

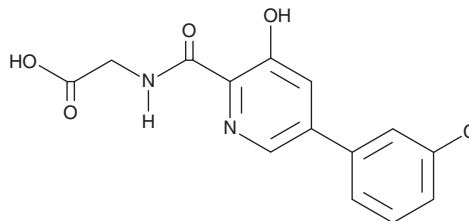
# PRODUCT INFORMATION



## Vadadustat

Item No. 15295

**CAS Registry No.:** 1000025-07-9  
**Formal Name:** N-[[5-(3-chlorophenyl)-3-hydroxy-2-pyridinyl]carbonyl]-glycine  
**Synonyms:** AKB6548, PG-1016548  
**MF:** C<sub>14</sub>H<sub>11</sub>ClN<sub>2</sub>O<sub>4</sub>  
**FW:** 306.7  
**Purity:** ≥98%  
**UV/Vis.:** λ<sub>max</sub>: 262, 317 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

Vadadustat is supplied as a crystalline solid. A stock solution may be made by dissolving the vadadustat in the solvent of choice, which should be purged with an inert gas. Vadadustat is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of vadadustat in ethanol and DMSO is approximately 2 mg/ml and approximately 5 mg/ml in DMF.

Vadadustat is sparingly soluble in aqueous buffers. For maximum solubility in aqueous buffers, vadadustat should first be dissolved in DMF and then diluted with the aqueous buffer of choice. Vadadustat has a solubility of approximately 0.04 mg/ml in a 1:20 solution of DMF:PBS (pH 7.2) using this method. We do not recommend storing the aqueous solution for more than one day.

### Description

Vadadustat is an inhibitor of hypoxia-inducible factor prolyl hydroxylase 2 (HIF-PH2) and HIF-PH3 (IC<sub>50</sub>s = 1.1 and 0.39 micromolar, respectively).<sup>1</sup> It stabilizes HIF-1α in mouse hepatocyte nuclear extracts. Vadadustat (50 and 100 mg/kg) increases erythropoietin levels in mouse serum.

### Reference

1. Kawamoto, R.M. Prolyl hydroxylase inhibitors and methods of use. *Procter & Gamble. US2007/0299086A1* (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, 12/09/2022

#### 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