

# PRODUCT INFORMATION

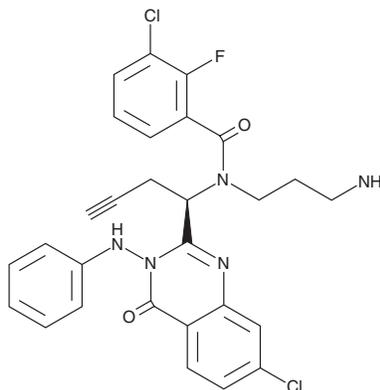


## ARQ 621

Item No. 34329

**CAS Registry No.:** 1095253-39-6  
**Formal Name:** N-(3-aminopropyl)-3-chloro-N-[(1R)-1-[7-chloro-3,4-dihydro-4-oxo-3-(phenylamino)-2-quinazoliny]-3-butyn-1-yl]-2-fluoro-benzamide

**MF:** C<sub>28</sub>H<sub>24</sub>Cl<sub>2</sub>FN<sub>5</sub>O<sub>2</sub>  
**FW:** 552.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 239 nm  
**Supplied as:** A 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

ARQ 621 is supplied as a solid. A stock solution may be made by dissolving the ARQ 621 in the solvent of choice, which should be purged with an inert gas. ARQ 621 is soluble in DMSO.

### Description

ARQ 621 is an Eg5 inhibitor.<sup>1</sup> It reduces tumor growth in a MIA PaCa-2 pancreatic cancer mouse xenograft model when administered at a dose of 6.25 mg/kg

### Reference

1. Jeay, S., Ali, S., Chen, C.-R., *et al.* Discovery of a novel Eg5 kinesin inhibitor, ARQ 621, with potent antitumor activity while sparing bone marrow-derived cells. *Cancer Res.* **68(9 Suppl.)**, 656 (2008).

#### 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, 12/08/2022

#### 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