

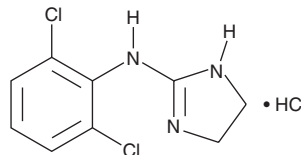
PRODUCT INFORMATION



Clonidine (hydrochloride)

Item No. 38558

CAS Registry No.: 4205-91-8
Formal Name: N-(2,6-dichlorophenyl)-4,5-dihydro-1H-imidazol-2-amine, monohydrochloride
Synonym: ST 155
MF: C₉H₉Cl₂N₃ • HCl
FW: 266.6
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

Clonidine (hydrochloride) (Item No. 38558) is an analytical reference standard categorized as an α_2 -adrenergic receptor agonist.¹ Formulations containing clonidine have been used in the treatment of opioid withdrawal.² This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 15949).

References

1. Jasper, J.R., Lesnick, J.D., Chang, L.K., *et al.* Ligand efficacy and potency at recombinant α_2 adrenergic receptors: Agonist-mediated [³⁵S]GTPγS binding. *Biochem. Pharmacol.* **55(7)**, 1035-1043 (1998).
2. Kuzmaul, A.K., Palmer, E.C., and Frederick, E.K. Lofexidine versus clonidine for mitigation of opioid withdrawal symptoms: A systematic review. *J. Am. Pharm. Assoc.* (2003) **60(1)**, 145-152 (2020).

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, 02/26/2025

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