

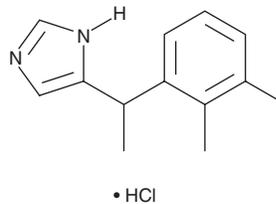
PRODUCT INFORMATION



Medetomidine (hydrochloride)

Item No. 38582

CAS Registry No.: 86347-15-1
Formal Name: 5-[1-(2,3-dimethylphenyl)ethyl]-1H-imidazole, monohydrochloride
Synonym: MPV 785
MF: C₁₃H₁₆N₂ • HCl
FW: 236.7
Purity: ≥98%
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.

Description

Medetomidine (hydrochloride) (Item No. 38582) is an analytical reference standard categorized as a sedative.¹ Medetomidine induces sedation in dogs. It has analgesic activity alone and an additive analgesic effect in combination with fentanyl (Item Nos. ISO60197 | 14719 | 23580) or morphine (Item Nos. ISO60147 | 15464) in rats.² Medetomidine has been found in fentanyl samples seized by law enforcement.³ Formulations containing medetomidine have been used as veterinary sedatives. This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 16454).

References

1. Vainio, O. and Vähä-Vahe, T. Reversal of medetomidine sedation by atipamezole in dogs. *J. Vet. Pharmacol. Ther.* **13(1)**, 15-22 (1990).
2. Ossipov, M.H., Harris, S., Lloyd, P., et al. Antinociceptive interaction between opioids and medetomidine: Systemic additivity and spinal synergy. *Anesthesiology* **73(6)**, 1227-1235 (1990).
3. Sisco, E. and Appley, M. Identification of the veterinary sedative medetomidine in combination with opioids and xylazine in Maryland. *J. Forensic Sci.* (2023).

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, 04/11/2025

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