

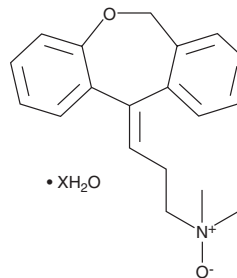
PRODUCT INFORMATION



Doxepin N-oxide (hydrate)

Item No. 9002302

Formal Name: 3-dibenz[b,e]oxepin-11(6H)-ylidene-N,N-dimethyl-1-propanamine, N-oxide, hydrate
MF: $C_{19}H_{21}NO_2 \cdot XH_2O$
FW: 295.4
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

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

Description

Doxepin N-oxide is a metabolite of the tricyclic antidepressant doxepin (Item No. 15888).¹ It is also a photodegradation product of doxepin.²

References

1. Maurer, H. and Pfleger, K. Screening procedure for detection of antidepressants and their metabolites in urine using a computerized gas chromatographic--Mass spectrometric technique. *J. Chromatogr.* **305**(2), 309-323 (1984).
2. Székely, P., Gyéresi, Á., Hancu, G., et al. Identification of photodegradation products of tricyclic antidepressants drugs clomipramine and doxepine. *Acta Medica Marisiensis* **57**(6), 745-747 (2011).

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