

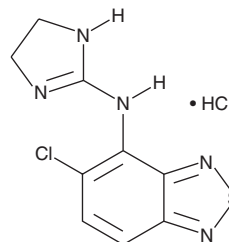
PRODUCT INFORMATION



Tizanidine (hydrochloride)

Item No. 38559

CAS Registry No.: 64461-82-1
Formal Name: 5-chloro-N-(4,5-dihydro-1H-imidazol-2-yl)-2,1,3-benzothiadiazol-4-amine, monohydrochloride
Synonym: DS 103-282
MF: C₉H₈ClN₅S • HCl
FW: 290.2
Purity: ≥98%
UV/Vis.: λ_{max}: 227, 320 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: ≥2 years



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

Description

Tizanidine (hydrochloride) (Item No. 38559) is an analytical reference standard categorized as a muscle relaxant.^{1,2} Tizanidine inhibits α - and γ -rigidity in rats, reflex muscle tone in rabbits, and the linguomandibular reflex in cats.² It has been associated with fatal overdoses.³ Formulations containing tizanidine have been used in the treatment of spasticity. This product is intended for analytical forensic applications. This product is also available as a general research tool (Item No. 16477).

References

1. Sakitani, K. The effects of centrally acting muscle relaxants on the intrathecal noradrenaline-induced facilitation of the flexor reflex mediated by group II afferent fibers in rats. *Jpn. J. Pharmacol.* **63(3)**, 369-376 (1993).
2. Sayers, A.C., Bürki, H.R., and Eichenberger, E. The pharmacology of 5-chloro-4-(2-imidazol-2-yl-amino)-2,1,3-benzothiadiazole (DS 103-282), a novel myotonolytic agent. *Arzneimittelforschung* **30(5)**, 793-803 (1980).
3. Sklerov, J.H., Cox, D.E., Moore, K.A., et al. Tizanidine distribution in a postmortem case. *J. Anal. Toxicol.* **30(5)**, 331-334 (2006).

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, 05/18/2023

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