

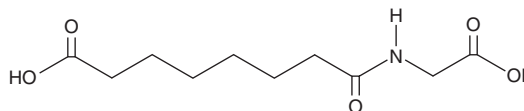
# PRODUCT INFORMATION



## Suberylglycine

Item No. 34619

CAS Registry No.: 60317-54-6  
Formal Name: 8-[(carboxymethyl)amino]-8-oxo-octanoic acid  
MF:  $C_{10}H_{17}NO_5$   
FW: 231.2  
Purity:  $\geq 95\%$   
Supplied as: A solid  
Storage:  $-20^{\circ}\text{C}$   
Stability:  $\geq 4$  years



Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Laboratory Procedures

Suberylglycine is supplied as a solid. A stock solution may be made by dissolving the suberylglycine in the solvent of choice, which should be purged with an inert gas. Suberylglycine is slightly soluble in DMSO and methanol.

### Description

Suberylglycine is a glycine-conjugated form of the dicarboxylic acid suberic acid.<sup>1</sup> Urinary levels of suberylglycine are increased in patients with medium-chain acyl-CoA dehydrogenase (MCAD) deficiency, an inborn error of fatty acid metabolism characterized by hypoketotic hypoglycemia, medium-chain dicarboxylic aciduria, and intolerance to fasting.<sup>2-4</sup>

### References

1. Truscott, R.J., Hick, L., Pullin, C., *et al.* Dicarboxylic aciduria: The response to fasting. *Clin. Chim. Acta* **94**(1), 31-39 (1979).
2. Bhuiyan, A.K., Watmough, N.J., Turnbull, D.M., *et al.* A new simple screening method for the diagnosis of medium chain acyl-CoA dehydrogenase deficiency. *Clin. Chim. Acta* **165**(1), 39-44 (1987).
3. Onkenhout, W., Venizelos, V., van der Poel, P.F.H., *et al.* Identification and quantification of intermediates of unsaturated fatty acid metabolism in plasma of patients with fatty acid oxidation disorders. *Clin. Chem.* **41**(10), 1467-1474 (1995).
4. Rinaldo, P., O'Shea, J.J., Coates, P.M., *et al.* Medium-chain acyl-CoA dehydrogenase deficiency. Diagnosis by stable-isotope dilution measurement of urinary n-hexanoylglycine and 3-phenylpropionylglycine. *N. Engl. J. Med.* **319**(20), 1308-1313 (1988).

#### 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, 11/10/2022

#### 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