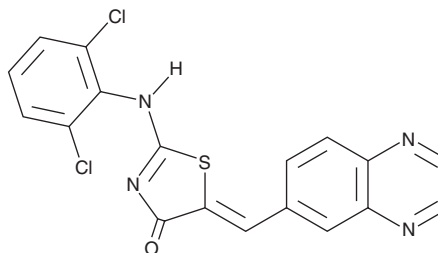


PRODUCT INFORMATION



GSK626616
Item No. 34621

CAS Registry No.: 1025821-33-3
Formal Name: (5Z)-2-[(2,6-dichlorophenyl)amino]-5-(6-quinoxalinylmethylene)-4(5H)-thiazolone
MF: C₁₈H₁₀Cl₂N₄OS
FW: 401.3
Purity: ≥98%
UV/Vis.: λ_{max}: 359 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

GSK626616 is supplied as a solid. A stock solution may be made by dissolving the GSK626616 in the solvent of choice, which should be purged with an inert gas. GSK626616 is soluble in ethanol.

Description

GSK626616 is an inhibitor of dual-specificity tyrosine phosphorylation-regulated kinase 3 (DYRK3; IC₅₀ = 0.7 nM).¹ It enhances erythropoietin-induced erythropoiesis in isolated human bone marrow. GSK626616 increases hemoglobin levels in a mouse model of carboplatin- and radiation-induced anemia.

Reference

1. Erickson-Miller, C.L., Creasy, C., Chadderton, A., *et al.* GSK626616: A DYRK3 inhibitor as a potential new therapy for the treatment of anemia. *Blood* **110**(11), 510 (2007).

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, 11/07/2022

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