

# PRODUCT INFORMATION



## Lasmiditan

Item No. 28369

**CAS Registry No.:** 439239-90-4  
**Formal Name:** 2,4,6-trifluoro-N-[6-[(1-methyl-4-piperidiny)carbonyl]-2-pyridinyl]-benzamide

**Synonyms:** COL-144, LY573144

**MF:** C<sub>19</sub>H<sub>18</sub>F<sub>3</sub>N<sub>3</sub>O<sub>2</sub>

**FW:** 377.4

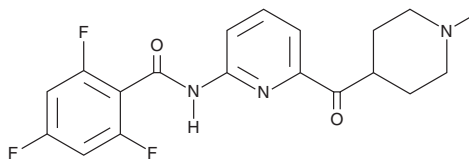
**Purity:** ≥98%

**UV/Vis.:** λ<sub>max</sub>: 225, 256, 296 nm

**Supplied as:** A crystalline solid

**Storage:** -20°C

**Stability:** ≥5 years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Description

Lasmiditan (Item No. 28369) is an analytical reference standard categorized as an agonist of the serotonin (5-HT) receptor subtype 5-HT<sub>1F</sub>.<sup>1</sup> Formulations containing lasmiditan have been used in the treatment of migraine headache. Lasmiditan has low abuse potential.<sup>2</sup> Lasmiditan is regulated as a Schedule V compound in the United States. This product is intended for research and forensic applications.

### References

1. Nelson, D.L., Phebus, L.A., Johnson, K.W., *et al.* Preclinical pharmacological profile of the selective 5-HT<sub>1F</sub> receptor agonist lasmiditan. *Cephalalgia* **30(10)**, 1159-1169 (2010).
2. Wilbraham, D., Berg, P.H., Tsai, M., *et al.* Abuse potential of lasmiditan: A phase 1 randomized, placebo- and alprazolam-controlled crossover study. *J. Clin. Pharmacol.* (2019).

#### 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, 07/24/2023

#### 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