

Bucladesine . sodium

Cell-permeable cAMP analog and PDE inhibitor

Bucladesine . sodium, is a cell-permeable cyclic AMP (cAMP) analog that mimics the action of endogenous cAMP. It activates protein kinase A (PKA) and inhibits phosphodiesterases (PDEs), leading to elevated intracellular cAMP levels. This compound is widely used in research to study cAMP-mediated signaling pathways, inflammation, and wound healing.

Key features and applications include:

- **Selective Inhibition:** Targets ALK5/4/7 with high potency, effectively blocking Smad2/3 phosphorylation.
- **cAMP Mimetic:** Activates cAMP-dependent protein kinase (PKA) pathways.
- **Phosphodiesterase Inhibitor:** Enhances intracellular cAMP levels by resisting enzymatic breakdown.
- **Cell-Permeable:** Easily crosses cell membranes to exert intracellular effects.

Research Applications:

- **Neuroscience:** Investigates memory, learning, and synaptic plasticity.
- **Cancer Biology:** Explores cAMP's role in cell proliferation and apoptosis.
- **Dermatology:** Used in models of acute skin inflammation.
- **Cell Signaling:** Studies second messenger systems and PKA activation.

Relevant disease states include:

- **Inflammatory Skin Conditions:** Demonstrates efficacy in reducing acute skin inflammation.
- **Cognitive Disorders:** Enhances spatial memory in preclinical models, supporting its use in neurodegenerative disease research.

- **Wound Healing Impairments:** Used to study and promote healing in damaged or diseased skin.
- **Pain and Sensory Disorders:** Investigated for its role in modulating pain perception via cAMP signaling.

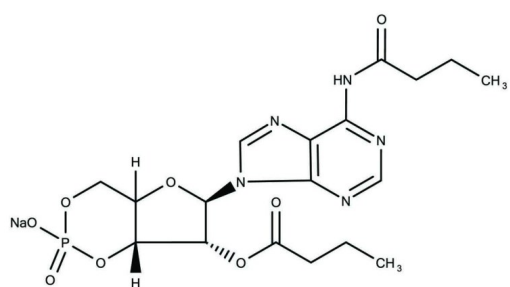
Ordering Information

Order Online »

ENZ-CHM363-0100	100mg
-----------------	-------

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Short Term Storage -20°C

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Dibutyl cAMP . sodium salt, DBcAMP . sodium salt

Appearance White solid.

CAS 16980-89-5

Couple Target PDE, PKA

Couple Type Activator, Inhibitor

Formula $C_{18}H_{23}N_5O_8P \cdot Na$

Identity Determined by NMR.

MW 491.37

Purity ≥98% (HPLC)

Solubility Soluble in DMSO (up to 50 mg/mL) or in water (up to 50 mg/mL).

Last modified: July 28, 2025



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-
fr@enzolifesciences.com](mailto:info-fr@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)