

8-pHPT-2'-O-Me-cAMP

Potent Epac agonist

Analogue of the natural signal molecule cyclic AMP in which the hydrogen in position 8 of the heterocyclic nucleobase is replaced by the lipophilic 4-hydroxyphenylthio moiety. In addition, the ribose 2'-hydroxy group has been methylated.

Potent stimulator of exchange factors directly activated by cAMP (Epac or cAMP-GEF), a newly discovered receptor for cyclic AMP. Since a free 2'-ribose hydroxyl group in cyclic AMP is essential for stimulation of cAMP-dependent protein kinase (PKA), the methylated structure of 8-pHPT-2'-O-Me-cAMP is an extremely poor PKA activator and allows for specific discrimination between both signalling pathways.

Its high lipophilicity (> dibutyryl-cAMP), allows for good membrane permeability in most biosystems, and its increased resistance towards phosphodiesterases prevents from rapid hydrolysis.

BLG-H010-25 (5 x 5 µmol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

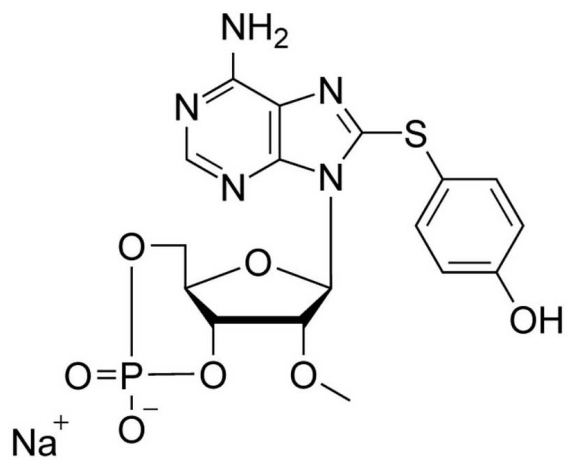
Ordering Information

[Order Online »](#)

BLG-H010-05	5µmol
BLG-H010-25	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-Hydroxyphenylthio)-2'-O-methyladenosine-3',5'-cyclicmonophosphate . sodiumsalt

Appearance Lyophilized or crystallized sodium salt.

CAS 612513-15-2

Couple Type Modified nucleotides

Formula $C_{17}H_{17}N_5O_7PS \cdot Na$

MW 489.4

Purity > 98% HPLC

Quantity 5µmol≈2.4mg

Solubility 8-pHPT-2'-O-Me-cAMP has limited solubility in water or buffers, however at least a 100 mM aqueous can be achieved. Please rinse tube walls carefully and preferably use ultrasonic or vortex to achieve total and uniform mixing.

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)