

# PRODUCT INFORMATION



**HC-067047**

Item No. 20927

**CAS Registry No.:** 883031-03-6  
**Formal Name:** 2-methyl-1-[3-(4-morpholinyl)propyl]-5-phenyl-N-[3-(trifluoromethyl)phenyl]-1H-pyrrole-3-carboxamide

**MF:** C<sub>26</sub>H<sub>28</sub>F<sub>3</sub>N<sub>3</sub>O<sub>2</sub>

**FW:** 471.5

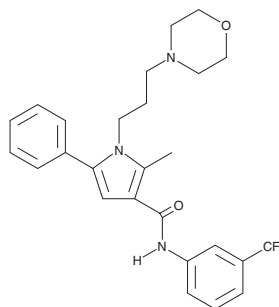
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 279 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

HC-067047 is supplied as a crystalline solid. A stock solution may be made by dissolving the HC-067047 in the solvent of choice. HC-067047 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF), which should be purged with an inert gas. The solubility of HC-067047 in these solvents is approximately 1, 15, and 20 mg/ml, respectively.

HC-067047 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, HC-067047 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. HC-067047 has a solubility of approximately 0.33 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

HC-067047 is a TRPV4 antagonist that reversibly inhibits currents through human, rat, and mouse TRPV4 orthologs (IC<sub>50</sub>s = 48, 133, and 17 nM, respectively).<sup>1</sup> It is selective for TRPV4 over TRPV1, TRPV2, TRPV3, and TRPM8 channels.<sup>1</sup> HC-067047 has been shown to increase functional bladder capacity and to reduce micturition frequency in wild-type mice and rats with cystitis.<sup>1</sup>

## Reference

1. Everaerts, W., Zhen, X., Ghosh, D., *et al.* Inhibition of the cation channel TRPV4 improves bladder function in mice and rats with cyclophosphamide-induced cystitis. *Proc. Natl. Acad. Sci. USA* **107**(44), 19084-19089 (2010).

### 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, 02/15/2024

## 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)