

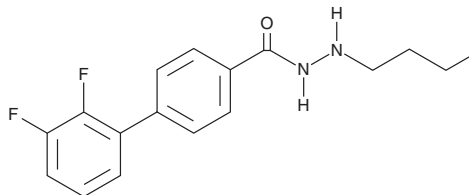
PRODUCT INFORMATION



SR 4370

Item No. 25543

CAS Registry No.: 1816294-67-3
Formal Name: 2',3'-difluoro-[1,1'-biphenyl]-4-carboxylic acid, 2-butylhydrazide
MF: C₁₇H₁₈F₂N₂O
FW: 304.3
Purity: ≥98%
UV/Vis.: λ_{max}: 261 nm
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

SR 4370 is supplied as a crystalline solid. A stock solution may be made by dissolving the SR 4370 in the solvent of choice, which should be purged with an inert gas. SR 4370 is soluble in the organic solvent DMSO.

Description

SR 4370 is an inhibitor of histone deacetylase 3 (HDAC3; IC₅₀ = 6 nM).¹ It is selective for HDAC3 over HDAC1, -2, -6, and -8 (IC₅₀s = 0.13, 0.58, 3.7, and 2.3 μM, respectively). SR 4370 decreases viability of MDA-MB-231 human breast cancer cells with an IC₅₀ value of 12.6 μM.

Reference

1. Liao, D., Roush, W.R., and Stowe, R.L. HDAC inhibitor compounds and methods of treatment. *University of Florida Research Foundation*. **WO 2015/153516 A1** (2015).

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, 10/10/2022

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