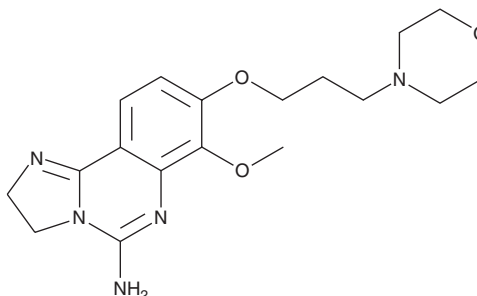


PRODUCT INFORMATION

7-Methoxy-8-(3-morpholin-4-ylpropoxy)-2,3-dihydroimidazo[1,2-c]quinazolin-5-amine

Item No. 30944

CAS Registry No.: 1032570-74-3
Formal Name: 2,3-dihydro-7-methoxy-8-[3-(4-morpholinyl)propoxy]-imidazo[1,2-c]quinazolin-5-amine
MF: C₁₈H₂₅N₅O₃
FW: 359.4
Purity: ≥95%
UV/Vis.: λ_{max}: 214, 264 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

7-Methoxy-8-(3-morpholin-4-ylpropoxy)-2,3-dihydroimidazo[1,2-c]quinazolin-5-amine is supplied as a crystalline solid. A stock solution may be made by dissolving the 7-methoxy-8-(3-morpholin-4-ylpropoxy)-2,3-dihydroimidazo[1,2-c]quinazolin-5-amine in the solvent of choice, which should be purged with an inert gas. 7-Methoxy-8-(3-morpholin-4-ylpropoxy)-2,3-dihydroimidazo[1,2-c]quinazolin-5-amine is soluble in the organic solvent chloroform at a concentration of approximately 10 mg/ml.

Description

7-Methoxy-8-(3-morpholin-4-ylpropoxy)-2,3-dihydroimidazo[1,2-c]quinazolin-5-amine is a building block.¹ It has been used in the synthesis of PI3K inhibitors.

Reference

1. Scott, W.J., Hentemann, M.F., Rowley, R.B., *et al.* Discovery and SAR of novel 2,3-dihydroimidazo[1,2-c]quinazoline PI3K inhibitors: Identification of copanlisib (BAY 80-6946). *ChemMedChem.* **11**(14), 1517-1530 (2016).

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, 12/09/2022

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