

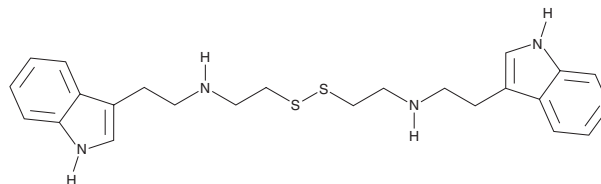
PRODUCT INFORMATION



G6PD Activator AG1

Item No. 39675

CAS Registry No.: 421581-52-4
Formal Name: N,N'-(dithiodi-2,1-ethanediyl)bis-1H-indole-3-ethanamine
Synonym: Glucose-6-phosphate Dehydrogenase Activator AG1
MF: C₂₄H₃₀N₄S₂
FW: 438.6
Purity: ≥95%
Supplied as: A neat oil
Storage: -20°C
Stability: ≥4 years
Special Conditions: Unstable in solutions



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

Laboratory Procedures

G6PD Activator AG1 is supplied as a neat oil. A stock solution may be made by dissolving the G6PD Activator AG1 in the solvent of choice, which should be purged with an inert gas. G6PD Activator AG1 is soluble in acetonitrile.

Description

G6PD activator AG1 is an activator of glucose-6-phosphate dehydrogenase (G6PDH; EC₅₀ = 3.4 μM for G6PDH^{R459L}).¹ It selectively induces the dimerization of G6PDH^{R459L} over 6-phosphogluconate dehydrogenase (6PGD), GAPDH, aldehyde dehydrogenase 2 (ALDH2), and ALDH3A1. G6PD activator AG1 decreases the levels of reactive oxygen species (ROS) in, and increases the viability of, primary human lymphocytes isolated from a patient carrying the G6PDH^{R459L} mutation. It inhibits hemolysis induced by chloroquine (Item No. 14194) or diamide in isolated human erythrocytes when used at a concentration of 5 μM. G6PD activator AG1 (1 μM) decreases chloroquine-induced pericardial edema in zebrafish embryos. It decreases tumor volume and weight and increases survival in an NCI H226 lung mesothelioma mouse xenograft model when administered at a dose of 40 mg/kg.²

References

1. Hwang, S., Mruk, K., Rahighi, S., *et al.* Correcting glucose-6-phosphate dehydrogenase deficiency with a small-molecule activator. *Nat. Commun.* **9(1)**, 4045 (2018).
2. Lu, C., Yang, D., Klement, J.D., *et al.* H3K9me3 represses G6PD expression to suppress the pentose phosphate pathway and ROS production to promote human mesothelioma growth. *Oncogene* **41(18)**, 2651-2662 (2022).

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, 11/17/2023

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