

# 8-pCPT-2'-O-Me-Ado

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

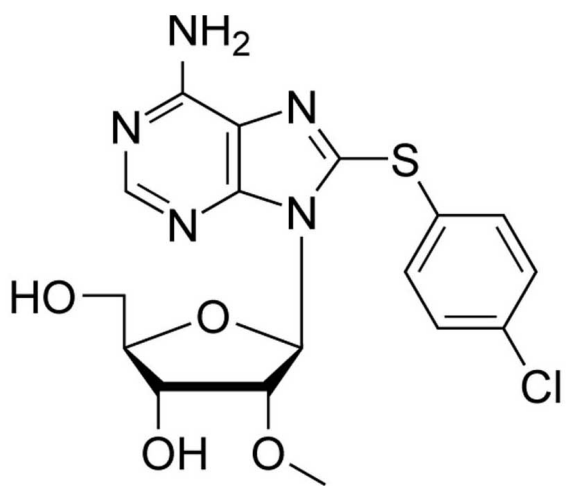
8-pCPT-2'-O-Me-Ado is a potential metabolite of the specific Epac activator 8-pCPT-2'-O-Me-cAMP (Prod. No. BLG-C041). The corresponding 5'-monophosphate 8-pCPT-2'-O-Me-5'-AMP (Prod. No. BLG-C078) as well as the nucleobase (Prod. No. BLG-C069) are available as well.

*BLG-C070-50 (5 x 10 µmol pack size) is not sold in the U.S. or Canada.  
Please [contact us](#) for available options.*

### Ordering Information [Order Online »](#)

BLG-C070-10	10µmol
BLG-C070-50	5x10µmol

### 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

CAS 1187179-96-9

Couple Type Modified nucleotides

Formula  $C_{17}H_{18}ClN_5O_4S$

MW 423.9

Purity > 98% HPLC

Quantity 10µmol≈4.2mg

Solubility Limited solubility in water and buffers. Soluble in DMSO or DMF

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)