

8-APT-cGMP

Isozyme-selective activator of PKG I α

Isozyme-selective activator of cGMP-dependent protein kinase I alpha which prefers type I α over I β by a factor of 200.

BLG-A032-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

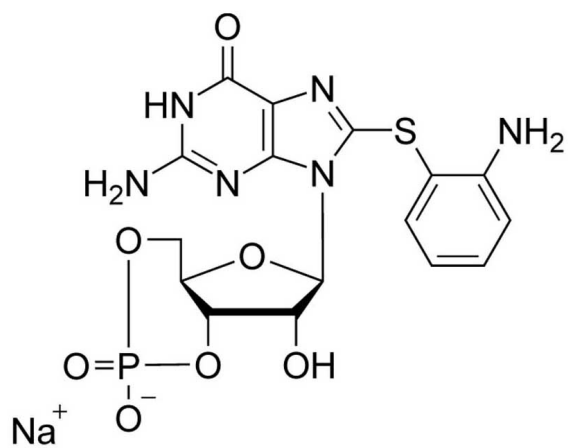
Ordering Information

[Order Online »](#)

BLG-A032-05	5 μ mol
BLG-A032-25	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-(2-Aminophenylthio)guanosine-3',5'-cyclic monophosphate . sodium salt

CAS 144509-87-5

Couple Type Modified nucleotides

Formula $C_{16}H_{16}N_6O_7PS \cdot Na$

MW 490.4

Purity > 98% HPLC

Quantity 5µmol≈2.4mg

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