

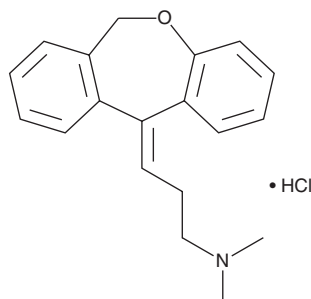
PRODUCT INFORMATION



Doxepin (hydrochloride)

Item No. 15888

CAS Registry No.: 1229-29-4
Formal Name: 3-(dibenz[b,e]oxepin-11(6H)-ylidene)-N,N-dimethyl-1-propanamine, monohydrochloride
MF: C₁₉H₂₁NO • HCl
FW: 315.8
Purity: ≥95%
UV/Vis.: λ_{max}: 296 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

Doxepin (hydrochloride) is supplied as a crystalline solid. A stock solution may be made by dissolving the doxepin (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Doxepin (hydrochloride) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of doxepin (hydrochloride) in these solvents is approximately 30, 25, and 20 mg/ml, respectively.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of doxepin (hydrochloride) can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of doxepin (hydrochloride) in PBS, pH 7.2, is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

Description

Doxepin is a tricyclic antidepressant that binds to the serotonin (5-HT) transporter (SERT) and norepinephrine transporter (NET; K_ds = 68 and 29.5 nM, respectively).¹ It is a histamine H₁ receptor antagonist (K_i = 1.23 nM).² Doxepin selectively binds to SERT and NET over the dopamine transporter (DAT; K_d = 12,100 nM) and inhibits histamine H₁ over H₂, H₃, and H₄ receptors (K_is = 170, 39,810, and 15,135 nM, respectively).^{1,2} It also binds to the 5-HT₂ receptor, as well as to muscarinic acetylcholine and α₁-adrenergic receptors (α₁-ARs; K_ds = 27, 23, and 23.5 nM, respectively).³ Doxepin (10 mg/kg, i.p.) decreases allodynia and hyperalgesia in a mouse model of chronic constriction injury-induced neuropathic pain.⁴ It increases the distance traveled in the center of the open field test and reduces immobility time in the forced swim test in a mouse model of depression induced by chronic stress when administered orally at a dose of 15 mg/kg.⁵ Formulations containing doxepin have been used in the treatment of depression and insomnia.

References

1. Tatsumi, M., Groshan, K., Blakely, R.D., et al. *Eur. J. Pharmacol.* **340(2-3)**, 249-258 (1997).
2. Appl, H., Holzammer, T., Dove, S., et al. *Maunyn-Schmiedeberg's Arch. Pharmacol.* **385**, 145-170 (2012).
3. Cusack, B., Nelson, A., and Richelson, E. *Psychopharmacology (Berl.)* **114(4)**, 559-565 (1994).
4. Mika, J., Jurga, A.M., Starnowska, J., et al. *Neuroscience* **294**, 38-50 (2015).
5. Kim, Y.R., Park, B.-K., Kim, Y.H., et al. *BioMed. Res. Int.* 8249563 (2018).

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, 11/28/2022

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