

# PRODUCT INFORMATION



## VU0453379 (hydrochloride)

Item No. 41392

**Formal Name:** 2-cyclopentyl-2,9-dihydro-9-methyl-N-[[[(2S)-1-(1-methylethyl)-2-pyrrolidinyl]methyl]-1-oxo-1H-pyrido[3,4-b]indole-4-carboxamide, monohydrochloride

**MF:** C<sub>26</sub>H<sub>34</sub>N<sub>4</sub>O<sub>2</sub> • HCl

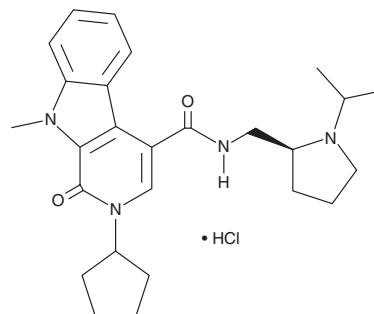
**FW:** 471.0

**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

VU0453379 (hydrochloride) is supplied as a solid. A stock solution may be made by dissolving the VU0453379 (hydrochloride) in the solvent of choice, which should be purged with an inert gas. VU0453379 (hydrochloride) is sparingly soluble (1-10 mg/ml) in DMSO and slightly soluble (0.1-1 mg/ml) in acetonitrile.

### Description

VU0453379 is an allosteric agonist and positive allosteric modulator (ago-PAM) of glucagon-like peptide 1 receptor (GLP-1R).<sup>1</sup> It induces calcium mobilization in 9-3-H cells expressing human GLP-1R (EC<sub>50</sub> = 1.3 μM) and enhances GLP-1-, exenatide-, or liraglutide-induced calcium mobilization in the same cells (EC<sub>50</sub>s = 1.8, 8.4, and 30 μM, respectively). VU0453379 (30 μM) potentiates exenatide-induced insulin secretion in primary mouse pancreatic islets. It reverses catalepsy induced by haloperidol (Item No. 12014) in a rat model of Parkinson's disease motor symptoms when administered at a dose of 30 mg/kg.

### Reference

1. Morris, L.C., Nance, K.D., Gentry, P.R., *et al.* Discovery of (S)-2-cyclopentyl-N-((1-isopropylpyrrolidin-2-yl)-9-methyl-1-oxo-2,9-dihydro-1H-pyrido[3,4-b]indole-4-carboxamide (VU0453379): A novel, CNS penetrant glucagon-like peptide 1 receptor (GLP-1R) positive allosteric modulator (PAM). *J. Med. Chem.* **57**(23), 10192-10197 (2014).

#### 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, 09/27/2024

#### 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