

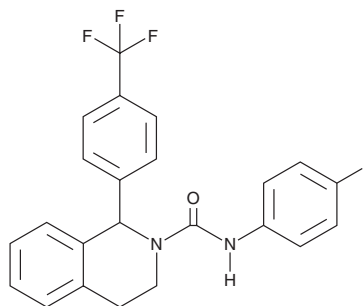
# PRODUCT INFORMATION



## TC-I 2000

Item No. 35463

**CAS Registry No.:** 1159996-20-9  
**Formal Name:** N-(4-fluorophenyl)-3,4-dihydro-1-[4-(trifluoromethyl)phenyl]-2(1H)-isoquinolinecarboxamide  
**MF:** C<sub>23</sub>H<sub>18</sub>F<sub>4</sub>N<sub>2</sub>O  
**FW:** 414.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 234 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

TC-I 2000 is supplied as a solid. A stock solution may be made by dissolving the TC-I 2000 in the solvent of choice, which should be purged with an inert gas. TC-I 2000 is soluble in the organic solvent ethanol at a concentration of approximately 1 mg/ml.

### Description

TC-I 2000 is an antagonist of transient receptor potential melastatin 8 (TRPM8; IC<sub>50</sub> = 53 nM for the rat receptor).<sup>1</sup>

### Reference

1. Tamayo, N.A., Bo, Y., Gore, V., *et al.* Fused piperidines as a novel class of potent and orally available transient receptor potential melastatin type 8 (TRPM8) antagonists. *J. Med. Chem.* **55**(4), 1593-1611 (2012).

#### 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, 12/12/2022

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