

L-Biopterin

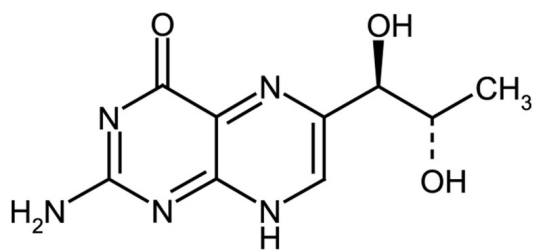
Ordering Information

[Order Online »](#)

ALX-440-008-M100	100mg
------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Stable for 2 years after receipt when stored at -20°C.
Handling	Protect from light.
Long Term Storage	-20°C
Shipping	Ambient Temperature

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	6-Biopterin, 2-Amino-6-[(1R,2S)-1,2-dihydroxypropyl]-4(1H)-pteridinone, 2-Amino-4-hydroxy-6-(1,2-dihydroxypropyl)pteridine, L-erythro-6-(1,2-Dihydroxypropyl)pterin
Appearance	White to pale yellow powder.
CAS	22150-76-1
Formula	$C_9H_{11}N_5O_3$
MI	14: 1229
MW	237.2
Purity	≥98% (HPLC)
RTECS	UO3506000
Solubility	Soluble in 1 N sodium hydroxide or 1N hydrogen chloride; very slightly soluble in water (0.7mg/ml) (ultrasonication). To obtain concentrated solutions, ammonium hydroxide should be added.

Last modified: May 29, 2024

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