

PRODUCT INFORMATION



TP-0463518

Item No. 33788

CAS Registry No.: 1558021-37-6

Formal Name: N-[[1-[[6-(4-chlorophenoxy)-3-pyridinyl]methyl]-1,2,5,6-tetrahydro-4-hydroxy-2-oxo-3-pyridinyl]carbonyl]-glycine

MF: C₂₀H₁₈ClN₃O₆

FW: 431.8

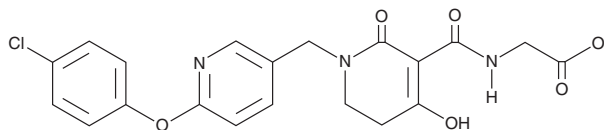
Purity: ≥98%

UV/Vis.: λ_{max}: 231, 274 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

TP-0463518 is supplied as a solid. A stock solution may be made by dissolving the TP-0463518 in the solvent of choice, which should be purged with an inert gas. TP-0463518 is soluble in organic solvents such as DMSO and dimethyl formamide (DMF). The solubility of TP-0463518 in these solvents is approximately 25 and 30 mg/ml, respectively.

TP-0463518 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, TP-0463518 should first be dissolved in DMF and then diluted with the aqueous buffer of choice. TP-0463518 has a solubility of approximately 0.3 mg/ml in a 1:2 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

TP-0463518 is a pan hypoxia-inducible factor prolyl hydroxylase (HIF-PH) inhibitor.¹ It inhibits human HIF-PH1 (K_i = 5.3 nM), as well as human HIF-PH2 and HIF-PH3 (IC₅₀s = 18 and 63 nM, respectively). TP-0463518 (5 mg/kg) increases serum levels of erythropoietin (EPO) in mice. It also increases serum EPO levels in a 5/6 nephrectomy-induced rat model of chronic kidney disease when administered at a dose of 10 mg/kg.

Reference

1. Kato, S., Takayama, N., Takano, H., *et al.* TP0463518, a novel inhibitor for hypoxia-inducible factor prolyl hydroxylases, increases erythropoietin in rodents and monkeys with a good pharmacokinetics-pharmacodynamics correlation. *Eur. J. Pharmacol.* **838**, 138-144 (2018).

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, 10/25/2022

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