

(6R)-5,6,7,8-Tetrahydro-L-biopterin . dihydrochloride

NOS cofactor

Cofactor of nitric oxide synthases (NOS).

Citations: 13

[View Online »](#)

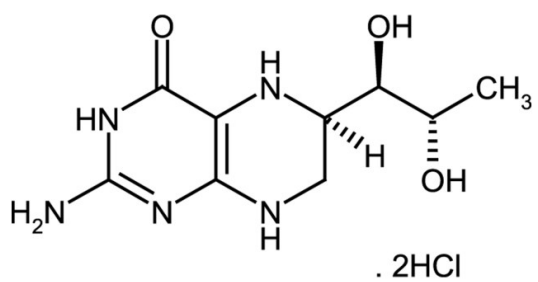
Ordering Information

[Order Online »](#)

ALX-440-001-M005	5mg
ALX-440-001-M025	25mg
ALX-440-001-M100	100mg

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for several years when stored at -20°C. Sensitive to oxidation, especially in neutral and alkaline solutions. Solutions must be stored at -20°C or colder immediately after preparation. Use oxygen free water with the lowest possible pH. Ascorbic acid, DTT, or other antioxidants may be added in order to increase stability.
Handling	Hygroscopic.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	(6R)-BH4 . 2HCl, [6R-[6R*(1R*,2S*)]]-2-Amino-6-(1,2-dihydroxypropyl)-5,6,7,8-tetrahydro-4(1H)-pteridinone . 2HCl, Sapropterin . dihydrochloride
Appearance	White to off-white powder.
CAS	69056-38-8
Formula	$C_9H_{15}N_5O_3 \cdot 2HCl$
MI	14: 8368
MW	314.2
Purity	≥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-
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)