

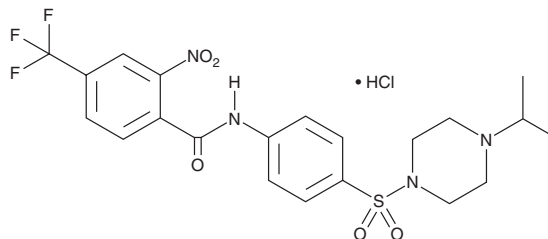
# PRODUCT INFORMATION



## RN-9893 (hydrochloride)

Item No. 29172

**CAS Registry No.:** 2109450-40-8  
**Formal Name:** N-[4-[[4-(1-methylethyl)-1-piperazinyl]sulfonyl]phenyl]-2-nitro-4-(trifluoromethyl)-benzamide, monohydrochloride  
**MF:** C<sub>21</sub>H<sub>23</sub>F<sub>3</sub>N<sub>4</sub>O<sub>5</sub>S • HCl  
**FW:** 537.0  
**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

RN-9893 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the RN-9893 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. RN-9893 (hydrochloride) is soluble in DMSO.

### Description

RN-9893 is an antagonist of transient receptor potential vanilloid 4 (TRPV4; IC<sub>50</sub>s = 0.42 and 0.66 μM, respectively, for the human and rat receptors).<sup>1</sup> It is selective for TRPV4 over TRPV1, TRPV3, and TRPM8 (IC<sub>50</sub>s = 10, >30, and 30 μM, respectively). RN-9893 reduces rat TRPV4 activity induced by 4α-phorbol 12,13-didecanoate (4α-PDD; Item No. 20446) or hypotonicity (IC<sub>50</sub>s = 0.57 and 2.1 μM, respectively, in cell free assays).

### Reference

1. Nguyen, M.T., O'Mahony, D.J., *et al.* Identification of orally-bioavailable antagonists of the TRPV4 ion-channel. *Bioorg. Med. Chem. Lett.* **25(18)**, 4011-4015 (2015).

#### 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, 12/07/2022

#### 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