

Sp-6-Phe-cAMPS

PDE-resistant form of N6-Phenyl-cAMP

Sp-6-Phe-cAMPS is a hydrolysis-resistant form of 6-Phe-cAMP (Prod. No. [BLG-P006](#)), a potent site-selective and membrane-permeant activator of cAMP-dependent protein kinase. Sp-6-Phe-cAMPS does not activate the exchange factors directly activated by cAMP (Epac or cAMP-GEF) and thus can be used as an Epac-negative control.

BLG-P018-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

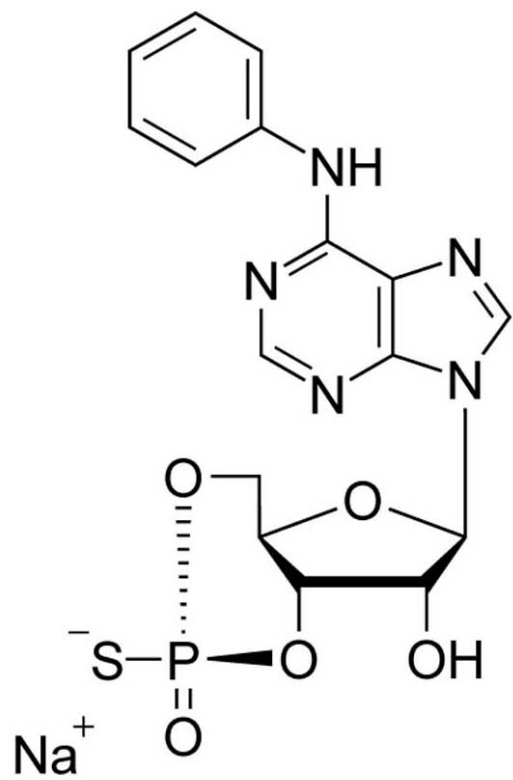
Ordering Information

[Order Online »](#)

BLG-P018-05	5 μ mol
BLG-P018-25	5x5 μ mol

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 N6-Phenyladenosine-3',5'-cyclic monophosphorothioate, Sp-isomer, sodium salt

CAS 169335-92-6

Couple Type Modified nucleotides

Formula $C_{16}H_{15}N_5O_5PS \cdot Na$

MW 443.4

Purity > 98% HPLC

Quantity 5µmol≈2.2mg

Solubility Soluble in water.

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