

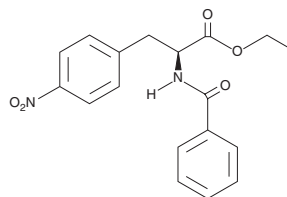
PRODUCT INFORMATION



SB 297006

Item No. 11964

CAS Registry No.: 58816-69-6
Formal Name: N-benzoyl-4-nitro-L-phenylalanine, ethyl ester
MF: C₁₈H₁₈N₂O₅
FW: 342.4
Purity: ≥99%
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 297006 is supplied as a solid. A stock solution may be made by dissolving the SB 297006 in the solvent of choice, which should be purged with an inert gas. SB 297006 is soluble in organic solvents such as ethanol and DMSO. The solubility of SB 297006 in these solvents is approximately 25 and 10 mM, respectively.

Description

SB 297006 is an antagonist of C-C chemokine receptor 3 (CCR3; IC₅₀ = 39 nM), which normally is activated by eotaxin, eotaxin-3, MCP-3, MCP-4, RANTES, and MIP-1δ.¹ It is at least 250-fold selective for CCR3 over a panel of other chemokine receptors. SB 297006 blocks CCR3-mediated calcium mobilization induced by eotaxin, eotaxin-2, and MCP-4 in transfected cells.¹ It suppresses antigen-induced accumulation of Th2 lymphocytes and eosinophils in lungs of mice when delivered subcutaneously (100 mg/kg).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. Mori, A., Ogawa, K., Someya, K., *et al.* Selective suppression of T_H2-mediated airway eosinophil infiltration by low-molecular weight CCR3 antagonists. *Int. Immunol.* **19**(8), 913-921 (2007).

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, 11/15/2022

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