

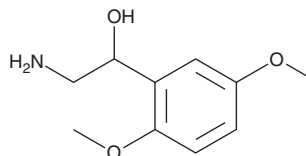
PRODUCT INFORMATION



Desglymidodrine

Item No. 35608

CAS Registry No.: 3600-87-1
Formal Name: α -(aminomethyl)-2,5-dimethoxy-benzenemethanol
Synonym: ST 1059
MF: $C_{10}H_{15}NO_3$
FW: 197.2
Purity: $\geq 98\%$
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

Desglymidodrine is supplied as a solid. A stock solution may be made by dissolving the desglymidodrine in the solvent of choice, which should be purged with an inert gas. Desglymidodrine is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of desglymidodrine in these solvents is approximately 2 and 1 mg/ml, respectively.

Description

Desglymidodrine is the active metabolite of the prodrug midodrine (Item No. 17349).¹ It is formed from midodrine by enzymatic hydrolysis.² Desglymidodrine is an agonist of α_1 -adrenergic receptors (α_1 -ARs).¹ It binds to α_1 -ARs (K_i s = 9.77, 9.55, and 8.51 μM for α_{1A} , α_{1B} , and α_{1D} -ARs, respectively) and increases intraurethral pressure (IUP) and mean arterial pressure (MAP) by 5 and 20 mm Hg, respectively, in anesthetized dogs at doses of 205 and 250 nmol/kg, respectively.

References

1. Altenbach, R.J., Khilevich, A., Kolasa, T., *et al.* Synthesis and structure-activity studies on *N*-[5-(1*H*-imidazol-4-yl)-5,6,7,8-tetrahydro-1-naphthalenyl]methanesulfonamide, an imidazole-containing α_{1A} -adrenoceptor agonist. *J. Med. Chem.* **47**(12), 3220-3235 (2004).
2. 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).

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, 12/12/2022

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