

MANT-cUMP

Fluorescent, membrane-permeant cyclic UMP analogue for PDE studies

Fluorescent, membrane-permeant analogue of cUMP (Cat. No. U 001) with λ_{exc} 350 nm and λ_{em} 441 nm, which can be useful in phosphodiesterase studies. The compound shows an increase in fluorescence when cleaved by PDE.

*BLG-M068-25 (5 x 5 μmol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

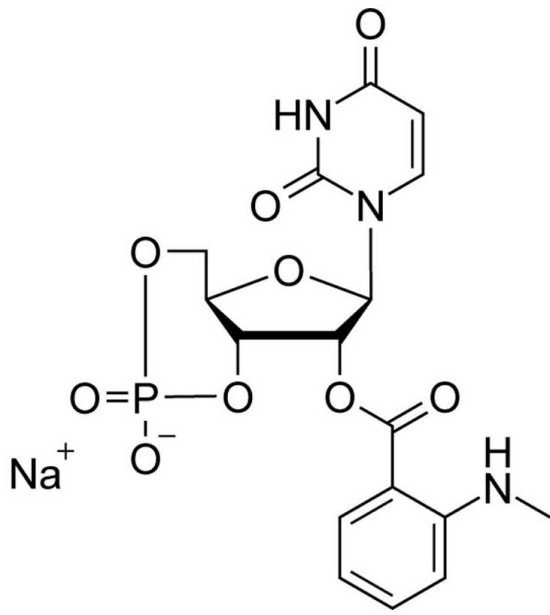
Ordering Information

[Order Online »](#)

BLG-M068-05	5 μmol
BLG-M068-25	5x5 μmol

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name 2'-O-(N'-Methylantraniloyl)uridine-3',5'-cyclic monophosphate . sodium salt

Appearance Solid.

CAS 1477518-35-6

Couple Type Modified nucleotides

Formula $C_{17}H_{17}N_3O_{19}P \cdot Na$

MW 461.3

Purity > 98% HPLC

Quantity 5µmol are equivalent to ~2.3 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