

EHNA

PDE inhibitor

Phosphodiesterase PDE II inhibitor, $IC_{50}=0.8\text{ }\mu\text{M}$. Also inhibits adenosine deaminase, $K_i=4\text{ nM}$.

Citations: 4

[View Online »](#)

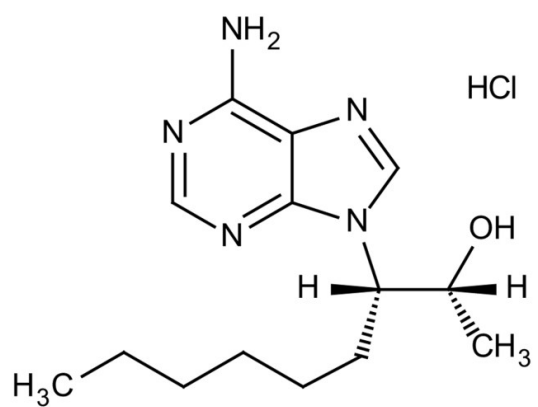
Ordering Information

[Order Online »](#)

BML-PD129-0010	10mg
BML-PD129-0050	50mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage Ambient

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name Erythro-9-(2-hydroxy-3-nonyl)adenine . HCl

Appearance White solid.

CAS 51350-19-7

Couple Target Adenosine Deaminase, PDE

Couple Type Inhibitor

Formula $C_{14}H_{23}N_5O \cdot HCl$

MW 313.8

Purity $\geq 98\%$ (HPLC)

Solubility Soluble in water (25mg/ml).

Last modified: May 29, 2024



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)