

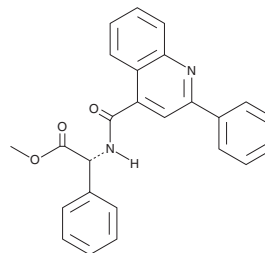
# PRODUCT INFORMATION



**SB-218795**

Item No. 42373

**CAS Registry No.:** 174635-53-1  
**Formal Name:** αR-[[[(2-phenyl-4-quinolinyl)carbonyl]amino]-benzeneacetic acid, methyl ester  
**MF:** C<sub>25</sub>H<sub>20</sub>N<sub>2</sub>O<sub>3</sub>  
**FW:** 396.4  
**Purity:** ≥98%  
**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-218795 is supplied as a solid. A stock solution may be made by dissolving the SB-218795 in the solvent of choice, which should be purged with an inert gas. SB-218795 is sparingly soluble (1-10 mg/ml) in ethanol and soluble (≥10 mg/ml) in DMSO.

## Description

SB-218795 is a neurokinin-3 (NK<sub>3</sub>) receptor antagonist (K<sub>i</sub> = 0.013 μM).<sup>1</sup> It is selective for NK<sub>3</sub> over NK<sub>1</sub> and NK<sub>2</sub> receptors (K<sub>s</sub> = >100 and 1.221 μM, respectively). SB-218795 inhibits senktide-induced contractions in isolated rabbit iris sphincter muscle preparations (K<sub>b</sub> = 43 nM). It also inhibits substance P-induced contraction of longitudinal and circular smooth muscles isolated from rabbit duodenum, jejunum, or ileum when used at a concentration of 1 μM.<sup>2</sup>

## References

- Giardina, G.A., Sarau, H.M., Farina, C., *et al.* Discovery of a novel class of selective non-peptide antagonists for the human neurokinin-3 receptor. 1. Identification of the 4-quinolinecarboxamide framework. *J. Med. Chem.* **40(12)**, 1794-1807 (1997).
- Valero, M.S., Fagundes, D.S., Grasa, L., *et al.* Contractile effect of tachykinins on rabbit small intestine. *Acta Pharmacol. Sin.* **32(4)**, 487-494 (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, 12/18/2024

## 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