

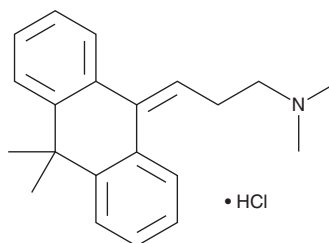
PRODUCT INFORMATION



Melitracen (hydrochloride)

Item No. 40117

CAS Registry No.: 10563-70-9
Formal Name: 3-(10,10-dimethyl-9(10H)-anthracenylidene)-N,N-dimethyl-1-propanamine, monohydrochloride
Synonyms: Melitracene, U-24973A
MF: C₂₁H₂₅N • HCl
FW: 327.9
Purity: ≥98%
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

Melitracen (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the melitracen (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Melitracen (hydrochloride) is soluble in the organic solvent DMSO. It is slightly soluble in acetonitrile.

Melitracen (hydrochloride) is slightly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

Description

Melitracen is a thymoleptic and tricyclic antidepressant.^{1,2} It inhibits serotonin (5-HT) uptake in rat brain synaptosomes (IC₅₀ = 670 nM).¹ Melitracen (20 and 40 mg/kg) increases the lethality of the α₂-adrenergic receptor (α₂-AR) antagonist yohimbine (Item No. 19869) in mice.²

References

1. Hyttel, J. Effect of a specific 5-HT uptake inhibitor, citalopram (Lu 10-171), on ³H-5-HT uptake in rat brain synaptosomes in vitro. *Psychopharmacol. (Berl)* **60(1)**, 13-18 (1978).
2. Lapin, I.P. Adrenergic nonspecific potentiation of yohimbine toxicity in mice by antidepressants and related drugs and antiyohimbine action of antiadrenergic and serotonergic drugs. *Psychopharmacol. (Berl)* **70(2)**, 179-185 (1980).

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, 02/12/2024

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