

# 8-pCPT-2'-O-Me-5'-AMP

Metabolite of the Epac agonist 8-pCPT-2'-O-Me-cAMP

8-pCPT-2'-O-Me-5'-AMP is a potential metabolite of the specific Epac activator 8-pCPT-2'-O-Me-cAMP (Prod. No. [BLG-C041](#)), released by action of phosphodiesterases.

*BLG-C078-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

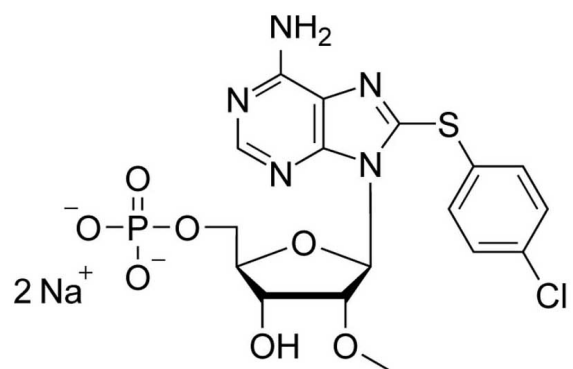
## Ordering Information

[Order Online »](#)

BLG-C078-05	5µmol
BLG-C078-25	5x5µmol

## 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** 8-(4-Chlorophenylthio)-2'-O-methyladenosine-5'-O-monophosphate . sodium salt

**CAS** 1187179-94-7

**Couple Type** Modified nucleotides

**Formula**  $C_{17}H_{19}ClN_5O_7PS$  (free acid)

**MW** 503.9

**Purity** > 97% HPLC

**Quantity** 5µmol≈2.5mg

**Solubility** Soluble in water or aqueous buffers.

**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)