

Oxamflatin

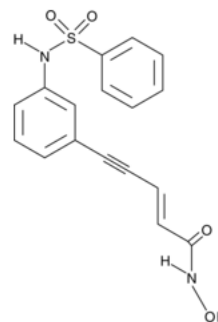
Catalog #: 27201

Lot #: 120203

Size: 1 mg

Structure:

CAS Registry #: 151720-43-3



Purity: ≥95%

Chemical Formula: C₁₇H₁₄N₂O₄S

Molecular Weight: 342.4

Description: Oxamflatin is a potent inhibitor of histone deacetylases. It has the ability to alter expression of genes involved in cell morphology, motility, apoptosis and cell cycle control, reducing the proliferation of cancer cells.

Appearance: White to off-white crystalline solid

Solubility: Solubility in ethanol, DMSO and DMF is 0.5 mg/ml, 5 mg/ml, and 10 mg/ml respectively. Sparingly soluble in aqueous buffers. For maximum solubility, dissolve first in DMF and then dilute with aqueous buffer of choice. Do not store in aqueous solution for more than one day.

Biological Activity: Oxamflatin inhibits histone deacetylases with IC₅₀ = 15.7 nM.

Storage/Stability: Store at or below -20 °C.

Quality Control: The purity was determined by HPLC analysis.

References:

1. Kim, Y.B., *et al.*, *Oncogene* 1999; **18**: 2461-2470.
2. Peart, M.J., *et al.*, *Cancer Res* 2003; **63**: 4460-4471.
3. Dear, A.E., *et al.*, *Biochem. Biophys. Acta* 2000; **1492**: 15-22.