

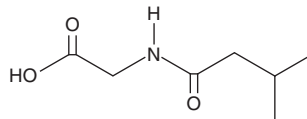
# PRODUCT INFORMATION



## Isovaleroylglycine

Item No. 17581

CAS Registry No.: 16284-60-9  
Formal Name: N-(3-methyl-1-oxobutyl)-glycine  
Synonym: N-Isovaleryl-glycine  
MF:  $C_7H_{13}NO_3$   
FW: 159.2  
Purity:  $\geq 95\%$   
Supplied as: A crystalline solid  
Storage:  $-20^{\circ}C$   
Stability:  $\geq 4$  years



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

### Laboratory Procedures

Isovaleroylglycine is supplied as a crystalline solid. A stock solution may be made by dissolving the isovaleroylglycine in the solvent of choice, which should be purged with an inert gas. Isovaleroylglycine is soluble in organic solvents such as methanol and DMSO. The solubility of isovaleroylglycine in these solvents is approximately 1 and 5 mg/ml, respectively.

### Description

Isovaleroylglycine is a compound found in the urine of patients with isovaleric acidemia, a disorder created by defective leucine metabolism due to a deficiency in the isovaleryl coenzyme A dehydrogenase.<sup>1</sup> It has been used as a diagnostic tool in the diagnosis of various acidemias and mitochondrial fatty acid  $\beta$ -oxidation defects.

### Reference

1. Tanaka, K. and Isselbacher, K.J. The isolation and identification of N-isovaleryl-glycine from urine of patients with isovaleric acidemia. *J. Biol. Chem.* **242**(12), 2966-2972 (1967).

#### 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, 11/11/2022

#### 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