

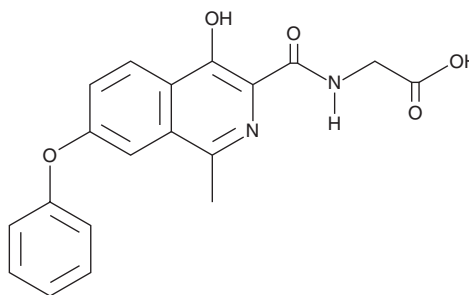
# PRODUCT INFORMATION



## Roxadustat

Item No. 15294

**CAS Registry No.:** 808118-40-3  
**Formal Name:** N-[(4-hydroxy-1-methyl-7-phenoxy-3-isoquinolinyl)carbonyl]-glycine  
**Synonyms:** ASP1517, FG-4592  
**MF:** C<sub>19</sub>H<sub>16</sub>N<sub>2</sub>O<sub>5</sub>  
**FW:** 352.4  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 223, 261, 314, 333 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

Roxadustat is supplied as a crystalline solid. A stock solution may be made by dissolving the roxadustat in the solvent of choice, which should be purged with an inert gas. Roxadustat is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide. The solubility of roxadustat in these solvents is approximately 0.1, 30, and 50 mg/ml, respectively.

### Description

Roxadustat is an analog of 2-OG that has been developed as an inhibitor of HIF-PHs. While little has been published regarding the actions of this compound, similar HIF-PH inhibitors suppress HIF degradation and in this way induce erythropoietin expression, promoting erythropoiesis or preventing anemia *in vivo*.<sup>1</sup>

### References

1. 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, 10/07/2022

#### 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