

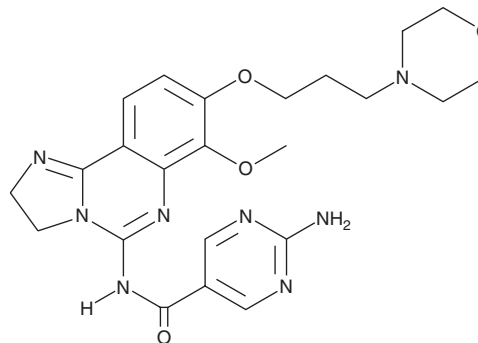
# PRODUCT INFORMATION



## Copanlisib

Item No. 20354

**CAS Registry No.:** 1032568-63-0  
**Formal Name:** 2-amino-N-[2,3-dihydro-7-methoxy-8-[3-(4-morpholinyl)propoxy]imidazo[1,2-c]quinazolin-5-yl]-5-pyrimidinecarboxamide  
**Synonym:** BAY 80-6946  
**MF:** C<sub>23</sub>H<sub>28</sub>N<sub>8</sub>O<sub>4</sub>  
**FW:** 480.5  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 222, 262, 275, 310 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

Copanlisib is supplied as a crystalline solid. A stock solution may be made by dissolving the copanlisib in the solvent of choice. Copanlisib is sparingly soluble in organic solvents such as acetonitrile and warmed DMSO.

### Description

Copanlisib is an inhibitor of class I phosphoinositide 3-kinase (PI3K) isoforms (IC<sub>50</sub>s = 0.5, 3.7, 6.4, and 0.7 nM for α, β, γ, and δ, respectively).<sup>1</sup> It is weakly active or inactive against numerous other kinases. Copanlisib is effective *in vivo* when administered intravenously, inhibiting PI3K signaling and inducing tumor regression *via* apoptosis in xenograft tumors in mice.<sup>1,2</sup>

### References

1. Liu, N., Rowley, B.R., Bull, C.O., *et al.* BAY 80-6946 is a highly selective intravenous PI3K inhibitor with potent p110α and p110δ activities in tumor cell lines and xenograft models. *Mol Cancer Ther.* **12(11)**, 2319-2330 (2013)
2. Will, M., Qin, A.C.R., Toy, W., *et al.* Rapid induction of apoptosis by PI3K inhibitors is dependent upon their transient inhibition of RAS -ERK signaling. *Cancer Discov.* **4(3)**, 334-347 (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, 12/19/2022

#### 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