

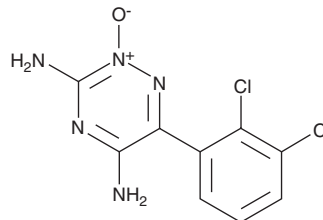
# PRODUCT INFORMATION



## Lamotrigine N2-oxide

Item No. 35524

**CAS Registry No.:** 136565-76-9  
**Formal Name:** 6-(2,3-dichlorophenyl)-1,2,4-triazine-3,5-diamine 2-oxide  
**Synonym:** LTG N2-oxide  
**MF:** C<sub>9</sub>H<sub>7</sub>Cl<sub>2</sub>N<sub>5</sub>O  
**FW:** 272.1  
**Purity:** ≥95%  
**Supplied as:** A 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 N2-oxide is supplied as a solid. A stock solution may be made by dissolving the lamotrigine N2-oxide in the solvent of choice, which should be purged with an inert gas. Lamotrigine N2-oxide is slightly soluble in methanol and DMSO.

### Description

Lamotrigine N2-oxide is a metabolite of the anticonvulsant lamotrigine (Item No. 15428).<sup>1</sup>

### Reference

1. Doig, M.V. and Clare, R.A. Use of thermospray liquid chromatography-mass spectrometry to aid in the identification of urinary metabolites of a novel antiepileptic drug, Lamotrigine. *J. Chromatogr.* **554**(1-2), 181-189 (1991).

#### 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, 11/18/2024

#### 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