

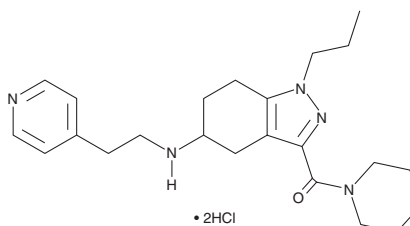
PRODUCT INFORMATION



NUCC-390 (hydrochloride)

Item No. 30957

CAS Registry No.: 2749281-71-6
Formal Name: 1-piperidinyl[4,5,6,7-tetrahydro-1-propyl-5-[[2-(4-pyridinyl)ethyl]amino]-1H-indazol-3-yl]-methanone, dihydrochloride
MF: $C_{23}H_{33}N_5O \cdot 2HCl$
FW: 468.5
Purity: $\geq 98\%$
Supplied as: A solid
Storage: $-20^{\circ}C$
Stability: ≥ 4 years



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

Laboratory Procedures

NUCC-390 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the NUCC-390 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. NUCC-390 (hydrochloride) is soluble in DMSO.

Description

NUCC-390 is a chemokine (C-X-C motif) receptor 4 (CXCR4) agonist.¹ It induces calcium mobilization in C8161 melanoma cells expressing CXCR4 when used at a concentration of 10 μM , an effect that can be prevented by the CXCR4 antagonist AMD3100 (plerixafor; Item No. 10011332). NUCC-390 (0.0625, 0.125, and 0.25 μM) increases axon length in cultured rat spinal cord motor neurons.² *In vivo*, NUCC-390 (26 mg/kg per day) restores gastrocnemius muscle compound muscle action potentials (CMAPs) in a mouse model of sciatic nerve crush injury. It also restores CMAPs in a mouse model of neuroparalysis induced by Papuan taipan (*O. scutellatus*) venom.³

References

1. Mishra, R.K., Shum, A.K., Platanias, L.C., *et al.* Discovery and characterization of novel small-molecule CXCR4 receptor agonists and antagonists. *Sci. Rep.* **6**:30155, 1-9 (2016).
2. Zanetti, G., Negro, S., Megighian, A., *et al.* A CXCR4 receptor agonist strongly stimulates axonal regeneration after damage. *Ann. Clin. Transl. Neurol.* **6**(12), 2395-2402 (2019).
3. Stazi, M., D'Este, G., Mattarei, A., *et al.* An agonist of the CXCR4 receptor accelerates the recovery from the peripheral neuroparalysis induced by Taipan snake envenomation. *PLoS Negl. Trop. Dis.* **14**(9), e0008547 (2020).

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, 07/24/2023

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