

Buthionine sulfoximine

γ -Glutamyl transpeptidase inhibitor

L-Buthionine-sulfoximine (BSO) induces oxidative stress in cells by irreversibly inhibiting γ -glutamylcysteine synthetase, an essential enzyme for the synthesis of glutathione.

Citations: 4

[View Online »](#)

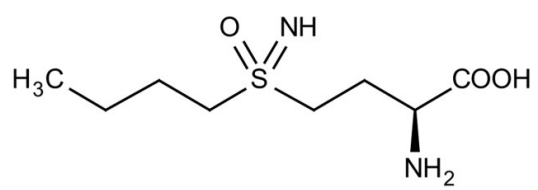
Ordering Information

[Order Online »](#)

BML-FR117-0500	500mg
-----------------------	-------

Manuals, SDS & CofA

[View Online »](#)



Handling & Storage

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

Long Term Storage -20°C

Shipping Ambient Temperature

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name	BSO
Appearance	White solid.
CAS	83730-53-4
Couple Target	gamma-Glutamyl transpeptidase
Couple Type	Inhibitor
Formula	$C_8H_{18}N_2O_3S$
MW	222.3
Purity	≥97% (TLC)
Solubility	Soluble in water.

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)