

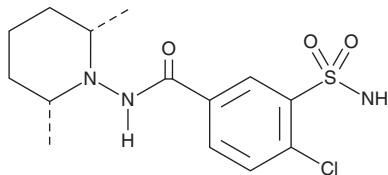
PRODUCT INFORMATION



Clopamide

Item No. 19356

CAS Registry No.: 636-54-4
Formal Name: *rel*-3-(aminosulfonyl)-4-chloro-N-(2*R*,6*S*-dimethyl-1-piperidiny)-benzamide
MF: C₁₄H₂₀ClN₃O₃S
FW: 345.8
Purity: ≥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

Clopamide is supplied as a solid. A stock solution may be made by dissolving the clopamide in the solvent of choice, which should be purged with an inert gas. Clopamide is soluble in organic solvents such as ethanol and DMSO. The solubility of clopamide in these solvents is approximately 10 and 50 mM, respectively.

Description

Clopamide is a thiazide-like diuretic that inhibits sodium/chloride cotransporters in the proximal convoluted tubule of the kidney, which prevents the reabsorption of sodium and chloride.^{1,2} This compound is intended for research purposes only.

References

1. Odland, B. and Lönnerholm, G. Renal tubular secretion and effects of chlorothiazide, hydrochlorothiazide and clopamide: A study in the avian kidney. *Acta. Pharmacol. Toxicol (Copenh)* **51(3)**, 187-197 (1982).
2. McNeil, J.J., Conway, E.L., Drummer, O.H., *et al.* Clopamide: Plasma concentrations and diuretic effect in humans. *Clin. Pharmacol. Ther.* **42(3)**, 299-304 (1987).

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, 10/25/2022

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