

# Roxadustat

## HIF-PHI inhibitor

Roxadustat is a small molecule inhibitor of hypoxia-inducible factor prolyl hydroxylase (HIF-PH), with an IC50 value of 591.4 nM. This compound is primarily used to treat anemia associated with chronic kidney disease (CKD) by stimulating erythropoiesis.

Key features and applications include:

- **HIF-PH Inhibition:** Inhibits HIF-PH enzymes, leading to increased levels of hypoxia-inducible factors (HIFs) and promoting erythropoietin production.
- **Erythropoiesis Stimulation:** Increases hemoglobin levels and red blood cell production. Used to treat symptomatic anemia in patients with chronic kidney disease, both dialysis-dependent and non-dialysis-dependent.
- **Cancer Research:** Investigated for its potential to reduce tumor growth by modulating the tumor microenvironment.

Relevant disease states include:

- **Chronic Kidney Disease (CKD):** Approved for the treatment of anemia associated with CKD.
- **Cancer:** Potential applications in reducing tumor growth and improving cancer treatment outcomes.

## Ordering Information

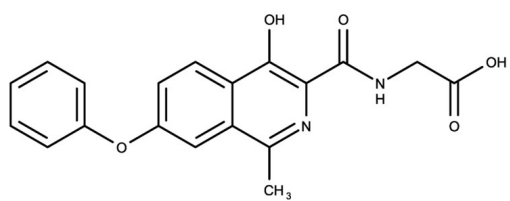
[Order Online »](#)

ENZ-CHM356-0025	25mg
-----------------	------

## Manuals, SDS & CofA

[View Online »](#)

- GMP format available



## Handling & Storage

Use/Stability	As indicated on product label or CoA when stored as recommended. Solutions in DMSO may be stored at -20°C for up to 2 months.
Handling	Keep container tightly closed in a dry and well-ventilated place.
Short Term Storage	-20°C
Long Term Storage	-20°C
Shipping	Ambient Temperature

## Regulatory Status

RUO - Research Use Only

## Product Details

Alternative Name	2-[(4-Hydroxy-1-methyl-7-phenoxyisoquinoline-3-carbonyl)amino]acetic acid
Appearance	Off-white or white solid.
CAS	808118-40-3
Couple Target	HIF-PH
Couple Type	Inhibitor
Formula	$C_{19}H_{16}N_2O_5$
Identity	Determined by NMR.
MW	352.35
Purity	≥99% (HPLC)
Solubility	Soluble in DMSO (up to at least 25 mg/ml).



ENZO LIFE SCIENCES,  
INC.  
Phone: 800.942.0430  
[info-  
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office  
ENZO LIFE SCIENCES  
(ELS) AG  
Phone: +41 61 926 8989  
[info-  
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands  
& Luxembourg  
Phone: +32 3 466 0420  
[info-  
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France  
Phone: +33 472 440 655  
[info-  
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany  
Phone: +49 7621 5500 526  
[info-  
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland  
Phone (UK customers):  
0845 601 1488  
Phone: +44 1392 825900  
[info-  
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)