# **PRODUCT** INFORMATION



## MOMIPP

Item No. 39671

CAS Registry No.: Formal Name:	1363421-46-8 (2E)-3-(5-methoxy-2-methyl- 1H-indol-3-yl)-1-(4-pyridinyl)-2- propen-1-one	N O
MF:	$C_{18}H_{16}N_2O_2$	
FW:	292.3	
Purity:	≥98%	/
Supplied as:	A solid	N
Storage:	-20°C	/ ~
Stability:	≥4 years	Н

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

#### Laboratory Procedures

MOMIPP is supplied as a solid. A stock solution may be made by dissolving the MOMIPP in the solvent of choice, which should be purged with an inert gas. MOMIPP is soluble in organic solvents such as DMSO. MOMIPP is slightly soluble in acetonitrile and methanol.

### Description

 $\label{eq:MOMIPP} \begin{array}{l} \text{MOMIPP} \mbox{ is an inhibitor of PIKfyve (IC_{50} = 5.05 \mbox{ nM}).^1 It selectively binds to PIKfyve over phosphatidylinositol 4-phosphate 5-kinase type-1 <math display="inline">\gamma$  (PIP5K1C; K\_ds = 5.3 and >15,000 \mbox{ nM}, respectively). MOMIPP (10  $\mu\text{M})$ induces vacuolization and cell lysis, markers of methuosis, and decreases the extracellular acidification rate (ECAR), indicating inhibition of glycolytic flux, but not the oxygen consumption rate (OCR), in U251 glioblastoma cells.<sup>2</sup> Intraperitoneal administration of MOMIPP (80 mg/kg) decreases tumor growth in a U251 orthotopic mouse xenograft model.

#### References

- 1. Cho, H., Geno, E., Patoor, M., et al. Indolyl-pyridinyl-propenone-induced methuosis through the inhibition of PIKFYVE. ACS Omega 3(6), 6097-6103 (2018).
- 2. Li, Z., Mbah, N.E., Overmeyer, J.H., et al. The JNK signaling pathway plays a key role in methuosis (non-apoptotic cell death) induced by MOMIPP in glioblastoma. BMC Cancer 19(1), 77 (2019).

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.

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, 11/06/2023

## 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