

# PRODUCT INFORMATION

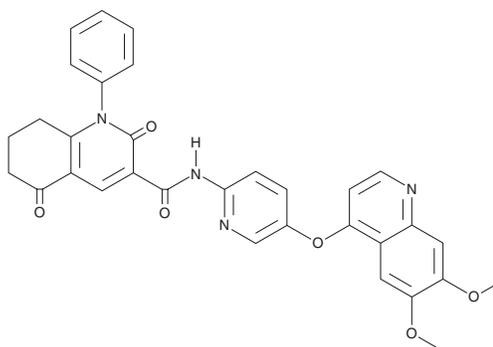


**ONO-7475**

Item No. 33258

**CAS Registry No.:** 1646839-59-9  
**Formal Name:** N-[5-[(6,7-dimethoxy-4-quinolinyl)oxy]-2-pyridinyl]-1,2,5,6,7,8-hexahydro-2,5-dioxo-1-phenyl-3-quinolinecarboxamide

**MF:** C<sub>32</sub>H<sub>26</sub>N<sub>4</sub>O<sub>6</sub>  
**FW:** 562.6  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 240, 279, 325 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

ONO-7475 is supplied as a crystalline solid. A stock solution may be made by dissolving the ONO-7475 in the solvent of choice, which should be purged with an inert gas. ONO-7475 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of ONO-7475 in these solvents is approximately 30 mg/ml.

ONO-7475 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, ONO-7475 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. ONO-7475 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

## Description

ONO-7475 is an inhibitor of the TAM family kinases Axl and Mer (IC<sub>50</sub>s = 0.7 and 1 nM, respectively).<sup>1</sup> It is selective for Axl and Mer over FLT3 (IC<sub>50</sub> = 147 nM). ONO-7475 (10 nM) reduces viability of, and induces apoptosis in, MOLM-13 and MV4-11 leukemia cells. *In vivo*, ONO-7475 (6 mg/kg) reduces leukemic burden and increases the median survival time in a MOLM-13 mouse xenograft model.

## Reference

1. Ruvolo, P.P., Ma, H., Ruvolo, V.R., *et al.* Anexelekt/MER tyrosine kinase inhibitor ONO-7475 arrests growth and kills FMS-like tyrosine kinase 3-internal tandem duplication mutant acute myeloid leukemia cells by diverse mechanisms. *Haematologica* **102(12)**, 2048-2057 (2017).

### 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, 10/05/2022

## 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