

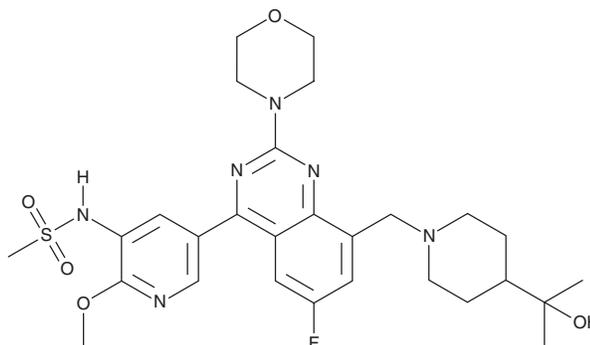
# PRODUCT INFORMATION



## Linperlisib

Item No. 44441

**CAS Registry No.:** 1702816-75-8  
**Formal Name:** N-[5-[6-fluoro-8-[[4-(1-hydroxy-1-methyl ethyl)-1-piperidinyl]methyl]-2-(4-morpholinyl)-4-quinazoliny]-2-methoxy-3-pyridinyl]-methanesulfonamide  
**Synonym:** YY-20394  
**MF:** C<sub>28</sub>H<sub>37</sub>FN<sub>6</sub>O<sub>5</sub>S  
**FW:** 588.7  
**Purity:** ≥95%  
**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

Linperlisib is supplied as a solid. A stock solution may be made by dissolving the linperlisib in the solvent of choice, which should be purged with an inert gas. Linperlisib is slightly soluble (0.1-1 mg/ml) in DMSO.

### Description

Linperlisib is a PI3K $\delta$  inhibitor (IC<sub>50</sub> = 4.6 nM).<sup>1</sup> It is selective for PI3K $\delta$  over PI3K $\alpha$ , PI3K $\beta$ , and PI3K $\gamma$  (IC<sub>50</sub>s = 1,159, 138, and 5,175 nM, respectively).

### Reference

1. Jiang, B., Qi, J., Song, Y., *et al.* Phase 1 clinical trial of the PI3K $\delta$  inhibitor YY-20394 in patients with B-cell hematological malignancies. *J. Hematol. Oncol.* **14**(1), 130 (2021).

#### 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, 10/21/2025

#### 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