

2',2''-Di-AHC-c-diAMP

Functionalized c-diAMP for labelling or immobilization

Analogue of the bacterial second messenger c-diAMP (Cat. No. BLG-C088) which is suitable as a ligand in affinity chromatography or for coupling of various labelling structures including fluorophores.

BLG-D121-005 (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-D121-001	1µmol
BLG-D121-005	5x1µmol

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"-O-(Di-[6-aminohexylcarbamoyl])-cyclic diadenosine monophosphate . sodium salt

Appearance Crystallized or lyophilized sodium salt.

Couple Type Modified nucleotides

Formula $C_{34}H_{52}N_{14}O_{14}P_2$ (free acid)

MW 942.8 (free acid)

Purity > 95% HPLC

Quantity 0.1 μmol ≈ 94 μg

Solubility 2',2"-Di-AHC-c-diAMP is difficult to dissolve in water. However, the compound is soluble to approx. 1 mM in 20:80% (v/v) MeOH/H₂O. Please rinse tube walls carefully and preferably use ultrasonic or vortex to achieve total and uniform mixing. When opening the tube please make sure that no substance is lost within the cap.

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