

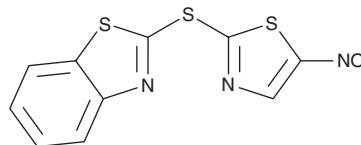
PRODUCT INFORMATION



BI-87G3

Item No. 42045

CAS Registry No.: 2207-44-5
Formal Name: 2-[(5-nitro-2-thiazolyl)thio]-benzothiazole
Synonyms: JNK Inhibitor XI, c-Jun N-terminal Kinase Inhibitor XI
MF: C₁₀H₅N₃O₂S₃
FW: 295.4
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

BI-87G3 is supplied as a solid. A stock solution may be made by dissolving the BI-87G3 in the solvent of choice, which should be purged with an inert gas. BI-87G3 is sparingly soluble (1-10 mg/ml) in DMSO.

Description

BI-87G3 is an inhibitor of JNK1 (IC₅₀ = 1.804 μM).¹ It is selective for JNK1 over p38α MAPK, Akt, furin, and lethal factor (IC₅₀s = >100, >100, >100, and >50 μM, respectively). BI-87G3 inhibits TNF-α-induced JNK1 phosphorylation in HeLa cells (IC₅₀ = ~15 μM).

Reference

1. De, S.K., Stebbins, J.L., Chen, L.-H., *et al.* Design, synthesis, and structure-activity relationship of substrate competitive, selective, and in vivo active triazole and thiadiazole inhibitors of the c-Jun N-terminal kinase. *J. Med. Chem.* **52**(7), 1943-1952 (2009).

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, 10/22/2024

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