

PRODUCT INFORMATION

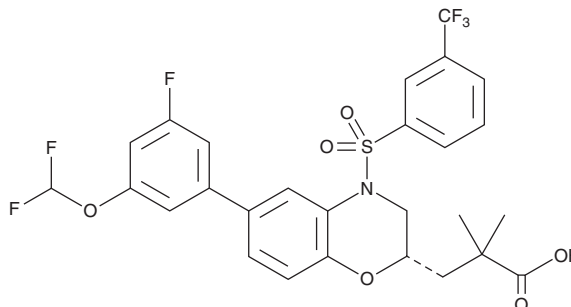


Cintirorgon

Item No. 28425

CAS Registry No.: 2055536-64-4
Formal Name: (2S)-6-[3-(difluoromethoxy)-5-fluorophenyl]-3,4-dihydro- α,α -dimethyl-4-[[3-(trifluoromethyl)phenyl]sulfonyl]-2H-1,4-benzoxazine-2-propanoic acid

Synonym: LYC-55716
MF: C₂₇H₂₃F₆NO₆S
FW: 603.5
Purity: \geq 98%
UV/Vis.: λ_{\max} : 214, 267 nm
Supplied as: A crystalline solid
Storage: -20°C
Stability: \geq 4 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

Laboratory Procedures

Cintirorgon is supplied as a crystalline solid. A stock solution may be made by dissolving the cintirorgon in the solvent of choice, which should be purged with an inert gas. Cintirorgon is soluble in the organic solvent DMSO at a concentration of approximately 100 mg/ml.

Description

Cintirorgon is a retinoic acid receptor-related orphan receptor γ (ROR γ) agonist with EC₅₀ values of less than 0.5 μ M in a reporter assay and time-resolved FRET (TR-FRET) assay.¹

Reference

1. Aicher, T.D., Taylor, C.B., Vanhuis, C.A. Aryl dihydro-2h-benzo[b][1,4]oxazine sulfonamide and related compounds for use as agonists of ROR γ and the treatment of disease. *Lycern Corporation*. WO2016/201225A1, (2016).

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, 12/05/2022

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