

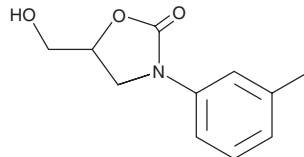
# PRODUCT INFORMATION



## Toloxatone

Item No. 33413

**CAS Registry No.:** 29218-27-7  
**Formal Name:** 5-(hydroxymethyl)-3-(3-methylphenyl)-2-oxazolidinone  
**Synonyms:** Humoryl, MD 69276  
**MF:** C<sub>11</sub>H<sub>13</sub>NO<sub>3</sub>  
**FW:** 207.2  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 239 nm  
**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

Toloxatone is supplied as a crystalline solid. A stock solution may be made by dissolving the tolloxatone in the solvent of choice, which should be purged with an inert gas. Tolloxatone is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of tolloxatone in these solvents is approximately 30 mg/ml.

### Description

Tolloxatone is an inhibitor of monoamine oxidase A (MAO-A; K<sub>i</sub> = 1.8 μM).<sup>1</sup> It is selective for MAO-A over MAO-B (K<sub>i</sub> = 44 μM). Tolloxatone (100 mg/kg) inhibits deamination of serotonin (5-HT) and norepinephrine in rat brain.<sup>2</sup> It also decreases immobility time in the forced swim test in mice when administered at a dose of 256 mg/kg.<sup>3</sup>

### References

1. Cesura, A.M. and Pletscher, A. The new generation of monoamine oxidase inhibitors. *Prog. Drug Res.* **38**, 171-297 (1992).
2. Keane, P.E., Kan, J.P., Sontag, N., *et al.* Monoamine oxidase inhibition and brain amine metabolism after oral treatment with tolloxatone in the rat. *J. Pharm. Pharmacol.* **31(11)**, 752-754 (1979).
3. Van der Meersch-Mougeot, V., da Rocha, M., Jr., Monier, C., *et al.* Benzodiazepines reverse the anti-immobility effect of antidepressants in the forced swimming test in mice. *Neuropharmacology* **32(5)**, 439-446 (1993).

#### 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, 11/14/2022

#### 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](http://WWW.CAYMANCHEM.COM)