

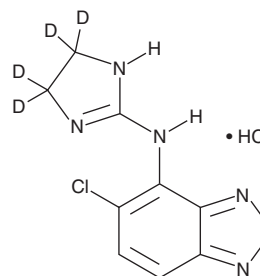
PRODUCT INFORMATION



Tizanidine-d₄ (hydrochloride)

Item No. 26521

CAS Registry No.: 1188263-51-5
Formal Name: 5-chloro-N-(4,5-dihydro-1H-imidazol-2-yl-4,4,5,5-d₄)-2,1,3-benzothiadiazol-4-amine, monohydrochloride
MF: C₉H₄ClD₄N₅S • HCl
FW: 294.2
Chemical Purity: ≥98% (Tizanidine)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₄); ≤1% d₀
Supplied as: A solid
Storage: -20°C
Stability: ≥4 years



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

Laboratory Procedures

Tizanidine-d₄ (hydrochloride) is intended for use as an internal standard for the quantification of tizanidine (Item No. 16477) by GC- or LC-MS. The accuracy of the sample weight in this vial is between 5% over and 2% under the amount shown on the vial. If better precision is required, the deuterated standard should be quantitated against a more precisely weighed unlabeled standard by constructing a standard curve of peak intensity ratios (deuterated versus unlabeled).

Tizanidine-d₄ (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the tizanidine-d₄ (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Tizanidine-d₄ (hydrochloride) is soluble in organic solvents such as methanol.

Description

Tizanidine is an α_2 -adrenergic receptor (α_2 -AR) agonist.¹ Administration of tizanidine reduces neuronal excitation induced by noxious stimuli and depresses spontaneous neuronal firing in anesthetized cats. Tizanidine-induced neuronal depression is reversed by the selective α_2 -AR antagonist RX781094 but not the α_1 -AR antagonists prazosin (Item No. 15023) and WB-4101. It acts as a muscle relaxant that inhibits α - and γ -rigidity in rats, reflex muscle tone in rabbits, and the linguomandibular reflex in cats without sedative, hemodynamic, or neurochemical effects.² Formulations containing tizanidine have been used in the treatment of spasticity.

References

1. Davies, J. and Quinlan, J.E. *Neuroscience* **16**(3), 673-682 (1985).
2. Sayers, A.C., Bürki, H.R., and Eichenberger, E. *Arzneimittelforschung* **30**(5), 793-803 (1980).

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, 05/17/2023

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