

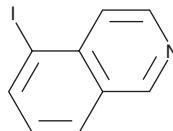
PRODUCT INFORMATION



DSRM-3716

Item No. 39517

CAS Registry No.: 58142-99-7
Formal Name: 5-iodo-isoquinoline
MF: C_9H_6IN
FW: 255.1
Purity: $\geq 95\%$
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

DSRM-3716 is supplied as a solid. A stock solution may be made by dissolving the DSRM-3716 in the solvent of choice, which should be purged with an inert gas. DSRM-3716 is soluble in methanol.

Description

DSRM-3716 is an inhibitor of the axonal degeneration mediator and NADase sterile alpha and TIR motif containing 1 (SARM1; $IC_{50} = 75 \text{ nM}$).¹ It decreases levels of cyclic ADP ribose (cADPR) and increases NAD^+ levels in axotomized primary mouse dorsal root ganglia (DRG) axons when used at concentrations ranging from 1.1 to 30 μM . DSRM-3716 (0.1-10 μM) prevents axonal degeneration in primary wild-type but not *Sarm1*^{-/-} axotomized DRG axons.

Reference

1. Hughes, R.O., Bosanac, T., Mao, X., *et al.* Small molecule SARM1 inhibitors recapitulate the SARM1^{-/-} phenotype and allow recovery of a metastable pool of axons fated to degenerate. *Cell Rep.* **34**(1), 108588 (2021).

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, 01/22/2024

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