

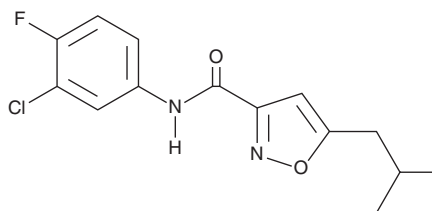
PRODUCT INFORMATION



SJ000063181

Item No. 42546

CAS Registry No.: 945189-68-4
Formal Name: N-(3-chloro-4-fluorophenyl)-5-(2-methylpropyl)-3-isoxazolecarboxamide
Synonyms: Ventromorphin 2, Vm2
MF: C₁₄H₁₄ClFN₂O₂
FW: 296.7
Purity: ≥98%
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

SJ000063181 is supplied as a solid. A stock solution may be made by dissolving the SJ000063181 in the solvent of choice, which should be purged with an inert gas. SJ000063181 is sparingly soluble (1-10 mg/ml) in ethanol and DMSO.

Description

SJ000063181 is a bone morphogenetic protein (BMP) signaling activator.¹ It activates BMP4 in C33A-2D2 human cervical carcinoma clonal reporter cells (EC₅₀ = ≤1 μM). SJ000063181 (6.25 μM) induces ventralization, as well as increases expression of the BMP target genes *bmp2b* and *szl* and the cooperatively induced BMP target genes *vent* and *vox* in zebrafish embryos. It also increases the stability of survival motor neuron protein (SMN; EC₅₀ = 1.5 μM) in a reporter assay using the *SMN2* gene, which produces an unstable truncated SMN protein.²

References

1. Genthe, J.R., Min, J., Farmer, D.M., *et al.* Ventromorphins: A new class of small molecule activators of the canonical BMP signaling pathway. *ACS Chem. Biol.* **12**(9), 2436-2447 (2017).
2. Rietz, A., Li, H., Quist, K.M., *et al.* Discovery of a small molecule probe that post-translationally stabilizes the survival motor neuron protein for the treatment of spinal muscular atrophy. *J. Med. Chem.* **60**(11), 4594-4610 (2017).

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, 02/27/2025

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