

## Daminozide

Epigenetic Enzyme Inhibitor

### Catalog # K421-904B

Lot # P3320-14  
CAS # 1596-84-5

### Product Description

Molecular Formula: C<sub>6</sub>H<sub>12</sub>N<sub>2</sub>O<sub>3</sub>  
Appearance: White powder  
Melting Point: 158-162°C  
Molecular Weight: 160.2  
Purity: 98%; NMR (Conforms)  
Solubilization: May be dissolved in DMSO (50 mg/ml); or Water (20 mg/ml, warm)

### Alias

1-(2,2-Dimethylhydrazide)butanedioic acid

### Specific Activity

KDM2/7 Demethylase Inhibitor

### Storage and Stability

Store desiccated as supplied at ambient temperature for up to 3 years. Store solutions at -20°C for up to 3 months.

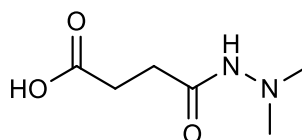
### Scientific Background

A selective inhibitor of the KDM2/7 JmjC subfamily of histone demethylases. IC<sub>50</sub>=1.5, 0.55 and 2.1 μM for KDM2A, PHF8 and KDM7A respectively. Highly selective over other subfamily members (KDM3, KDM4, KDM5 and KDM6) as well as other 2-oxoglutarate oxygenases (FIH, PHD2 and BBOX1), IC<sub>50</sub>>100 μM1.

### References

1. NR Rose et al. J. Med. Chem. 2012 55:6639

### Molecular Structure



## DAMINOZIDE

Epigenetic Enzyme Inhibitor

Catalog #	K421-904B
CAS #	1596-84-5
Lot #	P3320-14
Purity	98%
Format	White powder
Solubilization	May be dissolved in DMSO (50 mg/ml); or Water (20 mg/ml, warm)
Storage & Shipping	Store desiccated as supplied at ambient temperature for up to 3 years. Store solutions at -20°C for up to 3 months.

To place your order, please contact us by phone 1-(604)-232-4600, fax 1-604-232-4601 or by email: [orders@signalchem.com](mailto:orders@signalchem.com)  
[www.signalchem.com](http://www.signalchem.com)

**FOR IN VITRO RESEARCH PURPOSES ONLY. NOT INTENDED FOR USE IN HUMAN OR ANIMALS.**