

# Lamotrigine

Na<sup>+</sup> channel blocker

Lamotrigine blocks voltage-gated sodium channels. Inhibits the release of glutamate. Neuroprotective. Anticonvulsant.

Citations: 16

[View Online »](#)

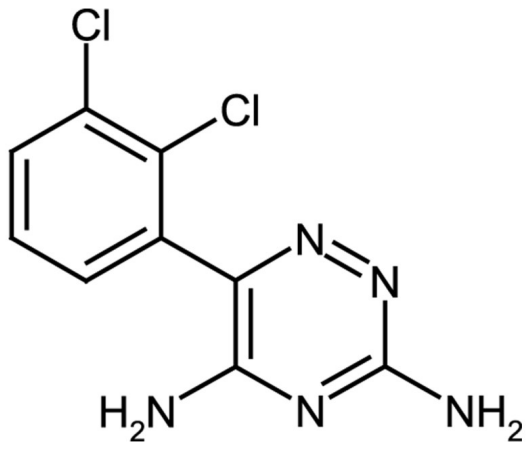
## Ordering Information

[Order Online »](#)

ALX-430-160-M100	100mg
ALX-430-160-M500	500mg

## Manuals, SDS & CofA

[View Online »](#)



## Handling & Storage

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

**Long Term Storage** +4°C

**Shipping** Ambient Temperature

**Regulatory Status** RUO - Research Use Only

## Product Details

**Alternative Name** 6-(2,3-Dichlorophenyl)-1,2,4-triazine-3,5-diamine, LTG, BW-430C

**Appearance** White to off-white powder.

**CAS** 84057-84-1

**Couple Target** Sodium channel

**Couple Type** Blocker

**Formula**  $C_9H_7Cl_2N_5$

**Identity** Determined by IR.

**MI** 14: 5353

**MW** 256.1

**Purity** ≥98% (Assay)

**RTECS** XY5850700

**Solubility** Soluble in DMSO (60°C, 20mg/ml) or dimethyl formamide; slightly soluble in anhydrous ethanol or dilute hydrochloric acid; almost insoluble in water.



ENZO LIFE SCIENCES,  
**Last updated May 29, 2024**  
Phone: 800.942.0430  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[usa@enzolifesciences.com](mailto:usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[eu@enzolifesciences.com](mailto:eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[be@enzolifesciences.com](mailto:be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[fr@enzolifesciences.com](mailto:fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[de@enzolifesciences.com](mailto:de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info@enzolifesciences.com](mailto:info@enzolifesciences.com)  
[uk@enzolifesciences.com](mailto:uk@enzolifesciences.com)