

# 8-[ $\phi$ -575]-2'-O-Me-cAMP

## Fluorescent Epac agonist

In 8-[ $\phi$ -575]-2'-O-Me-cAMP the relatively small Pharos 575 dye is connected to position 8 of the heterocyclic nucleobase of cyclic AMP (cAMP) via a spacer. In addition, the ribose 2'-hydroxy group has been methylated.

8-[ $\phi$ -575]-2'-O-Me-cAMP is a fluorescent analogue of cAMP with  $\lambda_{exc}$  577 nm and  $\lambda_{em}$  605 nm. The fluorophore can be excited e.g. by a Kr/Ar laser. Due to its high lipophilicity/membrane permeability and bright red fluorescence, 8-[ $\phi$ -575]-2'-O-Me-cAMP is especially suitable for studies with intact cells. Since a free 2'-ribose hydroxy group in cAMP is essential for stimulation of cAMP-dependent protein kinase (PKA), the methylated structure of 8-[ $\phi$ -575]-2'-O-Me-cAMP does not activate PKA, but is expected to be a selective stimulator of the exchange proteins directly activated by cAMP (Epac or cAMP-GEF).

*BLG-P021-005 (5 x 0.1  $\mu$ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

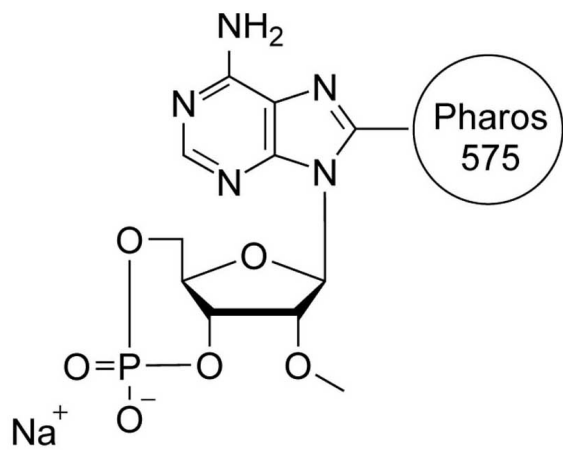
## Ordering Information

[Order Online »](#)

BLG-P021-001	0.1 $\mu$ mol
BLG-P021-005	5x0.1 $\mu$ mol

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 8-[Pharos-575]-2'-O-methyladenosine-3',5'-cyclic monophosphate . sodium salt

**Couple Type** Modified nucleotides

**MW** 731

**Purity** > 98% HPLC

**Quantity** 0.1µmol≈~74µg

**Solubility** Has limited solubility in water, but is better soluble in methanol.

**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)