

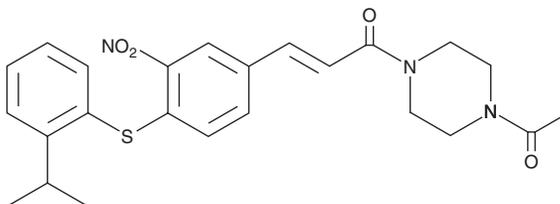
PRODUCT INFORMATION



A286982

Item No. 27646

CAS Registry No.: 280749-17-9
Formal Name: (2E)-1-(4-acetyl-1-piperazinyl)-3-[4-[[2-(1-methylethyl)phenyl]thio]-3-nitrophenyl]-2-propen-1-one
MF: C₂₄H₂₇N₃O₄S
FW: 453.6
Purity: ≥98%
UV/Vis.: λ_{max}: 316 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

A286982 is supplied as a crystalline solid. A stock solution may be made by dissolving the A286982 in the solvent of choice, which should be purged with an inert gas. A286982 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of A286982 in these solvents is approximately 30 mg/ml.

A286982 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, A286982 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. A286982 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

A286982 is an inhibitor that blocks the integrin-ligand interaction between leukocyte function-associated antigen-1 (LFA-1) and intercellular adhesion molecule-1 (ICAM-1; IC₅₀ = 35 nM).¹ It inhibits adhesion of JY-8 cells to ICAM-1-coated microtiter plates (IC₅₀ = 44 nM).

Reference

1. Liu, G., Link, J.T., Pei, Z., *et al.* Discovery of novel p-arylthio cinnamides as antagonists of leukocyte function-associated antigen-1/intracellular adhesion molecule-1 interaction. 1. Identification of an additional binding pocket based on an anilino diaryl sulfide lead. *J. Med. Chem.* **43(21)**, 4025-4040 (2000).

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, 12/02/2022

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