

# PRODUCT INFORMATION

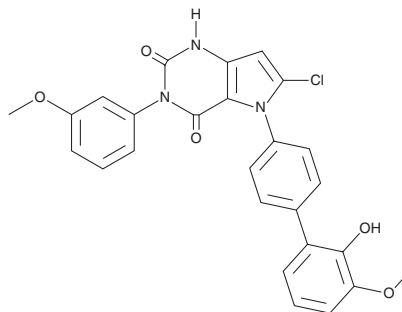


## GSK621

Item No. 19155

CAS Registry No.: 1346607-05-3  
Formal Name: 6-chloro-5-(2'-hydroxy-3'-methoxy[1,1'-biphenyl]-4-yl)-3-(3-methoxyphenyl)-1H-pyrrolo[3,2-d]pyrimidine-2,4(3H,5H)-dione

MF:  $C_{26}H_{20}ClN_3O_5$   
FW: 489.9  
Purity:  $\geq 98\%$   
UV/Vis.:  $\lambda_{max}$ : 273 nm  
Supplied as: A crystalline solid  
Storage:  $-20^\circ\text{C}$   
Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

GSK621 is supplied as a crystalline solid. A stock solution may be made by dissolving the GSK621 in the solvent of choice, which should be purged with an inert gas. GSK621 is soluble in organic solvents such as DMSO and dimethyl formamide. The solubility of GSK621 in these solvents is approximately 30 mg/ml.

GSK621 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, GSK621 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. GSK621 has a solubility of approximately 0.25 mg/ml in a 1:3 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

GSK621 is an activator of AMP-activated protein kinase (AMPK). It has been shown to increase phosphorylation at AMPK $\alpha$  T172, a marker of AMPK activation, in both acute myeloid leukemia (AML) cell lines and primary AML samples.<sup>1</sup> GSK621 is reported to reduce the proliferation of AML cell lines ( $IC_{50}$ s = 13-30  $\mu\text{M}$ ), triggering autophagy via the eIF2 $\alpha$ /ATF4 signaling pathway.<sup>1</sup>

### Reference

1. Sujobert, P., Poulain, L., Paubelle, E., *et al.* Co-activation of AMPK and mTORC1 induces cytotoxicity in acute myeloid leukemia. *Cell Rep.* **11**(9), 1446-1457 (2015).

#### 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, 11/29/2022

#### 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