

# BCMCM-caged cAMP

BCMCM-caged cAMP is an only weakly fluorescent, caged precursor of the second messenger cyclic AMP, which upon irradiation, releases both cAMP and a fluorescent coumarin analogue.

*BLG-B016-005 (5 x 0.1  $\mu$ mol pack size) is not sold in the U.S. or Canada.  
Please [contact us](#) for available options.*

## Ordering Information

[Order Online »](#)

<b>BLG-B016-001</b>	0.1 $\mu$ mol
<b>BLG-B016-005</b>	5x0.1 $\mu$ mol

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

Handling	Protect from light.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

Alternative Name	[6,7-Bis(carboxymethoxy)coumarin-4-yl]methyl adenosine-3',5'-cyclic monophosphate
Couple Type	Modified nucleotides
Formula	$C_{24}H_{22}N_5O_{14}P \cdot 2.5H_2O$
MW	635.5 . 45.0
Purity	>98% by HPLC
Quantity	0.1 μmol≈~68 μg
Solubility	Excellent solubility in water or buffer.
Technical Info / Product Notes	For the Original Manufacturer's data sheet please <a href="#">click here</a> .



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)