

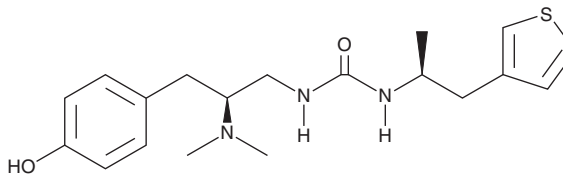
# PRODUCT INFORMATION



## PZM21

Item No. 20576

**CAS Registry No.:** 1997387-43-5  
**Formal Name:** N-[(2S)-2-(dimethylamino)-3-(4-hydroxyphenyl)propyl]-N'-[(1S)-1-methyl-2-(3-thienyl)ethyl]-urea  
**MF:** C<sub>19</sub>H<sub>27</sub>N<sub>3</sub>O<sub>2</sub>S  
**FW:** 361.5  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 229, 281 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

PZM21 is supplied as a crystalline solid. A stock solution may be made by dissolving the PZM21 in the solvent of choice, which should be purged with an inert gas. PZM21 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of PZM21 in ethanol and DMF is approximately 10 mg/ml and approximately 1 mg/ml in DMSO.

PZM21 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PZM21 should first be dissolved in ethanol and then diluted with the aqueous buffer of choice. PZM21 has a solubility of approximately 0.2 mg/ml in a 1:4 solution of ethanol:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

PZM21 is a  $\mu$ -opioid receptor agonist ( $K_i = 1.1$  nM).<sup>1</sup> It is selective for  $\mu$ - over  $\kappa$ - and  $\delta$ -opioid receptors ( $K_{\mu} = 18$  and  $506$  nM, respectively). PZM21 induces  $G_{i/o}$  signaling ( $EC_{50} = 4.6$  nM in HEK293T cells expressing human receptors) but has no detectable activity in a  $\beta$ -arrestin2 recruitment assay. *In vivo*, PZM21 (10, 20, and 40 mg/kg) induces analgesia in the hot plate test and reduces formalin-induced paw licking in mice when administered at a dose of 40 mg/kg.

### Reference

1. Manglik, A., Lin, H., Aryal, D.K., *et al.* Structure-based discovery of opioid analgesics with reduced side effects. *Nature* **537(7619)**, 185-190 (2016).

#### 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, 10/07/2022

#### 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