

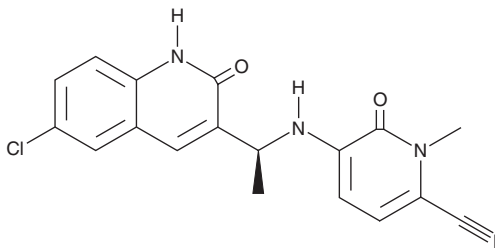
PRODUCT INFORMATION



Olutasidenib

Item No. 30370

CAS Registry No.: 1887014-12-1
Formal Name: 5-[[[(1S)-1-(6-chloro-1,2-dihydro-2-oxo-3-quinolinyl)ethyl]amino]-1,6-dihydro-1-methyl-6-oxo-2-pyridinecarbonitrile
Synonym: FT-2102
MF: C₁₈H₁₅ClN₄O₂
FW: 354.8
Purity: ≥98%
UV/Vis.: λ_{max}: 236, 335 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

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

Description

Olutasidenib is an inhibitor of mutant isocitrate dehydrogenase 1 (IDH1; IC₅₀s = 0.021 and 0.114 μM for IDH1^{R132H} and IDH1^{R132C}, respectively).¹ It is selective for mutant IDH1 over wild-type IDH1 and mutant IDH2 (IC₅₀s = 22.4, 27.3, and >100 μM for wild-type IDH1, IDH2^{R172K}, and IDH2^{R140Q}, respectively). Olutasidenib inhibits production of α-hydroxyglutaric acid (2-HG; Item Nos. 25894 | 16374) in U87 cells expressing mutant IDH1^{R132H}, IDH1^{R132C}, IDH1^{R132L}, IDH1^{R132G}, and IDH1^{R132S} with IC₅₀ values of 0.009, 0.039, 0.042, 0.006, and 0.009 μM, respectively. It reduces intratumor 2-HG levels in HCT116-IDH1^{R132H} and HCT116-IDH1^{R132C} mouse xenograft models when administered at doses of 12.5, 25, and 50 mg/kg.

Reference

1. Caravella, J.A., Lin, J., Diebold, R.B., *et al.* Structure-based design and identification of FT-2102 (olutasidenib), a potent mutant-selective IDH1 inhibitor. *J. Med. Chem.* **63**(4), 1612-1623 (2020).

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, 11/23/2022

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