

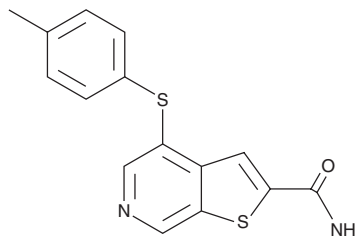
PRODUCT INFORMATION



A-205804

Item No. 21252

CAS Registry No.: 251992-66-2
Formal Name: 4-[(4-methylphenyl)thio]-thieno[2,3-c]pyridine-2-carboxamide
MF: C₁₅H₁₂N₂OS₂
FW: 300.4
Purity: ≥95%
UV/Vis.: λ_{max}: 229, 340 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

A-205804 is supplied as a crystalline solid. A stock solution may be made by dissolving the A-205804 in the solvent of choice, which should be purged with an inert gas. A-205804 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of A-205804 in these solvents is approximately 15 and 25 mg/ml, respectively.

A-205804 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, A-205804 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. A-205804 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

A-205804 is a potent inhibitor of E-selectin and ICAM-1 expression in human vascular endothelial cells stimulated with TNF-α (IC₅₀s = 20 and 25 nM, respectively).¹ It does not block expression of VCAM-1 (IC₅₀ > 1,000 nM).

Reference

1. Stewart, A.O., Bhatia, P.A., McCarty, C.M., *et al.* Discovery of inhibitors of cell adhesion molecule expression in human endothelial cells. 1. Selective inhibition of ICAM-1 and E-selectin expression. *J. Med. Chem.* **44**(6), 988-1002 (2001).

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, 10/05/2022

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