

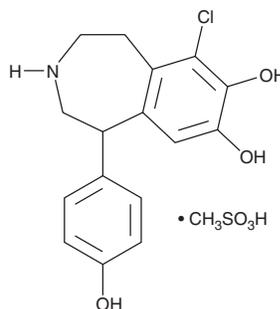
# PRODUCT INFORMATION



## Fenoldopam (mesylate)

Item No. 17629

**CAS Registry No.:** 67227-57-0  
**Formal Name:** 6-chloro-2,3,4,5-tetrahydro-1-(4-hydroxyphenyl)-1H-3-benzazepine-7,8-diol, monomethanesulfonate  
**Synonym:** SKF 82526J  
**MF:** C<sub>16</sub>H<sub>16</sub>ClNO<sub>3</sub> • CH<sub>3</sub>SO<sub>3</sub>H  
**FW:** 401.9  
**Purity:** ≥95%  
**Supplied as:** A crystalline solid  
**Storage:** -20°C  
**Stability:** ≥2 years



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

### Laboratory Procedures

Fenoldopam (mesylate) is supplied as a crystalline solid. A stock solution may be made by dissolving the fenoldopam (mesylate) in the solvent of choice, which should be purged with an inert gas. Fenoldopam (mesylate) is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of fenoldopam (mesylate) in these solvents is approximately 0.25, 1, and 2 mg/ml, respectively.

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

### Description

Fenoldopam is an agonist of dopamine D<sub>1A</sub> (D1R) and D<sub>1B</sub> (D5R) receptors (K<sub>d</sub>s = 17 and 11 nM, respectively).<sup>1,2</sup> Fenoldopam is used to study the roles of these receptors, in cells and *in vivo*, and to alter hemodynamic properties, including hypertension, in animals.<sup>3,4</sup>

### References

- Carey, R.M., Stote, R.M., Dubb, J.W., *et al.* Selective peripheral dopamine-1 receptor stimulation with fenoldopam in human essential hypertension. *J. Clin. Invest.* **74**(6), 2198-2207 (1984).
- Tiberi, M., Jarvie, K.R., Silvia, C., *et al.* Cloning, molecular characterization, and chromosomal assignment of a gene encoding a second D<sub>1</sub> dopamine receptor subtype: Differential expression pattern in rat brain compared with the D<sub>1A</sub> receptor. *Proc. Natl. Acad. Sci. USA* **88**(17), 7491-7495 (1991).
- Gildea, J.J., Shah, I.T., Van Sciver, R., *et al.* The cooperative roles of the dopamine receptors, D<sub>1</sub>R and D<sub>5</sub>R, on the regulation of renal sodium transport. *Kidney Int.* **86**(1), 118-126 (2014).
- Swanson, T.A., Conte, T., Deeley, B., *et al.* Hemodynamic correlates of drug-induced vascular injury in the rat using high-frequency ultrasound imaging. *Toxicol. Pathol.* **42**(4), 784-791 (2014).

#### 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, 06/12/2024

#### 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