

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

8-pCPT-2'-O-Me-Guo is a potential metabolite of the PKG- and Epac-inactive cGMP analogue 8-pCPT-2'-O-Me-cGMP (Prod. No. [BLG-C048](#)).

*BLG-C102-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.*

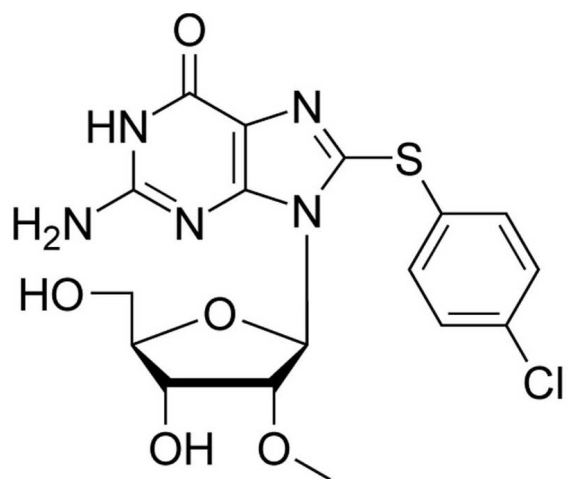
## Ordering Information

[Order Online »](#)

BLG-C102-05	5µmol
BLG-C102-25	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-methylguanosine

CAS 956935-50-5

Couple Type Modified nucleotides

Formula  $C_{17}H_{18}ClN_5O_5S$

MW 439.9

Purity > 97% HPLC

Quantity 5µmol≈2.2mg

**Solubility** Due to its increased lipophilicity, the solubility of 8-pCPT-2'-O-Me-Guo in water or buffer is limited. It is recommended to use a small amount of anhydrous organic solvent such as DMSO or DMF for dissolution, and to dilute with water or buffer down to the concentrations required.

**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)