

Sp-8-Br-cAMPS

Membrane-permeant and PDE-resistant PKA agonist

Metabolically resistant activator of protein kinase A (cyclic AMP agonist). No metabolic side effects as observed with dibutyryl-cAMP or 8-Br-cAMP. Significantly more lipophilic and membrane permeant compared to Sp-cAMPS.

BLG-B002-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

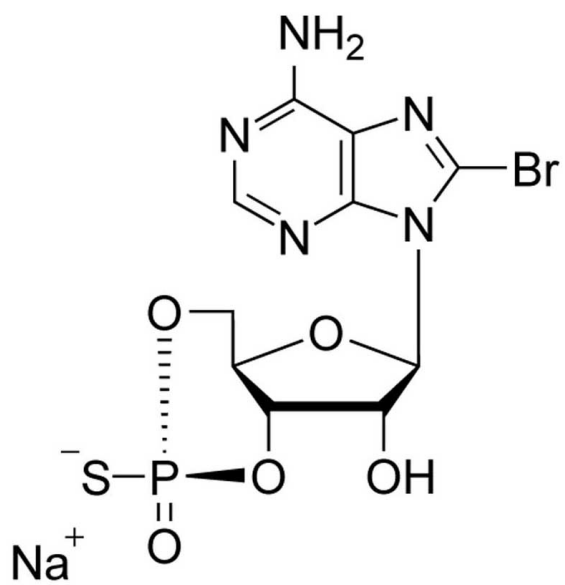
Ordering Information

[Order Online »](#)

BLG-B002-25	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 8-Bromoadenosine-3',5'-cyclic monophosphorothioate . sodium salt, Sp-isomer

CAS 127634-20-2

Couple Type Modified nucleotides

Formula $C_{10}H_{10}BrN_5O_5PS \cdot Na$

MW 446.2

Purity > 99% HPLC

Quantity 5µmol≈2.2mg

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