

DB-cCMP

Highly membrane permeant, metabolically activatable prodrug form of cyclic CMP

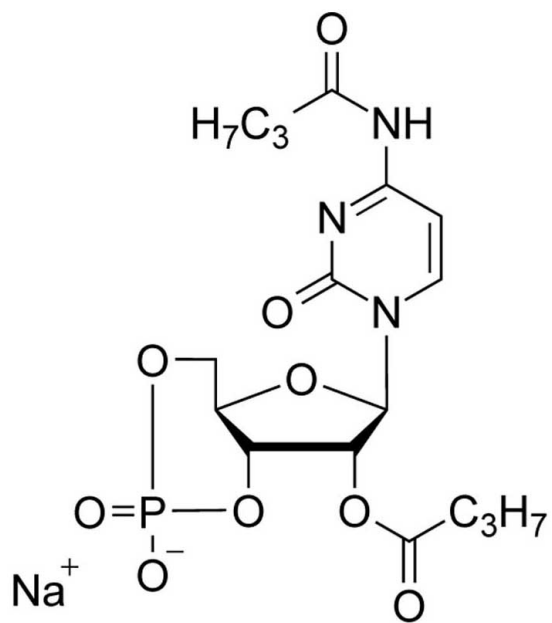
DB-cCMP is a membrane-permeant, metabolically activatable prodrug form of the potential second messenger cCMP (Prod. No. [BLG-C001](#)). Please note that the butyrate that is split off from DB-cCMP by esterases induces a lot of effects which are often reported to interfere with second messenger pathways. That means, that for each experiment with DB-cCMP, the influence of butyrate should be determined in control runs with sodium butyrate or tributyrin.

BLG-D075-50 (5 x 10 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

Ordering Information [Order Online »](#)

BLG-D075-10	10µmol
BLG-D075-50	5x10µmol

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 N4,2'-O-Dibutyrylcytidine-3',5'-cyclic monophosphate . sodium salt

CAS 64649-87-2

Couple Type Modified nucleotides

Formula $C_{17}H_{23}N_3O_9P \cdot Na$

MW 467.3

Purity > 97% HPLC

Quantity 10µmol≈4.7mg

Solubility Soluble in water (≥25mM).

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