

8-pCPT-Ade

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

8-pCPT-Ade is a potential metabolite of the specific Epac activator 8-pCPT-2'-O-Me-cAMP (Prod. No. [BLG-C041](#)). The corresponding nucleoside 8-pCPT-2'-O-Me-Ado (Prod. No. [BLG-C070](#)) and the 5'-O-monophosphate 8-pCPT-2'-O-Me-5'-AMP (Prod. No. [BLG-C078](#)) are available as well.

BLG-C069-50 (5 x 10 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

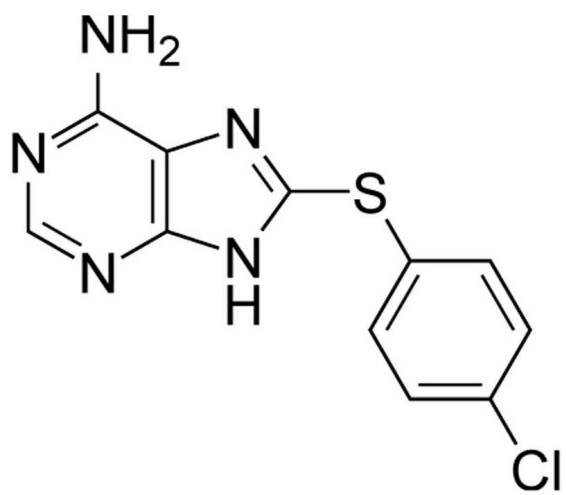
Ordering Information

[Order Online »](#)

BLG-C069-10	10µmol
BLG-C069-50	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)adenine

CAS 696574-61-5

Couple Type Modified nucleotides

Formula $C_{11}H_8ClN_5S$

MW 277.7

Purity > 95% HPLC

Quantity 10µmol≈2.8mg

Solubility 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

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-be@enzolifesciences.com

France
Phone: +33 472 440 655
info-fr@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-de@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com