

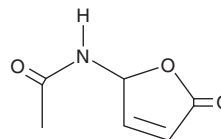
# PRODUCT INFORMATION



## 5-Acetamide-Butenolide

Item No. 40867

**CAS Registry No.:** 16275-44-8  
**Formal Name:** N-(2,5-dihydro-5-oxo-2-furanyl)-acetamide  
**Synonyms:** 4-acetamido-4-hydroxy-2-Butenoic Acid  $\gamma$ -lactone, NSC 114350  
**MF:**  $C_6H_7NO_3$   
**FW:** 141.1  
**Purity:**  $\geq 70\%$   
**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

5-Acetamide-butenolide is supplied as a solid. A stock solution may be made by dissolving the 5-acetamide-butenolide in the solvent of choice, which should be purged with an inert gas. 5-Acetamide-butenolide is soluble in ethanol, methanol, dichloromethane, and DMSO

### Description

5-Acetamide-butenolide is a mycotoxin that has been found in *Fusarium* and has pro-oxidant activity.<sup>1-3</sup> It disrupts the mitochondrial membrane potential in primary neonatal rat cardiomyocytes when used at a concentration of 50  $\mu\text{g}/\text{ml}$ .<sup>2</sup> 5-Acetamide-butenolide also induces the production of thiobarbituric acid reactive substances (TBARS) in isolated rat myocardial mitochondria in a concentration-dependent manner, an effect that can be reduced by glutathione (GSH). It increases the production of reactive oxygen species (ROS) and decreases the levels of GSH in, as well as reduces the viability of, HepG2 cells when used at concentrations ranging from 25 to 100  $\mu\text{g}/\text{ml}$ .<sup>3</sup> 5-Acetamide-butenolide is toxic to mice ( $\text{LD}_{50}$  = 43.6 mg/kg, i.p.).<sup>1</sup>

### References

1. Yates, A.S., Tookey, H.L., Ellis, J.J., *et al.* Mycotoxins produced by *Fusarium nivale* isolated from tall fescue (*Festuca arundinacea* Schreb.). *Phytochemistry* **7**(1), 139-146 (1968).
2. Wang, Y.-M., Liu, J.-B., and Peng, S.-Q. Effects of *Fusarium* mycotoxin butenolide on myocardial mitochondria in vitro. *Toxicol. Mech. Methods* **19**(2), 79-85 (2009).
3. Wang, Y.-M., Peng, S.-Q., Zhou, Q., *et al.* Depletion of intracellular glutathione mediates butenolide-induced cytotoxicity in HepG2 cells. *Toxicol. Lett.* **164**(3), 231-239 (2006).

#### 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, 04/23/2024

#### 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](http://WWW.CAYMANCHEM.COM)