

2'-AHC-cUMP

Functionalized cUMP for immobilization or conjugation with markers and dyes

Analogue of cUMP (Cat. No. U 001), in which a hexyl spacer with a terminal amino group has been attached to the ribose 2'-hydroxy group by a carbamate bond, suitable as a ligand for affinity chromatography of cUMP binding proteins or for coupling of various labelling structures including fluorophores. This ligand is also available immobilized on agarose gel (2'-AHC-cUMP-Agarose; Cat. No. A 164).

*BLG-A162-25 (5 x 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-A162-05	5 μ mol
BLG-A162-25	5x5 μ mol

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'-O-(6-Aminohexylcarbamoyl)uridine-3',5'-cyclic monophosphate . sodium salt

Couple Type Modified nucleotides

Formula $C_{16}H_{24}N_4O_9P \cdot Na$

MW 470.4

Purity > 98% HPLC

Quantity 5 µmol = ~2.4 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