

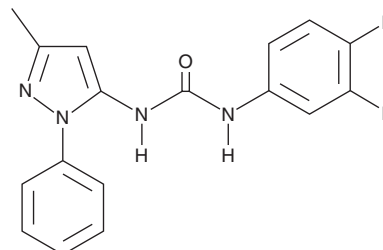
# PRODUCT INFORMATION



**ML-297**

Item No. 14792

**CAS Registry No.:** 1443246-62-5  
**Formal Name:** N-(3,4-difluorophenyl)-N'-(3-methyl-1-phenyl-1H-pyrazol-5-yl)-urea  
**Synonyms:** CID-56642816, VU0456810  
**MF:** C<sub>17</sub>H<sub>14</sub>F<sub>2</sub>N<sub>4</sub>O  
**FW:** 328.3  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 244 nm  
**Supplied as:** A crystalline 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

ML-297 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-297 in the solvent of choice, which should be purged with an inert gas. ML-297 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of ML-297 in ethanol is approximately 20 mg/ml and approximately 30 mg/ml in DMSO and DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of ML-297 can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of ML-297 in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

## Description

G protein-regulated inwardly rectifying potassium (GIRK1-4) channels are a family of K<sub>ir</sub>3.1-K<sub>ir</sub>3.4 ion channels that modulate cell excitability. The four different GIRK subunits are composed in different homo- and heterotetrameric combinations, which are expressed with regional specificity throughout the central nervous system and in the periphery.<sup>1</sup> ML-297 is a selective GIRK1/2 activator (EC<sub>50</sub>s = 0.16 and 1.8 μM for GIRK1/2 and GIRK1/4, respectively, and is completely inactive at GIRK2/3).<sup>1,2</sup> In two different mouse models of epilepsy, ML-297 at 60 mg/kg was shown to delay seizure onset and to prevent convulsions.<sup>1</sup>

## References

1. Kaufmann, K., Romaine, I.M., Days, E., *et al.* ML297 (VU0456810), the first potent and selective activator of the GIRK potassium channel, displays antiepileptic properties in mice. *ACS Chem. Neurosci.* **4**(9), 1278-1286 (2013).
2. Wen, W., Wu, W., Romaine, I.M., *et al.* Discovery of 'molecular switches' within a GIRK activator scaffold that afford selective GIRK inhibitors. *Bioorg. Med. Chem. Lett.* **23**(16), 4562-4566 (2013).

### 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, 01/31/2024

## 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](http://WWW.CAYMANCHEM.COM)