

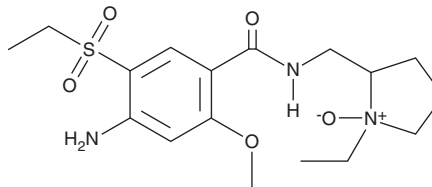
PRODUCT INFORMATION



Amisulpride N-oxide

Item No. 35447

CAS Registry No.: 71676-01-2
Formal Name: 4-amino-N-[(1-ethyl-1-oxido-2-pyrrolidiny)methyl]-5-(ethylsulfonyl)-2-methoxy-benzamide
MF: $C_{17}H_{27}N_3O_5S$
FW: 385.5
Purity: $\geq 95\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

Amisulpride N-oxide is supplied as a solid. A stock solution may be made by dissolving the amisulpride N-oxide in the solvent of choice, which should be purged with an inert gas. Amisulpride N-oxide is slightly soluble in chloroform and methanol.

Description

Amisulpride N-oxide is a degradation product of the dopamine and serotonin (5-HT) receptor antagonist amisulpride (Item No. 14619).¹ It is formed from amisulpride via oxidation by ozonation or photodegradation.² Amisulpride N-oxide has been found as a contaminant in surface water.³

References

1. Skibiński, R. Identification of photodegradation product of amisulpride by ultra-high-pressure liquid chromatography-DAD/ESI-quadrupole time-of-flight-mass spectrometry. *J. Pharm. Biomed. Anal.* **56**(5), 904-910 (2011).
2. Bollmann, A.F., Seitz, W., Prasse, C., *et al.* Occurrence and fate of amisulpride, sulpiride, and lamotrigine in municipal wastewater treatment plants with biological treatment and ozonation. *J. Hazard. Mater.* **320**, 204-215 (2016).
3. Diamanti, K.S., Alygizakis, N.A., Nika, M.-C., *et al.* Assessment of the chemical pollution status of the Dniester River Basin by wide-scope target and suspect screening using mass spectrometric techniques. *Anal. Bioanal. Chem.* **412**(20), 4893-4907 (2020).

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