

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

Monophosphate sodium salt

Specific, potent and membrane permeant activator of the Epac cAMP receptor. Activates Epac1 and Epac2, the guanine nucleotide-exchange factors (GEFs) for the small GTPases Rap1 and Rap2. Does not activate protein kinase A (PKA), neither *in vitro* nor *in vivo* and is therefore a valuable tool to specifically discriminate between PKA- and Epac-mediated effects. Since a free 2'-ribose hydroxyl group in cyclic AMP is essential for stimulation of PKA, the methylated structure of 8-CPT-2'-O-Me-cAMP is an extremely poor PKA activator. Its high lipophilicity allows for good membrane permeability in most biosystems and its increased resistance towards phosphodiesterases prevents rapid hydrolysis.

Citations: 3

[View Online »](#)

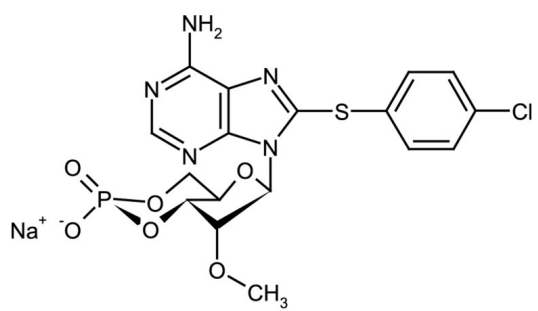
Ordering Information

[Order Online »](#)

ALX-480-093-M001	1mg
ALX-480-093-M005	5mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended.
Handling	Protect from light. Hygroscopic.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	8-CPT-2'-O-Me-cAMP . Na
Appearance	White to off-white crystalline powder.
CAS	510774-50-2
Couple Target	cAMP receptor, GEF
Couple Type	Activator
Formula	$C_{17}H_{16}ClN_5O_6PS \cdot Na$
Identity	Determined by ESI-MS and UV.
MW	484.8 . 23.0
Purity	≥98% (HPLC)
Purity Detail	Product is not sterile.
Solubility	Soluble in water, aqueous buffers, DMSO, dimethyl formamide or methanol.

Last modified: May 29, 2024



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)