

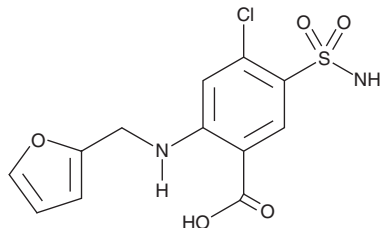
# PRODUCT INFORMATION



## Furosemide

Item No. 17273

**CAS Registry No.:** 54-31-9  
**Formal Name:** 5-(aminosulfonyl)-4-chloro-2-[(2-furanylmethyl)amino]-benzoic acid  
**Synonym:** Frusemide  
**MF:** C<sub>12</sub>H<sub>11</sub>ClN<sub>2</sub>O<sub>5</sub>S  
**FW:** 330.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 234, 273, 343 nm  
**Supplied as:** A crystalline 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

Furosemide is supplied as a crystalline solid. A stock solution may be made by dissolving the furosemide in the solvent of choice, which should be purged with an inert gas. Furosemide is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of furosemide in ethanol is approximately 10 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Furosemide is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, furosemide should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Furosemide has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Furosemide is a loop diuretic and an inhibitor of the Na<sup>+</sup>/K<sup>+</sup>/2Cl<sup>-</sup> (NKCC) cotransporters, NKCC1 and NKCC2 (K<sub>s</sub> = ~10 μM for both).<sup>1,2</sup> *In vivo*, furosemide (0.1 mg/kg, p.o.) increases diuresis in beagle dogs.<sup>3</sup> Furosemide (30 mg/kg) reduces ventricular collagen deposition and fibrosis in a rat model of dilated cardiomyopathy.<sup>4</sup> It is also an inhibitor of carbonic anhydrase I (CAI), CAII, and CAIII (K<sub>s</sub> = 0.052-0.065 μM) and organic ion transporter 1 (OAT1; K<sub>i</sub> = 9.5 μM), as well as a GABA<sub>A</sub> receptor antagonist.<sup>5-7</sup> This product is also available as an analytical reference standard (Item No. 26298).

### References

1. Gillen, C.M., Brill, S., Payne, J.A., et al. *J. Biol. Chem.* **271**(27), 16237-16244 (1996).
2. Markadieu, N. and Delpire, E. *Pflugers. Arch.* **466**(1), 91-105 (2014).
3. Potter, B.M., Ames, M.K., Hess, A., et al. *J. Vet. Cardiol.* **26**, 51-62 (2019).
4. Watanabe, K., Sreedhar, R., Thandavarayan, R.A., et al. *Biofactors* **43**(2), 187-194 (2017).
5. Temperini, C., Cecchi, A., Scozzafava, A., et al. *Bioorg. Med. Chem. Lett.* **18**(8), 2567-2573 (2008).
6. Rajan, S.T., Kumar, M.K., and Makrishna, M. *PCT/IN2014/000406* (2014).
7. Siess, W., Siegel, F.L., and Lapetina, E.G. *Biochemica et Biophysica Acta* **801**, 265-276 (1984).

#### 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, 11/17/2022

#### 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