

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

## NOS cofactor

Cofactor of nitric oxide synthases (NOS).

Citations: 13

[View Online »](#)

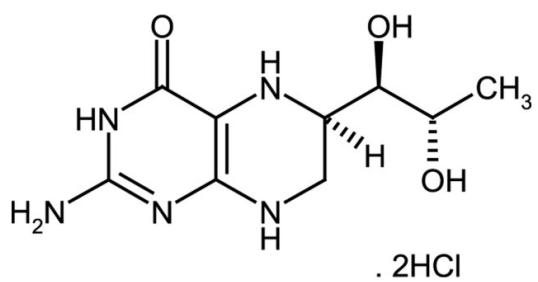
## Ordering Information

[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)