

# PRODUCT INFORMATION



**SB-452533**

Item No. 37379

**CAS Registry No.:** 459429-39-1

**Formal Name:** N-(2-bromophenyl)-N'-[2-[ethyl(3-methylphenyl)amino]ethyl]-urea

**MF:** C<sub>18</sub>H<sub>22</sub>BrN<sub>3</sub>O

**FW:** 376.3

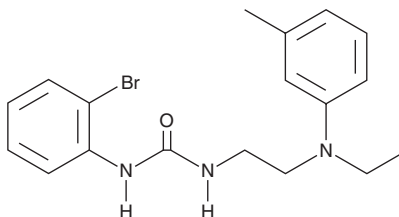
**Purity:** ≥95%

**UV/Vis.:** λ<sub>max</sub>: 252 nm

**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

SB-452533 is supplied as a solid. A stock solution may be made by dissolving the SB-452533 in the solvent of choice, which should be purged with an inert gas. SB-452533 is soluble in methanol and DMSO.

## Description

SB-452533 is a transient receptor potential vanilloid 1 (TRPV1) antagonist.<sup>1</sup> It inhibits heat-induced activation of TRPV1 in HEK293 cells expressing the human channel when used at a concentration of 1 μM.

## Reference

1. Rami, H.K., Thompson, M., Wyman, P., *et al.* Discovery of small molecule antagonists of TRPV1. *Bioorg. Med. Chem. Lett.* **14**(14), 3631-3634 (2004).

### 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, 09/28/2023

## 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)