

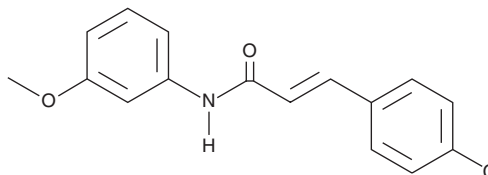
# PRODUCT INFORMATION



**SB-366791**

Item No. 11019

**CAS Registry No.:** 472981-92-3  
**Formal Name:** 3-(4-chlorophenyl)-N-(3-methoxyphenyl)-2-propenamamide  
**MF:** C<sub>16</sub>H<sub>14</sub>ClNO<sub>2</sub>  
**FW:** 287.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 297 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

SB-366791 is supplied as a crystalline solid. A stock solution may be made by dissolving the SB-366791 in the solvent of choice, which should be purged with an inert gas. SB-366791 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of SB-366791 in these solvents is approximately 10 mg/ml.

SB-366791 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, SB-366791 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. SB-366791 has a solubility of approximately 0.15 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

The transient receptor potential vanilloid type 1 (TRPV1) receptor is a nonselective cation channel gated by noxious heat, protons, and capsaicin. TRPV1 is present in primary sensory neurons and in both central and peripheral sensory nerve terminals and plays a role in thermal and mechanical hyperalgesia. SB-366791 is a selective TRPV1 antagonist that is widely used in pain research.<sup>1</sup> In cultured trigeminal ganglion neurons, SB-366791 inhibits capsaicin-evoked Ca<sup>2+</sup> influx with an IC<sub>50</sub> value of 0.7 μM.<sup>1</sup> A 1 nmol injection of SB-366791 in mouse paw reduces capsaicin-induced nociceptive responses in a model of pain.<sup>2</sup> Intrathecal pretreatment of 10 μl SB-366791 in rats suppresses tolerance to the analgesic effects of chronic morphine administration.<sup>3</sup> Co-administration of morphine and SB-366791 at 0.1 mg/kg has potent analgesic effects in a mouse model of bone cancer pain.<sup>4</sup>

## References

1. Varga, A., Németh, J., Szabó, A., *et al.* Effects of the novel TRPV1 receptor antagonist SB366791 in vitro and in vivo in the rat. *Neurosci. Lett.* **385(2)**, 137-142 (2005).
2. Andrade, E.L., Luiz, A.P., Ferreira, J., *et al.* Pronociceptive response elicited by TRPA1 receptor activation in mice. *Neuroscience* **152(2)**, 511-520 (2008).
3. Chen, Y., Geis, C., and Sommer, C. Activation of TRPV1 contributes to morphine tolerance: Involvement of the mitogen-activated protein kinase signaling pathway. *J. Neurosci.* **28(22)**, 5836-5845 (2008).
4. Niiyama, Y., Kawamata, T., Yamamoto, J., *et al.* SB366791, a TRPV1 antagonist, potentiates analgesic effects of systemic morphine in a murine model of bone cancer pain. *Br. J. Anaesth.* **102(2)**, 251-258 (2009).

### 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, 02/21/2024

## 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