

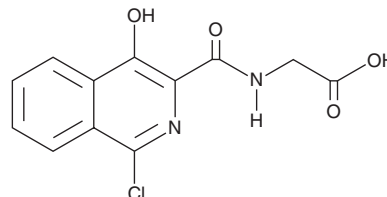
PRODUCT INFORMATION



FG-2216

Item No. 18382

CAS Registry No.: 223387-75-5
Formal Name: N-[(1-chloro-4-hydroxy-3-isoquinolinyl) carbonyl]-glycine
Synonyms: IOX3, YM-311
MF: C₁₂H₉ClN₂O₄
FW: 280.7
Purity: ≥98%
UV/Vis.: λ_{max}: 221, 257, 327, 341 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

FG-2216 is supplied as a crystalline solid. A stock solution may be made by dissolving the FG-2216 in the solvent of choice, which should be purged with an inert gas. FG-2216 is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of FG-2216 is approximately 0.5 mg/ml in ethanol and approximately 20 mg/ml in DMSO and DMF.

FG-2216 is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, FG-2216 should first be dissolved in DMSO and then diluted with the aqueous buffer of choice. FG-2216 has a solubility of approximately 0.5 mg/ml in a 1:1 solution of DMSO:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

Description

FG-2216 is an inhibitor of hypoxia-inducible factor prolyl hydroxylase 2 (HIF-PH2; IC₅₀ = 3.9 micromolar).¹ It has been shown to reversibly stimulate erythropoietin secretion at 100 μM *in vitro* and to increase hematocrit, red blood cell count, and hemoglobin levels at 50 mg/kg in mice.^{1,2}

References

1. Hong, Y.R., Kim, H.T., Lee, S.-L., *et al.* [(4-Hydroxyl-benzo[4,5]thieno[3,2-c]pyridine-3-carbonyl)-amino]-acetic acid derivatives; HIF prolyl 4-hydroxylase inhibitors as oral erythropoietin secretagogues. *Bioorg. Med. Chem.* **23**(21), 5953-5957 (2013).
2. Hsieh, M.M., Linde, N.S., Wynter, A., *et al.* HIF-prolyl hydroxylase inhibition results in endogenous erythropoietin induction, erythrocytosis, and modest fetal hemoglobin expression in rhesus macaque. *Blood* **110**(6), 2140-2147 (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, 10/31/2022

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