

2-CI-8-HA-cAMP

Cyclic AMP analogue with high selectivity for site B I of PKA and good membrane permeability. Combined with AI-selective analogues synergistic activation of PKA type I is achieved.

*BLG-C071-25 (5 x 5 μ mol pack size) is not sold in the U.S. or Canada.
Please [contact us](#) for available options.*

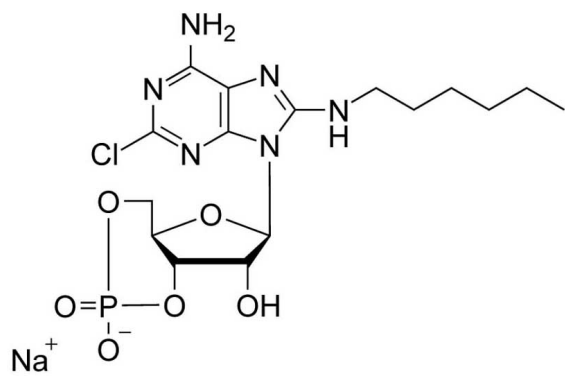
Ordering Information

[Order Online »](#)

BLG-C071-05	5 μ mol
BLG-C071-25	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 2-Chloro-8-n-hexylaminoadenosine-3',5'-cyclic monophosphate . sodium salt

Couple Type Modified nucleotides

Formula $C_{16}H_{23}ClN_6O_6P \cdot Na$

MW 484.8

Purity > 98% HPLC



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