

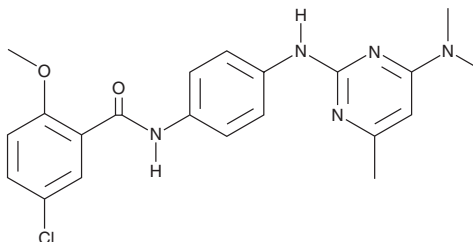
PRODUCT INFORMATION



TG53

Item No. 42514

CAS Registry No.: 946369-04-6
Formal Name: 5-chloro-N-[4-[[4-(dimethylamino)-6-methyl-2-pyrimidinyl]amino]phenyl]-2-methoxy-benzamide
MF: C₂₁H₂₂ClN₅O₂
FW: 411.9
Purity: ≥98%
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

TG53 is supplied as a crystalline solid. A stock solution may be made by dissolving the TG53 in the solvent of choice, which should be purged with an inert gas. TG53 is slightly soluble (0.1-1 mg/ml) in ethanol and soluble (≥10 mg/ml) in DMSO.

Description

TG53 is an inhibitor of the protein-protein interaction between transglutaminase 2 (TG2) and fibronectin (IC₅₀ = 10 μM in a cell-free assay).¹ It inhibits the adhesion of SKOV3 ovarian cancer cells to fibronectin-coated wells when used at a concentration of 25 μM. TG53 (10 μM) inhibits migration and invasion in a scratch assay using SKOV3 cells.

Reference

1. Yakubov, B., Chen, L., Belkin, A.M., *et al.* Small molecule inhibitors target the tissue transglutaminase and fibronectin interaction. *PLoS One* **9**(2), e89285 (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, 02/07/2025

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