

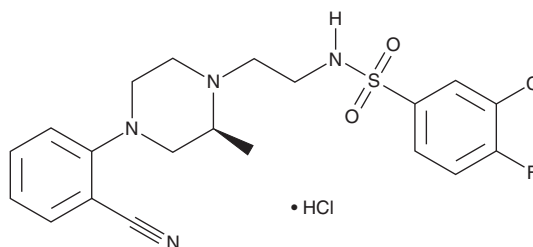
# PRODUCT INFORMATION



## VU6036720 (hydrochloride)

Item No. 40286

**Formal Name:** (S)-3-chloro-N-(2-(4-(2-cyanophenyl)-2-methylpiperazin-1-yl)ethyl)-4-fluorobenzenesulfonamide, monohydrochloride  
**MF:** C<sub>20</sub>H<sub>22</sub>ClFN<sub>4</sub>O<sub>2</sub>S • HCl  
**FW:** 473.4  
**Purity:** ≥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

VU6036720 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the VU6036720 (hydrochloride) in the solvent of choice which should be purged with an inert gas. VU6036720 (hydrochloride) is soluble in DMSO and is slightly soluble in acetonitrile.

VU6036720 (hydrochloride) is sparingly soluble in aqueous solutions. To enhance aqueous solubility, dilute the organic solvent solution into aqueous buffers or isotonic saline. If performing biological experiments, ensure the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. We do not recommend storing the aqueous solution for more than one day.

### Description

VU6036720 is an inhibitor of heteromeric inward-rectifier potassium channels (K<sub>ir</sub>) composed of K<sub>ir</sub>4.1 and K<sub>ir</sub>5.1 subunits (IC<sub>50</sub> = 0.24 μM).<sup>1</sup> It is selective for K<sub>ir</sub>4.1/5.1 channels over a variety of homomeric K<sub>ir</sub> channels, including K<sub>ir</sub>4.1 (IC<sub>50</sub> = >10 μM), as well as heteromeric K<sub>ir</sub>3.1/3.2, K<sub>ir</sub>3.1/3.4, K<sub>ir</sub>6.2/SUR1, and K<sub>ir</sub>6.1/SUR2b channels at 30 μM but does inhibit the voltage-gated potassium channel K<sub>v</sub>11.1, also known as human ether-a-go-go (hERG; IC<sub>50</sub> = 6.4 μM). VU6036720 (30 and 100 mg/kg) does not inhibit diuresis in mice due to low oral bioavailability and high hepatic clearance.

### Reference

1. McClenahan, S.J., Kent, C.N., Kharade, S.V., *et al.* VU6036720: The first potent and selective in vitro inhibitor of heteromeric Kir4.1/5.1 inward rectifier potassium channels. *Mol. Pharmacol.* **101(5)**, 357-370 (2022).

#### 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, 02/14/2024

#### 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