

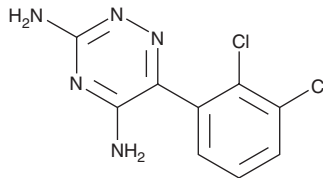
PRODUCT INFORMATION



Lamotrigine

Item No. 15428

CAS Registry No.: 84057-84-1
Formal Name: 6-(2,3-dichlorophenyl)-1,2,4-triazine-3,5-diamine
Synonyms: BW 430C, LTG
MF: C₉H₇Cl₂N₅
FW: 256.1
Purity: ≥98%
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Lamotrigine is supplied as a crystalline solid. A stock solution may be made by dissolving the lamotrigine in the solvent of choice, which should be purged with an inert gas. Lamotrigine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of lamotrigine in ethanol is approximately 2 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Lamotrigine is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, lamotrigine should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Lamotrigine has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

Lamotrigine is an anticonvulsant.¹ It inhibits voltage-gated sodium channels (Na_v) in HEK293 cells expressing recombinant human Na_v1.2, Na_v1.5, or Na_v1.8 (IC₅₀s = 10, 62, and 96 μM, respectively), as well as high voltage-activated calcium currents in isolated rat cortical neurons (IC₅₀ = 12.3 μM), an effect that can be reversed by the N-type calcium channel blocker ω-conotoxin GVIA (Item No. 24114) and P-type calcium channel blocker ω-agatoxin IVA (Item No. 21605).^{1,2} Lamotrigine protects against seizures induced by maximal electroshock (MES) in mice and rats (ED₅₀s = 10.1 and 7.4 μmol/kg, respectively).³ It also decreases mechanical allodynia in a rat model of neuropathic pain induced by spinal nerve ligation (ED₅₀ = 47 μmol/kg).¹ Formulations containing lamotrigine have been used in the treatment of epilepsy and bipolar disorder.

References

1. Drizin, I., Gregg, R.J., Scanio, M.J., *et al.* Discovery of potent furan piperazine sodium channel blockers for treatment of neuropathic pain. *Bioorg. Med. Chem.* **16(12)**, 6379-6386 (2008).
2. Stefani, A., Spadoni, F., Siniscalchi, A., *et al.* Lamotrigine inhibits Ca²⁺ currents in cortical neurons: Functional implications. *Eur. J. Pharmacol.* **307(1)**, 113-116 (1996).
3. Miller, A.A., Wheatley, P., Sawyer, D.A., *et al.* Pharmacological studies on lamotrigine, a novel potential antiepileptic drug: I. Anticonvulsant profile in mice and rats. *Epilepsia* **27(5)**, 483-489 (1986).

WARNING

THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA

This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the complete Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY

Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 09/29/2022

CAYMAN CHEMICAL

1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA

PHONE: [800] 364-9897
[734] 971-3335

FAX: [734] 971-3640

CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM