

PRODUCT INFORMATION

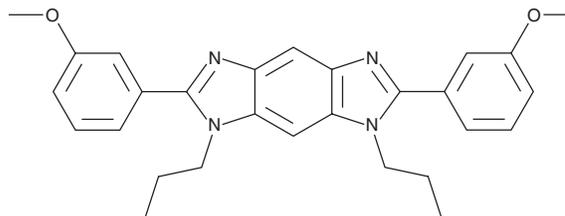


STAT3-IN-12

Item No. 44651

CAS Registry No.: 2980758-31-2
Formal Name: 2,6-bis(3-methoxyphenyl)-1,7-dipropyl-1,7-dihydrobenzo[1,2-d:4,5-d']diimidazole
Synonyms: Signal Transducer and Activator of Transcription 3-IN-12, Signal Transducer and Activator of Transcription 3 Inhibitor 12, STAT3 Inhibitor 12

MF: C₂₈H₃₀N₄O₂
FW: 454.6
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

STAT3-IN-12 is supplied as a solid. A stock solution may be made by dissolving the STAT3-IN-12 in the solvent of choice, which should be purged with an inert gas. STAT3-IN-12 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

STAT3-IN-12 is an inhibitor of STAT3 (IC₅₀ = 2.61 μM).¹ It induces apoptosis in HepG2 hepatocellular carcinoma and EC109 esophageal carcinoma cells when used at concentrations of 5, 10, and 20 μM. *In vivo*, STAT3-IN-12 (20 and 40 mg/kg) reduces tumor volume in a HepG2 mouse xenograft model.

Reference

1. Liu, Y.-C., Yang, Y.-D., Liu, W.-Q., *et al.* Benzobis(imidazole) derivatives as STAT3 signal inhibitors with antitumor activity. *Bioorg. Med. Chem.* **65**, 116757 (2022).

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, 12/08/2025

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