

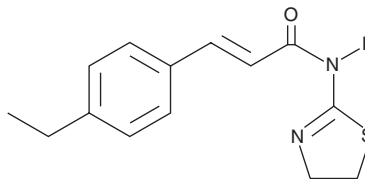
# PRODUCT INFORMATION



## ML-243

Item No. 15189

**CAS Registry No.:** 1426576-80-8  
**Formal Name:** (2E)-N-(4,5-dihydro-2-thiazolyl)-3-(4-ethylphenyl)-2-propenamide  
**Synonym:** CID-50910523  
**MF:** C<sub>14</sub>H<sub>16</sub>N<sub>2</sub>OS  
**FW:** 260.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 225, 308 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-243 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-243 in the solvent of choice, which should be purged with an inert gas. ML-243 is soluble in DMSO at a concentration of approximately 5 mg/ml.

### Description

ML-243 is an inhibitor of cancer stem cells, displaying 32-fold selective inhibition of a breast cancer stem cell-like cell line (EC<sub>50</sub> = 2.0 μM) over a control (mammary epithelial) cell line (EC<sub>50</sub> = 64 μM).<sup>1</sup> It weakly antagonizes the adenosine A<sub>2A</sub> receptor (IC<sub>50</sub> = 10 μM) and has no effect on 68 other targets commonly used in drug discovery for lead profiling.<sup>1</sup>

### Reference

1. Carmody, L., Germain, A., Morgan, B., *et al.* Identification of a selective small-molecule inhibitor of breast cancer stem cells - Probe 3, in Probe Reports from the NIH Molecular Libraries Program, 3, 1 (2011).

#### 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, 11/16/2022

#### 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