

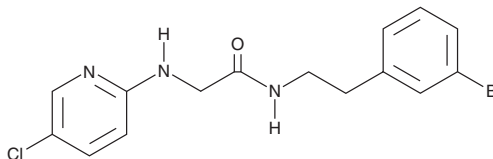
PRODUCT INFORMATION



SR 12343

Item No. 31031

CAS Registry No.: 2055101-86-3
Formal Name: N-[2-(3-bromophenyl)ethyl]-2-[(5-chloro-2-pyridinyl)amino]-acetamide
MF: C₁₅H₁₅BrClN₃O
FW: 368.7
Purity: ≥98%
UV/Vis.: λ_{max}: 249 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 12343 is supplied as a crystalline solid. A stock solution may be made by dissolving the SR 12343 in the solvent of choice, which should be purged with an inert gas. SR 12343 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of SR 12343 in these solvents is approximately 30 mg/ml.

Description

SR 12343 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 selectively disrupts the association of NEMO/IKKγ with IKKβ over TNF receptor-associated factor 2 (TRAF2), IκBα, and IKKα in co-immunoprecipitation assays using RAW 264.7 cells at concentrations ranging from 25 to 150 μM. SR 12343 inhibits TNF-α-induced NF-κB activation in a reporter assay. It reduces increased serum IL-6 levels in a mouse model of LPS-induced endotoxemia when administered at a dose of 10 mg/kg. SR 12343 (30 mg/kg) improves grip strength and reduces muscle fibrosis 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