

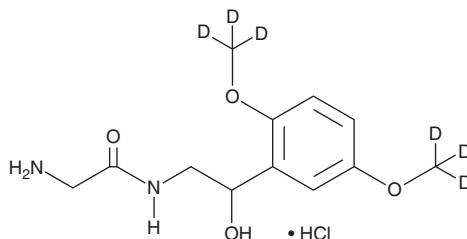
PRODUCT INFORMATION



Midodrine-d₆ (hydrochloride)

Item No. 33546

CAS Registry No.: 1188265-43-1
Formal Name: 2-amino-N-[2-[2,5-di(methoxy-d₃)phenyl]-2-hydroxyethyl]-acetamide, monohydrochloride
MF: C₁₂H₁₂D₆N₂O₄ • HCl
FW: 296.8
Chemical Purity: ≥95% (Midodrine)
Deuterium Incorporation: ≥99% deuterated forms (d₁-d₆); ≤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

Midodrine-d₆ (hydrochloride) is intended for use as an internal standard for the quantification of midodrine (Item No. 17349) 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).

Midodrine-d₆ (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the midodrine-d₆ (hydrochloride) in the solvent of choice, which should be purged with an inert gas. Midodrine-d₆ (hydrochloride) is soluble in methanol and DMSO.

Description

Midodrine is a prodrug form of the α₁-adrenergic receptor (α₁-AR) agonist desglymidodrine.¹ It is converted to desglymidodrine by enzymatic hydrolysis. Midodrine (5 mg/kg) increases mean arterial pressure (MAP) and decreases heart rate in normotensive rats.² Formulations containing midodrine have been used in the treatment of orthostatic hypotension.

References

1. McClellan, K.J., Wiseman, L.R., and Wilde, M.I. Midodrine. A review of its therapeutic use in the management of orthostatic hypotension. *Drugs Aging* **12**(1), 76-86 (1998).
2. Dabasaki, T., Shimojo, M., Ishikawa, H., et al. Anti-hypotensive effects of M6434, an orally active α₁-adrenoceptor agonist, in rats. *Jpn. J. Pharmacol.* **59**(2), 145-150 (1992).

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, 11/15/2022

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