

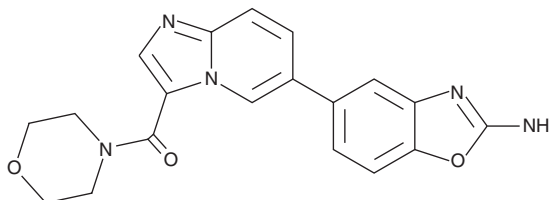
PRODUCT INFORMATION



INK1117

Item No. 19514

CAS Registry No.: 1268454-23-4
Formal Name: [6-(2-amino-5-benzoxazolyl)imidazo[1,2-a]pyridin-3-yl]-4-morpholinyl-methanone
Synonyms: MLN1117, TAK-117
MF: C₁₉H₁₇N₅O₃
FW: 363.4
Purity: ≥98%
UV/Vis.: λ_{max}: 234, 298 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

INK1117 is supplied as a crystalline solid. A stock solution may be made by dissolving the INK1117 in the solvent of choice, which should be purged with an inert gas. INK1117 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of INK1117 in these solvents is approximately 1 and 0.25 mg/ml, respectively.

INK1117 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, INK1117 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. INK1117 has a solubility of approximately 0.09 mg/ml in a 1:10 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

INK1117 is an inhibitor of phosphoinositide 3-kinase α (PI3K α) that is selective for p110 α *in vitro* (IC₅₀ = 15 nM for PI3K α vs. >1 μ M for other isoforms, as well as for mTOR) and in cells when used at 1 μ M.^{1,2} It blocks signaling to Akt and inhibits growth of cancer cells harboring wild-type or mutated p110 α .¹ INK1117 does not interfere with B cell proliferation or NK cell maturation and survival.^{1,2}

References

1. So, L., Yea, S.S., Oak, J.S., *et al.* Selective inhibition of phosphoinositide 3-kinase p110 α preserves lymphocyte function. *J. Biol. Chem.* **288**(8), 5718-5731 (2013).
2. Yea, S.S., So, L., Mallya, S., *et al.* Effects of novel isoform-selective phosphoinositide 3-kinase inhibitors on natural killer cell function. *PLoS One* **9**(6), (2014).

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, 12/02/2022

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