

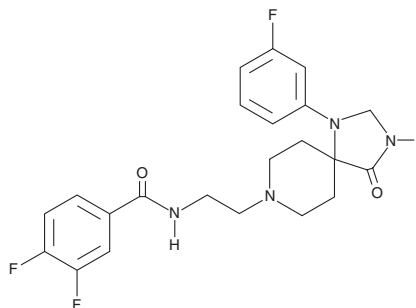
# PRODUCT INFORMATION



## ML-298

Item No. 15349

**CAS Registry No.:** 1426916-02-0  
**Formal Name:** 3,4-difluoro-N-[2-[1-(3-fluorophenyl)-4-oxo-1,3,8-triazaspiro[4.5]dec-8-yl]ethyl]-benzamide  
**MF:** C<sub>22</sub>H<sub>23</sub>F<sub>3</sub>N<sub>4</sub>O<sub>2</sub>  
**FW:** 432.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 203, 249 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-298 is supplied as a crystalline solid. A stock solution may be made by dissolving the ML-298 in the solvent of choice. ML-298 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide, which should be purged with an inert gas. The solubility of ML-298 in these solvents is approximately 1, 20, and 10 mg/ml, respectively.

### Description

ML-298 is an inhibitor of phospholipase D<sub>2</sub> (PLD<sub>2</sub>; IC<sub>50</sub> = 355 nM).<sup>1</sup> It is selective for PLD<sub>2</sub> over PLD<sub>1</sub> (IC<sub>50</sub> = 20,000 nM). It decreases invasive migration of U87-MG glioblastoma cells when used at a concentration of 10 μM without inducing cytotoxicity.

### Reference

- O'Reilly, M.C., Scott, S.A., Brown, K.A., *et al.* Development of dual PLD1/2 and PLD2 selective inhibitors from a common 1,3,8-triazaspiro[4.5]decane core: Discovery of ML298 and ML299 that decrease invasive migration in U87-MG glioblastoma cells. *J. Med Chem.* **56(6)**, 2695-2699 (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, 12/04/2017

#### 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