

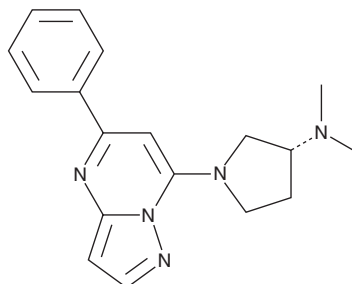
# PRODUCT INFORMATION



## (R)-ZINC3573

Item No. 41312

**CAS Registry No.:** 2089389-15-9  
**Formal Name:** N,N-dimethyl-1-(5-phenylpyrazolo[1,5-a]pyrimidin-7-yl)-3R-pyrrolidinamine  
**Synonym:** (R)-ZINC72453573  
**MF:** C<sub>18</sub>H<sub>21</sub>N<sub>5</sub>  
**FW:** 307.4  
**Purity:** ≥95%  
**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

(R)-ZINC3573 is supplied as a solid. A stock solution may be made by dissolving the (R)-ZINC3573 in the solvent of choice, which should be purged with an inert gas. (R)-ZINC3573 is soluble (≥10 mg/ml) in DMSO and sparingly soluble (1-10 mg/ml) in ethanol.

### Description

(R)-ZINC3573 is an agonist of MAS-related G protein-coupled receptor family member X2 (MRGPRX2).<sup>1</sup> It induces β-arrestin recruitment in a reporter assay using MRGPRX2-expressing HTLA cells and calcium mobilization in LAD 2 mast cells (EC<sub>50</sub>s = 0.74 and 1 μM, respectively). (R)-ZINC3573 is selective for MRGPRX2 over 21 other G protein-coupled receptors (GPCRs). It induces degranulation of LAD 2 mast cells in a concentration-dependent manner.

### Reference

1. Lansu, K., Karpiak, J., Liu, J., *et al.* *In silico* design of novel probes for the atypical opioid receptor MRGPRX2. *Nat. Chem. Biol.* **13**(5), 529-536 (2017).

#### 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, 09/09/2024

#### 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