

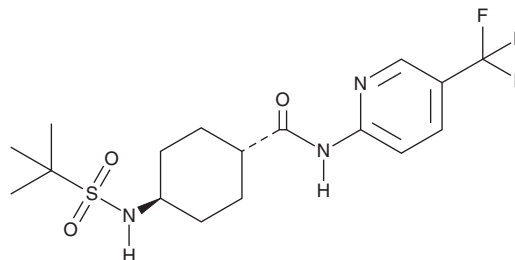
# PRODUCT INFORMATION



## Velneperit

Item No. 33850

CAS Registry No.: 342577-38-2  
Formal Name: *trans*-4-[[[1,1-dimethylethyl)sulfonyl]amino]-N-[5-(trifluoromethyl)-2-pyridinyl]-cyclohexanecarboxamide  
Synonym: S-2367  
MF: C<sub>17</sub>H<sub>24</sub>F<sub>3</sub>N<sub>3</sub>O<sub>3</sub>S  
FW: 407.5  
Purity: ≥98%  
UV/Vis.: λ<sub>max</sub>: 244, 279 nm  
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

Velneperit is supplied as a solid. A stock solution may be made by dissolving the velneperit in the solvent of choice, which should be purged with an inert gas. Velneperit is soluble in the organic solvent DMSO.

### Description

Velneperit is a nonpeptide antagonist of neuropeptide Y (NPY) receptor Y<sub>5</sub> (K<sub>i</sub> = 5.3 nM).<sup>1</sup> It is selective for Y<sub>5</sub> over Y<sub>1</sub>, Y<sub>2</sub>, and Y<sub>4</sub> receptors, as well as a panel of 44 receptors, ion channels, transporters and enzymes, at 10 μM. Velneperit inhibits NPY-induced decreases in forskolin-induced cAMP production in CHO-K1 cells expressing the human Y<sub>5</sub> receptor (IC<sub>50</sub> = 169 nM). *In vivo*, velneperit (50 mg/kg) suppresses Y<sub>5</sub> agonist-induced increases in food intake in high-fat diet-induced obese (DIO) mice. It also reduces body weight gain and suppresses food intake in DIO mice.

### Reference

1. Fukasaka, Y., Nambu, H., Tanioka, H., *et al.* An insurmountable NPY Y5 receptor antagonist exhibits superior anti-obesity effects in high-fat diet-induced obese mice. *Neuropeptides* **70**, 55-63 (2018).

#### 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, 11/17/2022

#### 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