

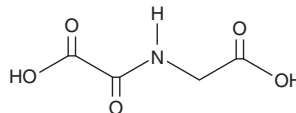
# PRODUCT INFORMATION



## N-Oxalylglycine

Item No. 13944

**CAS Registry No.:** 5262-39-5  
**Formal Name:** N-(carboxycarbonyl)-glycine  
**Synonym:** NOG  
**MF:** C<sub>4</sub>H<sub>5</sub>NO<sub>5</sub>  
**FW:** 147.1  
**Purity:** ≥98%  
**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

N-Oxalylglycine is supplied as a crystalline solid. A stock solution may be made by dissolving the N-oxalylglycine in the solvent of choice, which should be purged with an inert gas. N-Oxalylglycine is soluble in organic solvents such as ethanol, DMSO, and dimethyl formamide (DMF). The solubility of N-oxalylglycine in ethanol and DMSO is approximately 10 mg/ml and approximately 5 mg/ml in DMF.

Further dilutions of the stock solution into aqueous buffers or isotonic saline should be made prior to performing biological experiments. Ensure that the residual amount of organic solvent is insignificant, since organic solvents may have physiological effects at low concentrations. Organic solvent-free aqueous solutions of N-oxalylglycine can be prepared by directly dissolving the crystalline solid in aqueous buffers. The solubility of N-oxalylglycine in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

The jumonji domain-containing protein 2 (JMJD2) subfamily of histone demethylases have been shown to catalyze demethylation of the methylated forms of histone 3 lysine 9 (H3K9) and H3K36 *in vitro* and in cells.<sup>1</sup> Because histone demethylases are implicated in certain diseases, including cancer, selective inhibitors are candidate anticancer agents as well as potential tools for elucidating the biological functions of JMJDs.<sup>2</sup> N-Oxalylglycine, the amide analog of  $\alpha$ -ketoglutarate, is a cell permeable inhibitor of  $\alpha$ -ketoglutarate-dependent enzymes, including JMJD2A, JMJD2C, and JMJD2E (IC<sub>50</sub>s = 250, 500, and 24  $\mu$ M, respectively).<sup>3-6</sup> It can also inhibit the prolyl hydroxylase domain-containing proteins PHD1 and PHD2 with IC<sub>50</sub> values of 2.1 and 5.6  $\mu$ M, respectively.<sup>4-6</sup>

### References

1. Krishnan, S., Horowitz, S., and Trievel, R.C. *Chembiochem* **12**(2), 254-263 (2011).
2. Tian, X. and Fang, J. *Acta Biochim. Biophys. Sin. (Shanghai)* **39**(2), 81-88 (2007).
3. Hamada, S., Kim, T.D., Suzuki, T., et al. *Bioorg. Med. Chem. Lett.* **19**(10), 2852-2855 (2009).
4. Hamada, S., Suzuki, T., Mino, K., et al. *J. Med. Chem.* **53**(15), 5629-5638 (2010).
5. Rose, N.R., Ng, S.S., Mecinovic, J., et al. *J. Med. Chem.* **51**(22), 7053-7056 (2008).
6. Rose, N.R., Woon, E.C.Y., Kingham, G.L., et al. *J. Med. Chem.* **53**(4), 1810-1818 (2010).

#### 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, 02/03/2025

#### 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