & Storage

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Cyclic pentaadenosine monophosphate . sodium salt, Cyclic pentaadenylate, c-A5, cOA5

Appearance Crystallized or lyophilized solid.

CAS 163499-42-1

Excitation Maximum 258 nm

Formula $C_{50}H_{60}N_{25}O_{30}P_5$ (free acid)

MW 1646.0 (free acid)

Purity > 95% HPLC

Quantity 0.1 µmol / ~ 0.8 mg

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