

## 8-pCPT-2-O-Me-cAMP-AM

Kinase Activator

**Catalog # R46-905**

Lot # P3210-2

CAS # 1152197-23-3

### Product Description

Molecular Formula: C<sub>20</sub>H<sub>21</sub>ClN<sub>5</sub>O<sub>8</sub>PS

Appearance: White lyophilized solid

Melting Point: N/A

Molecular Weight: 557.9

Purity: >97% (HPLC)

Solubilization: May be dissolved in DMSO (50 mg/ml).

Poorly soluble in aqueous solution.

### Alias

8-(4-Chlorophenylthio)-2'-O-methyladenosine-3', 5'-cyclic monophosphate acetoxymethyl ester

### Specific Activity

Epac Pathway, RAP1 Pathway Activator

### Storage and Stability

Store desiccated as supplied at -20°C for up to 2 years.

Store solutions at -20°C for up to 1 month.

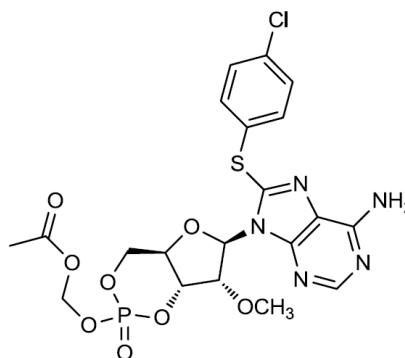
### Scientific Background

A potent, cell permeable Epac (exchange protein directly activated by cAMP) activator (1). Induces RAP1 activation and insulin secretion in pancreatic beta cell lines (2-5). Induces vascular relaxation in rate mesenteric artery (6). The acetoxymethyl ester confers increased cell-permeability and is cleaved by endogenous esterases to yield the active compound, 8-pCPT-2'-O-Me-cAMP. Addition to cell cultures should be done in serum-free media as esterases in the serum will cleave the acetoxymethyl ester and reduce cell permeability.

### References

1. MJ Vliem *et al.* *Chembiochem. Sci.* 2008 9:2052.
2. OG Chepurmy *et al.* *J. Biol. Chem.* 2009 284:10726.
3. GG Kelley *et al.* *Islets* 2009 1:260.
4. OG Chepurmy *et al.* *Am. J. Physiol. Endocrinol. Metab.* 2010 298:E622.
5. I Dzhura *et al.* *Islets* 2011 3:121.
6. OL Roberts *et al.* *J. Physiol.* 2013 591:5107.

### Molecular Structure



## 8-pCPT-2-O-Me-cAMP-AM

Kinase Activator

Catalog #	R46-905
Lot #	P3210-2
CAS #	1152197-23-3
Purity	>97% (HPLC)
Format	Lyophilized powder
Solubilization	May be dissolved in DMSO (50 mg/ml). Poorly soluble in aqueous solutions.
Storage & Shipping	Store desiccated as supplied at -20°C for up to 2 years. Store solutions at -20°C for up to 1 month.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**