

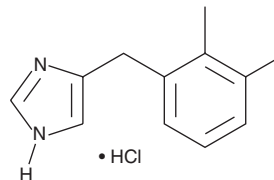
PRODUCT INFORMATION



Detomidine (hydrochloride)

Item No. 38583

CAS Registry No.: 90038-01-0
Formal Name: 5-[(2,3-dimethylphenyl)methyl]-1H-imidazole, monohydrochloride
Synonym: MPV 253All
MF: C₁₂H₁₄N₂ • HCl
FW: 222.7
Purity: ≥98%
UV/Vis.: λ_{max}: 214 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.

Description

Detomidine (hydrochloride) (Item No. 38583) is an analytical reference standard categorized as a sedative.¹ Detomidine has sedative activity in horses.² Formulations containing detomidine 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. 17853).

Reference

1. Grimsrud, K.N., Mama, K.R., Thomasy, S.M., *et al.* Pharmacokinetics of detomidine and its metabolites following intravenous and intramuscular administration in horses. *Equine Vet. J.* **41**(4), 361-365 (2009).
2. Freeman, S.L. and England, G.C.W. Investigation of romifidine and detomidine for the clinical sedation of horses. *Vet. Rec.* **147**(18), 507-511 (2000).

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, 03/10/2025

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