

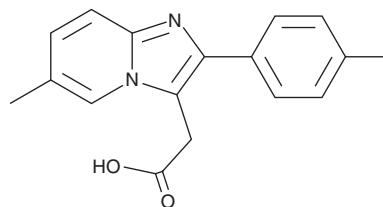
# PRODUCT INFORMATION



## Zolpidic Acid

Item No. 43000

CAS Registry No.: 189005-44-5  
Formal Name: 6-methyl-2-(4-methylphenyl)-imidazo[1,2-a]pyridine-3-acetic acid  
MF:  $C_{17}H_{16}N_2O_2$   
FW: 280.3  
Purity:  $\geq 98\%$   
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

Zolpidic acid is supplied as a solid. A stock solution may be made by dissolving the zolpidic acid in the solvent of choice, which should be purged with an inert gas. Zolpidic acid is sparingly soluble (1-10 mg/ml) in DMSO.

### Description

Zolpidic acid is an antioxidant and intermediate in the synthesis of the GABA<sub>A</sub> receptor positive allosteric modulator zolpidem.<sup>1</sup> It prevents iron sulfate-induced lipid peroxidation in rat liver and brain homogenates when used at a concentration of 1 mM.

### Reference

1. García-Santos, G. Herrera, F. Martín, V., *et al.* Antioxidant activity and neuroprotective effects of zolpidem and several synthesis intermediates. *Free Radic. Res.* **38(12)**, 1289-1299 (2004)

#### 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, 04/09/2025

#### 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