

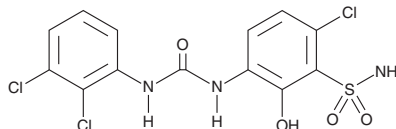
# PRODUCT INFORMATION



**SB-332235**

Item No. 32869

**CAS Registry No.:** 276702-15-9  
**Formal Name:** 6-chloro-3-[[[(2,3-dichlorophenyl)amino]carbonyl]amino]-2-hydroxy-benzenesulfonamide  
**MF:** C<sub>13</sub>H<sub>10</sub>Cl<sub>3</sub>N<sub>3</sub>O<sub>4</sub>S  
**FW:** 410.7  
**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

SB-332235 is supplied as a solid. A stock solution may be made by dissolving the SB-332235 in the solvent of choice, which should be purged with an inert gas. SB-332235 is soluble in the organic solvent DMSO.

## Description

SB-332235 is an antagonist of chemokine (C-X-C motif) receptor 2 (CXCR2; IC<sub>50</sub> = 9.3 nM in CHO cell membranes expressing the recombinant human receptor).<sup>1</sup> It is selective for CXCR2 over CXCR1 (IC<sub>50</sub> = 9,633 nM). SB-332235 (10 ng/ml) inhibits migration of isolated human peripheral blood monocytes induced by chemokine (C-X-C motif) ligand 1 (CXCL1), CXCL5, or CXCL8 by 92, 73, and 50%, respectively.<sup>2</sup> It decreases cigarette smoke-induced increases in bronchoalveolar lavage fluid (BALF) neutrophil infiltration when administered at a dose of 1 mg/kg.<sup>3</sup> SB-332235 (1 mg/kg) reduces increases in neuronal levels of 4-hydroxy nonenal (4-HNE; Item No. 32100) induced by amyloid-β (1-42) (Aβ42; Item No. 20574) in a rat model of Alzheimer's disease.<sup>4</sup>

## References

1. Podolin, P.L., Bolognese, B.J., Foley, J.J., *et al.* A potent and selective nonpeptide antagonist of CXCR2 inhibits acute and chronic models of arthritis in the rabbit. *J. Immunol.* **169(11)**, 6435-6444 (2002).
2. Traves, S.L., Smith, S.J., Barnes, P.J., *et al.* Specific CXC but not CC chemokines cause elevated monocyte migration in COPD: A role for CXCR<sub>2</sub>. *J. Leukoc. Biol.* **76(2)**, 441-450 (2004).
3. Stevenson, C.S., Coote, K., Webster, R., *et al.* Characterization of cigarette smoke-induced inflammatory and mucus hypersecretory changes in rat lung and the role of CXCR2 ligands in mediating this effect. *Am. J. Physiol. Lung Cell. Mol. Physiol.* **288(3)**, L514-L522 (2005).
4. Ryu, J.K., Cho, T., Choi, H.B., *et al.* Pharmacological antagonism of interleukin-8 receptor CXCR2 inhibits inflammatory reactivity and is neuroprotective in an animal model of Alzheimer's disease. *J. Neuroinflammation* **12**, 144 (2015).

### 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, 08/15/2023

## 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