

# MANT-5'-CMP

## Fluorescent analogue of 5'-CMP

Fluorescent analogue of cytidine-5'-monophosphate ( $\lambda_{exc}$  355 nm,  $\lambda_{em}$  448 nm). The MANT fluorophore has a certain sensitivity for its environment and can change its spectral properties upon binding.

*BLG-M069-50 (5 x 10  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

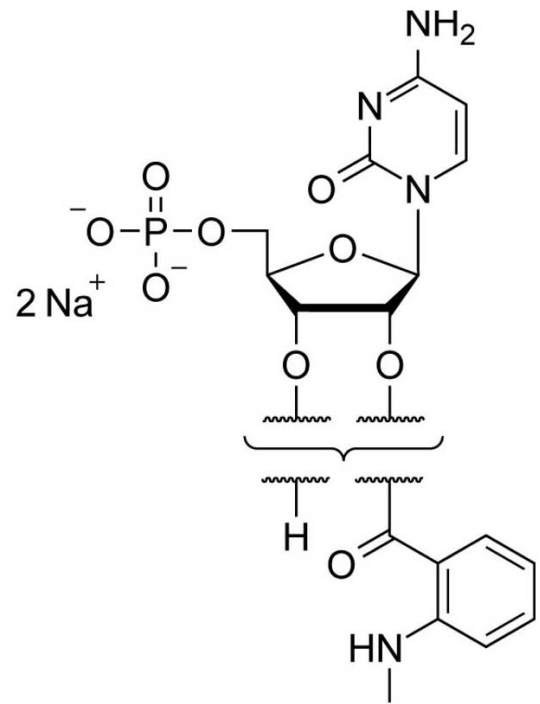
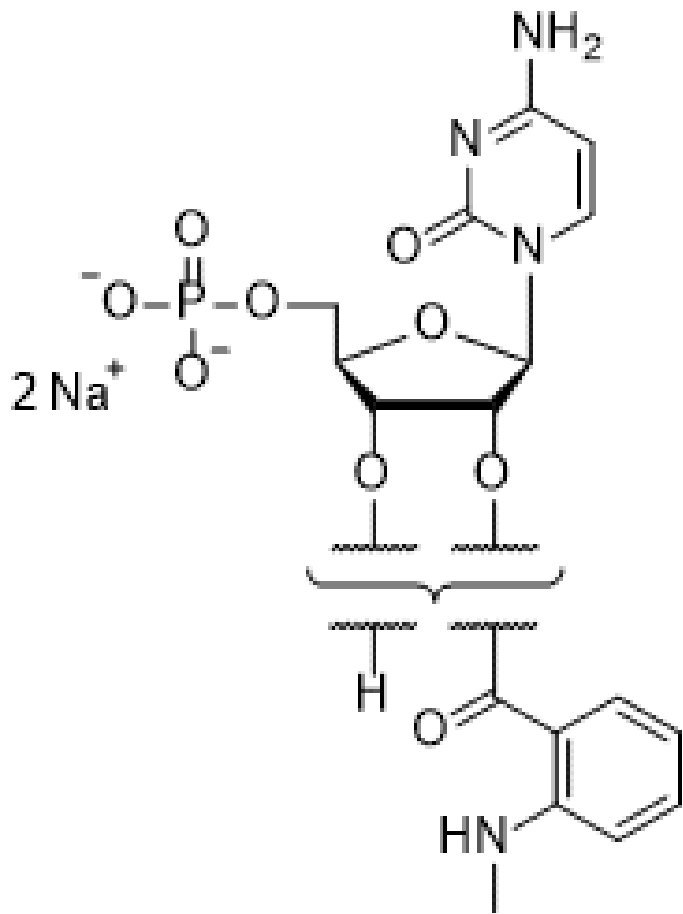
### Ordering Information

[Order Online »](#)

<b>BLG-M069-10</b>	10 $\mu$ mol
<b>BLG-M069-50</b>	5x10 $\mu$ 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'-/3'-O-(N'-Methylantraniloyl)cytidine-5'-O-monophosphate . sodium salt

**Appearance** Solid.

**CAS** 1477570-20-9

**Couple Type** Modified nucleotides

**Formula**  $C_{17}H_{21}N_4O_9P$  (free acid)

**MW** 456.4 (free acid)

**Purity** > 97% HPLC for mixture of isomers

**Quantity** 10µmol are equivalent to ~4.6 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](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)