

Daminozide

Catalog No: 14058, 14059

Format: 250 mg, 50 mg

Chemical Properties:

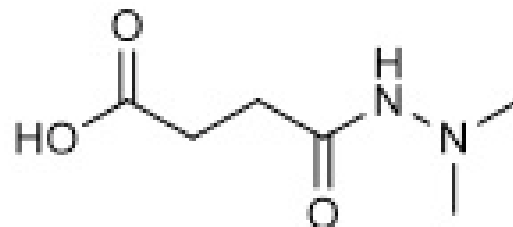
MW = 160.2

C₆H₁₂N₂O₃

CAS 1596-84-5

Physical Properties: White powder, MP = 158-162°C

Names: Daminozide, 1-(2,2-Dimethylhydrazide)butanedioic acid; N-(dimethylamino)succinamic acid



Pharmacology: A selective inhibitor of the KDM2/7 JmjC subfamily of histone demethylases. IC₅₀ = 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), IC₅₀ > 100 M (ref 1).

Solubilization: May be dissolved in DMSO (50 mg/ml); Water (20 mg/ml, warm)

Fluorescent Properties: N/A

Quality Control:

>98% (TLC); NMR (Conforms)

References:

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

Storage and Guarantee: Store desiccated as supplied at room temperature for up to 3 years. Store solutions at -20°C for up to 3 months. This product is guaranteed for 6 months from date of arrival.