

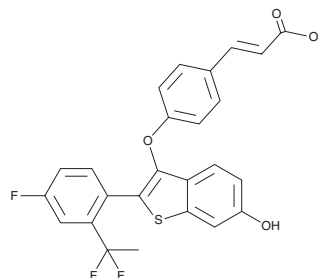
PRODUCT INFORMATION



LSZ-102

Item No. 41834

CAS Registry No.: 2135600-76-7
Formal Name: 3-[4-[[2E-[2-(1,1-difluoroethyl)-4-fluorophenyl]-6-hydroxybenzo[b]thien-3-yl]oxy]phenyl]-2-propenoic acid
MF: C₂₅H₁₇F₃O₄S
FW: 470.5
Purity: ≥95%
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

LSZ-102 is supplied as a solid. A stock solution may be made by dissolving the LSZ-102 in the solvent of choice, which should be purged with an inert gas. LSZ-102 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

LSZ-102 is an estrogen receptor α (ER α) degrader (IC₅₀ = 0.2 nM).¹ It reduces 17 β -estradiol-induced transcription in a reporter assay using MCF-7 breast cancer cells expressing ER α (IC₅₀ = 6 nM). LSZ-102 decreases the proliferation of MCF-7 cells (IC₅₀ = 1.7 nM). *In vivo*, LSZ-102 (20 mg/kg per day) reduces tumor volume in an MCF-7 mouse xenograft model.

Reference

1. Tria, G.S., Abrams, T., Baird, J., *et al.* Discovery of LSZ102, a potent, orally bioavailable selective estrogen receptor degrader (SERD) for the treatment of estrogen receptor positive breast cancer. *J. Med. Chem.* **61**(7), 2837-2864 (2018).

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, 10/09/2024

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