

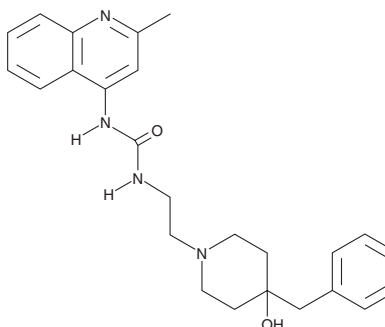
# PRODUCT INFORMATION



## Palosuran

Item No. 23460

**CAS Registry No.:** 540769-28-6  
**Formal Name:** N-[2-[4-hydroxy-4-(phenylmethyl)-1-piperidiny]ethyl]-N'-(2-methyl-4-quinoliny)-urea  
**Synonym:** ACT058362  
**MF:** C<sub>25</sub>H<sub>30</sub>N<sub>4</sub>O<sub>2</sub>  
**FW:** 418.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 216, 234, 338 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

Palosuran is supplied as a crystalline solid. A stock solution may be made by dissolving the palosuran in the solvent of choice. Palosuran is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of palosuran in ethanol is approximately 0.5 mg/ml and approximately 10 mg/ml in DMSO and DMF.

Palosuran is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, palosuran should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Palosuran 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

Palosuran is an antagonist of the urotensin II receptor (UTR; IC<sub>50</sub>s = 3.6 and 1,475 nM for CHO cell membranes expressing human and rat recombinant receptors, respectively).<sup>1</sup> It inhibits contraction of isolated rat aortic rings induced by urotensin II (Item Nos. 24711 | 24753) in a concentration-dependent manner. Palosuran (300 mg/kg twice per day) reduces plasma endothelin-1 (Item No. 24127), urotensin II, and TGF-β1 levels, as well as the main pulmonary arterial pressure and right ventricular hypertrophy index in a rat model of pulmonary arterial hypertension (PAH) induced by monocrotaline (Item No. 16666).<sup>2</sup> Palosuran (300 mg/kg per day) reduces serum glucose, cholesterol, and triglyceride levels and increases survival in a rat model of diabetes induced by streptozotocin (STZ; Item No. 13104).<sup>3</sup> It also decreases albuminuria and global kidney lesion scores in STZ-induced diabetic rats.

### References

1. Clozel, M., Binkert, C., Birker-Robaczewska, M., *et al.* Pharmacology of the urotensin-II receptor antagonist palosuran (ACT-058362; 1-[2-(4-benzyl-4-hydroxy-piperidin-1-yl)-ethyl]-3-(2-methyl-quinolin-4-yl)-urea sulfate salt): First demonstration of a pathophysiological role of the urotensin system. *J. Pharmacol. Exp. Ther.* **311**(1), 204-212 (2004).
2. Onat, A.M., Pehlivan, Y., Turkbeyler, I.H., *et al.* Urotensin inhibition with palosuran could be a promising alternative in pulmonary arterial hypertension. *Inflammation* **36**(2), 405-412 (2013).
3. Clozel, M., Hess, P., Qiu, C., *et al.* The urotensin-II receptor antagonist palosuran improves pancreatic and renal function in diabetic rats. *J. Pharmacol. Exp. Ther.* **316**(3), 1115-1121 (2006).

#### 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, 12/22/2022

#### 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](http://WWW.CAYMANCHEM.COM)