

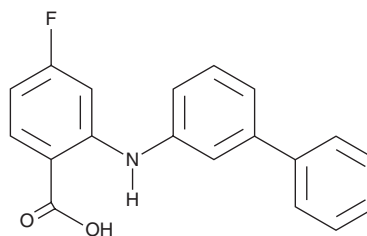
PRODUCT INFORMATION



LM-41

Item No. 45312

CAS Registry No.: 2996821-30-6
Formal Name: 2-([1,1'-biphenyl]-3-ylamino)-4-fluoro-benzoic acid
MF: C₁₉H₁₄FNO₂
FW: 307.3
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

LM-41 is supplied as a solid. A stock solution may be made by dissolving the LM-41 in the solvent of choice, which should be purged with an inert gas. LM-41 is soluble (≥10 mg/ml) in DMSO.

Description

LM-41 is a covalent TEA domain transcription factor 4 (TEAD4) inhibitor.¹ It reduces migration of MDA-MB-231 breast cancer cells when used at a concentration of 10 μM.

Reference

1. Fnaiche, A., Mélin, L., Suárez, N. G., *et al.* Development of LM-41 and AF-2112, two flufenamic acid-derived TEAD inhibitors obtained through the replacement of the trifluoromethyl group by aryl rings. **15:95**, 129488 (2023).

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, 03/16/2026

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