

PRODUCT INFORMATION

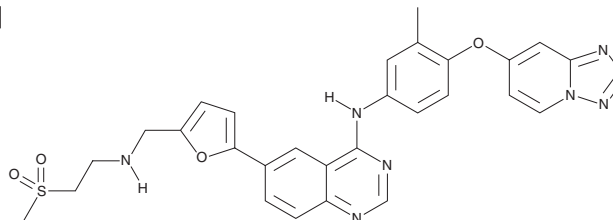


Arry-380 analog

Item No. 24186

CAS Registry No.: 937265-83-3
Formal Name: 6-[5-[[[2-(methylsulfonyl)ethyl]amino]methyl]-2-furanyl]-N-[3-methyl-4-([1,2,4]triazolo[1,5-a]pyridin-7-yloxy)phenyl]-4-quinazolinamine

MF: C₂₉H₂₇N₇O₄S
FW: 569.6
Purity: ≥98%
UV/Vis.: λ_{max}: 334, 361 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

Arry-380 analog is supplied as a crystalline solid. A stock solution may be made by dissolving the arry-380 analog in the solvent of choice. Arry-380 analog is soluble in organic solvents such as DMSO and dimethyl formamide, which should be purged with an inert gas. The solubility of arry-380 analog in these solvents is approximately 25 mg/ml.

Arry-380 analog is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, arry-380 analog should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. Arry-380 analog has a solubility of approximately 0.1 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

Arry-380 analog was designed as an inhibitor of EGFR (ErbB1).¹

Reference

1. Lyssikatos, J.P., Hicks, J.M., Marmsater, F.P., *et al.* N4-phenyl-quinazoline-4-amine derivatives and related compounds as ErbB type I receptor tyrosine kinase inhibitors for the treatment of hyperproliferative diseases. *Array Biopharma*. **US20170252317A1** (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, 10/21/2022

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