

PRODUCT INFORMATION

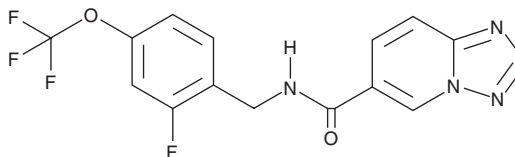


ONO-2920632

Item No. 43257

CAS Registry No.: 2230296-66-7
Formal Name: N-[[2-fluoro-4-(trifluoromethoxy)phenyl]methyl]-[1,2,4]triazolo[1,5-a]pyridine-6-carboxamide

Synonym: VU6011887
MF: C₁₅H₁₀F₄N₄O₂
FW: 354.3
Purity: ≥98%
Supplied as: A neat solid
Storage: 4°C
Stability: ≥4 years



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

Laboratory Procedures

ONO-2920632 is supplied as a neat solid. A stock solution may be made by dissolving the ONO-2920632 in the solvent of choice, which should be purged with an inert gas. ONO-2920632 is soluble (≥10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in ethanol.

Description

ONO-2920632 is an activator of two-pore domain potassium channel 2.1 (K_{2p}2.1/TREK-1) and K_{2p}10.1/TREK-2.¹ It induces K_{2p}2.1/TREK-1 and K_{2p}10.1/TREK-2 activity in thallium flux assays using CHO cells expressing K_{2p}2.1/TREK-1 and HEK293 cells expressing K_{2p}10.1/TREK-2 (EC₅₀s = 2.8 and 0.3 μM, respectively). ONO-2920632 is selective for K_{2p}2.1/TREK-1 and K_{2p}10.1/TREK-2 over 72 other ion channels, G protein-coupled receptors (GPCRs), transporters, and nuclear hormones at 10 μM. *In vivo*, ONO-2920632 (3 mg/kg) prevents acetic acid-induced writhing in mice.

Reference

1. Yashiro, K., Iwaki, Y., Urata, H., *et al.* Discovery of ONO-2920632 (VU6011887): A highly selective and CNS penetrant TREK-2 (TWIK-related K⁺ channel 2) preferring activator *in vivo* tool compound. *ACS Chem. Neurosci.* **16**(5), 960-967 (2025).

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, 05/07/2025

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