

cCMP-AM

Highly membrane-permeant prodrug form of the second messenger cyclic CMP

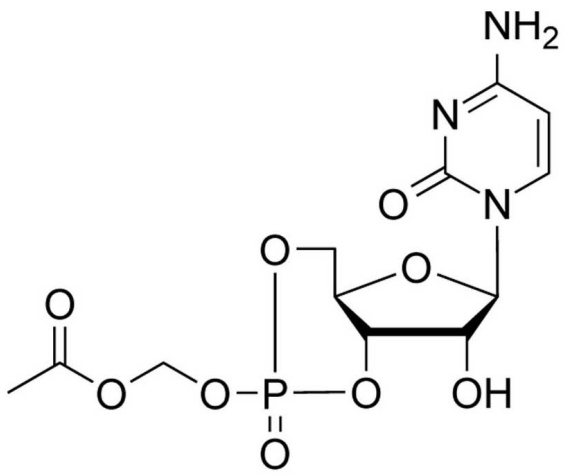
The acetoxymethyl group of cCMP-AM masks the charged polar phosphate and thus makes the molecule highly membrane-permeant. Inside the cell esterases release the polar cCMP which is trapped inside the cell and subsequently metabolized quickly, resulting in a pulse-type signal.

BLG-C067-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

Ordering Information [Order Online »](#)

BLG-C067-01	1 μ mol
BLG-C067-05	5x1 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 Cytidine-3',5'-cyclic monophosphate, acetoxymethyl ester

Couple Type Modified nucleotides

Formula $C_{12}H_{16}N_3O_9P$

MW 377.2

Purity > 95% HPLC

Quantity 1µmol≈0.4mg

Solubility Due to its rather high lipophilicity, the solubility of cCMP-AM in water or buffers is limited. It is recommended to use a small amount of organic solvent such as DMSO or DMF for dissolution and to dilute with water to the concentrations needed immediately before use.

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