

8-NBD-cAMP

Fluorescent cyclic AMP analogue

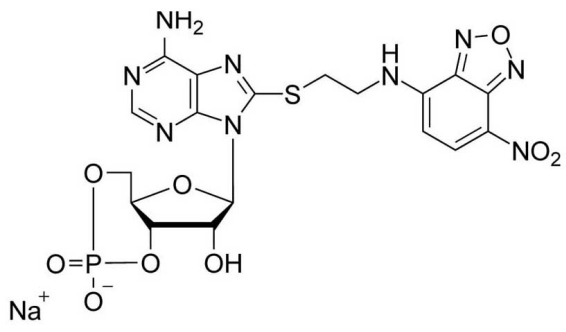
Potent membrane permeant fluorescent activator of cAMP-dependent protein kinase. While 8-NBD-cAMP is almost non-fluorescent in aqueous solutions it fluoresces strongly in hydrophobic environments such as in hydrophobic protein binding site.

BLG-N002-05 (5 x 1 μ mol pack size) is not sold in the U.S. or Canada. Please [contact us](#) for available options.

Ordering Information [Order Online »](#)

BLG-N002-01	1 μ mol
BLG-N002-05	5x1 μ mol

Manuals, SDS & CofA [View Online »](#)



Handling & Storage

Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	8-(2-[7-Nitro-4-benzofurazanyl]aminoethylthio)adenosine-3',5'-cyclic monophosphate . sodium salt
CAS	221905-51-7
Couple Type	Modified nucleotides
Formula	$C_{18}H_{17}N_9O_9PS \cdot Na$
MW	589.4
Purity	> 95% HPLC
Quantity	1µmol≈0.6mg
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)