

# PRODUCT INFORMATION



**SB-273005**

Item No. 34404

**CAS Registry No.:** 205678-31-5  
**Formal Name:** 2,3,4,5-tetrahydro-8-[2-[6-(methylamino)-2-pyridinyl]ethoxy]-3-oxo-2-(2,2,2-trifluoroethyl)-1H-2-benzazepine-4-acetic acid

**MF:** C<sub>22</sub>H<sub>24</sub>F<sub>3</sub>N<sub>3</sub>O<sub>4</sub>

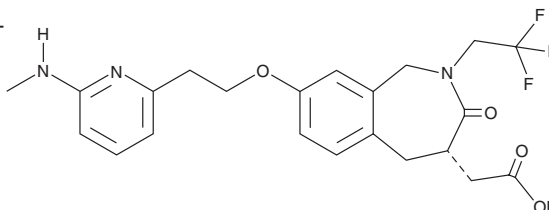
**FW:** 451.4

**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-273005 is supplied as a solid. A stock solution may be made by dissolving the SB-273005 in the solvent of choice, which should be purged with an inert gas. SB-273005 is slightly soluble (0.1-1 mg/ml) in acetonitrile.

## Description

SB-273005 is a nonpeptide antagonist of  $\alpha_V\beta_3$  and  $\alpha_V\beta_5$  integrins ( $K_i$ s = 1.2 and 0.3 nM, respectively).<sup>1</sup> *In vivo*, SB-273005 (30 and 60 mg/kg) reduces paw edema and increases bone density in a rat model of adjuvant-induced arthritis. It also reduces the number of T helper 2 (Th2) cells and production of IL-10 in pregnant mice when administered at a dose of 3 mg/kg.<sup>2</sup>

## References

1. Badger, A.M., Blake, S., Kapadia, R., *et al.* Disease-modifying activity of SB 273005, an orally active, nonpeptide  $\alpha_V\beta_3$  (vitronectin receptor) antagonist, in rat adjuvant-induced arthritis. *Arthritis Rheum.* **44**(1), 128-137 (2001).
2. Wang, S., Yang, J., Wang, C., *et al.* SB-273005, an antagonist of  $\alpha_V\beta_3$  integrin, reduces the production of Th2 cells and cytokine IL-10 in pregnant mice. *Exp. Ther. Med.* **7**(6), 1677-1682 (2014).

### 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, 05/05/2025

## 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