

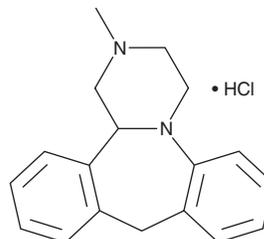
PRODUCT INFORMATION



Mianserin (hydrochloride)

Item No. 39378

CAS Registry No.: 21535-47-7
Formal Name: 1,2,3,4,10,14b-hexahydro-2-methyl-dibenzo[c,f]pyrazino[1,2-a]azepine, monohydrochloride
Synonyms: (±)-Mianserin, NSC 292267
MF: C₁₈H₂₀N₂ • HCl
FW: 300.8
Purity: ≥98%
Supplied as: A solid
Storage: -20°C
Stability: ≥3 years



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

Description

Mianserin is an analytical reference standard categorized as an antidepressant.¹ Formulations containing mianserin have been used in the treatment of depression and anxiety. This product is intended for research and forensic applications.

Reference

1. Pinder, R.M. Adrenoreceptor interactions of the enantiomers and metabolites of mianserin: Are they responsible for the antidepressant effect? *Acta Psychiatr. Scand. Suppl.* **320**, 1-9 (1985).

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, 12/13/2023

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