

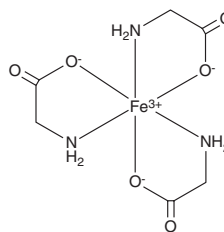
# PRODUCT INFORMATION



## Ferrous bisglycinate

Item No. 34128

**CAS Registry No.:** 34369-82-9  
**Formal Name:** tris(glycinato-κN,κO)-iron  
**Synonym:** Ferric glycinate  
**MF:** C<sub>6</sub>H<sub>12</sub>FeN<sub>3</sub>O<sub>6</sub>  
**FW:** 278.0  
**Purity:** ≥95%  
**Supplied as:** A 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

Ferrous bisglycinate is supplied as a solid. Aqueous solutions of ferrous bisglycinate can be prepared by directly dissolving the solid in aqueous buffers. The solubility of ferrous bisglycinate in PBS (pH 7.2) is approximately 10 mg/ml. We do not recommend storing the aqueous solution for more than one day.

### Description

Ferrous bisglycinate is an iron amino acid chelate.<sup>1,2</sup> Formulations containing ferrous bisglycinate have been used as dietary fortifying agents to prevent iron deficiency and anemia.

### References

1. Bovell-Benjamin, A.C., Viteri, R.E., and Allen, L.H. Iron absorption from ferrous bisglycinate and ferric trisglycinate in whole maize is regulated by iron status. *Am. J. Clin. Nutr.* **71(6)**, 1563-1569 (2000).
2. Pizzaro, F., Uicich, R., Olivaries, M., et al. Iron absorption of ferric glycinate is controlled by iron stores. *Nutr. Res.* **18(1)**, 3-9 (1998).

#### 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, 11/17/2022

#### 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