

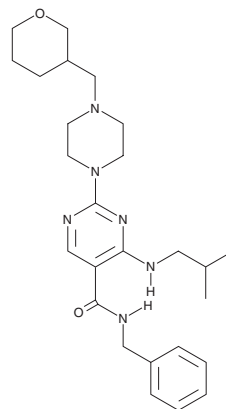
# PRODUCT INFORMATION



**PF-4840154**

Item No. 28915

**CAS Registry No.:** 1332708-14-1  
**Formal Name:** 4-[(2-methylpropyl)amino]-N-(phenylmethyl)-2-[4-[(tetrahydro-2H-pyran-3-yl)methyl]-1-piperazinyl]-5-pyrimidinecarboxamide  
**MF:** C<sub>26</sub>H<sub>38</sub>N<sub>6</sub>O<sub>2</sub>  
**FW:** 466.6  
**Purity:** ≥95%  
**UV/Vis.:** λ<sub>max</sub>: 234, 265, 311 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

PF-4840154 is supplied as a crystalline solid. A stock solution may be made by dissolving the PF-4840154 in the solvent of choice, which should be purged with an inert gas. PF-4840154 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of PF-4840154 in these solvents is approximately 10, 20, and 30 mg/ml, respectively.

PF-4840154 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, PF-4840154 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. PF-4840154 has a solubility of approximately 0.1 mg/ml in a 1:5 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

PF-4840154 is a transient receptor potential ankyrin 1 (TRPA1) agonist.<sup>1</sup> It induces calcium influx into HEK293 cells expressing recombinant human or rat TRPA1 (EC<sub>50</sub>s = 23 and 97 nM, respectively). Intraplantar injection of PF-4840154 (30 nmol) increases paw licking time in mice in a TRPA1-dependent manner.

## Reference

1. Ryckmans, T., Aubdool, A.A., Bodkin, J.V., *et al.* Design and pharmacological evaluation of PF-4840154, a non-electrophilic reference agonist of the TrpA1 channel. *Bioorg. Med. Chem.* **21(16)**, 4857-4859 (2011).

### 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, 12/08/2022

## 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