

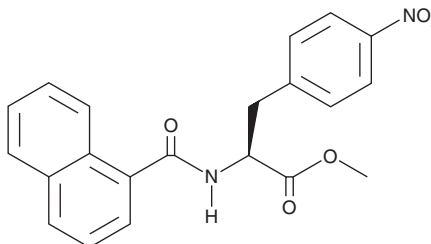
PRODUCT INFORMATION



SB-328437

Item No. 29677

CAS Registry No.: 247580-43-4
Formal Name: N-(1-naphthalenylcarbonyl)-4-nitro-L-phenylalanine, methyl ester
MF: C₂₁H₁₈N₂O₅
FW: 378.4
Purity: ≥98%
UV/Vis.: λ_{max}: 221, 278 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-328437 is supplied as a crystalline solid. A stock solution may be made by dissolving the SB-328437 in the solvent of choice, which should be purged with an inert gas. SB-328437 is soluble in the organic solvent DMSO at a concentration of approximately 100 mM.

Description

SB-328437 is a nonpeptide chemokine (C-C motif) receptor 3 (CCR3) antagonist (IC₅₀ = 4.5 nM in a radioligand binding assay).¹ It is selective for CCR3 over chemokine (C-X-C motif) receptor 1 (CXCR1), CXCR2, CCR7, complement 3a receptor (C3aR), and C5aR (IC₅₀s = >27, >27, >33, >10, and >10 μM, respectively). SB-328437 (3.3-1,000 nM) inhibits calcium mobilization induced chemokine (C-C motif) ligand 11 (CCL11), CCL24, and CCL13 in RBL-2H3-CCR3 cells and eosinophils. *In vivo*, SB-328437 reduces pulmonary *Ccr3* mRNA expression, the number of eosinophils and monocytes in bronchoalveolar lavage fluid (BALF), and pulmonary IL-4 and TNF-α levels in a mouse model of asthma.²

References

1. White, J.R., Lee, J.M., Dede, K., *et al.* Identification of potent, selective non-peptide CC chemokine receptor-3 antagonist that inhibits eotaxin-, eotaxin-2-, and monocyte chemotactic protein-4-induced eosinophil migration. *J. Biol. Chem.* **275**(47), 36626-36631 (2000).
2. Wang, T., Zhang, J.-Y., Wang, X.-D., *et al.* Effects of CCR3(chemokine receptor-3) on Muc5ac in airway of asthmatic mice. *Zhongguo Mianyixue Zazhi* **31**(8), 1019-1023 (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, 12/15/2022

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