

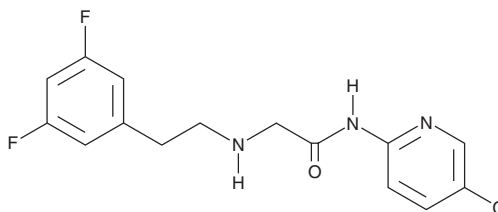
PRODUCT INFORMATION



SR 12460

Item No. 31066

CAS Registry No.: 2055101-66-9
Formal Name: N-(5-chloro-2-pyridinyl)-2-[[2-(3,5-difluorophenyl)ethyl]amino]-acetamide
MF: C₁₅H₁₄ClF₂N₃O
FW: 325.7
Purity: ≥98%
UV/Vis.: λ_{max}: 244, 288 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 12460 is supplied as a crystalline solid. A stock solution may be made by dissolving the SR 12460 in the solvent of choice, which should be purged with an inert gas. SR 12460 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of SR 12460 in these solvents is approximately 15 and 30 mg/ml, respectively.

Description

SR 12460 is a mimetic of the IκB kinase β (IKKβ) NEMO/IKKγ-binding domain (NBD) that inhibits the protein-protein interaction between the IKK complex subunits NF-κB essential modulator (NEMO/IKKγ) and IKKβ.¹ It inhibits TNF-α-induced NF-κB activation in a reporter assay when used at concentrations ranging from 25 to 150 μM. It reduces LPS-induced production of IL-6 in RAW 264.7 cells. SR 12460 (30 mg/kg) improves grip strength in the *mdx* mouse model of Duchenne muscular dystrophy.

Reference

1. Zhao, J., Zhang, L., Mu, X., *et al.* Development of novel NEMO-binding domain mimetics for inhibiting IKK/NF-κB activation. *PLoS Biol.* **16(6)**, e2004663 (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, 03/21/2024

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