Oxamflatin

HDAC inhibitor

Potent inhibitor of mammalian HDACs (histone deacetylases) (IC₅₀ =15.7nM). Acts as a ligand for the enzyme active site metal ion. Upregulates plasminogen activator inhibitor type 2 (PAI-2) expression with concomitant inhibition of urokinase (u-PA) gene and protein expression in HT-1080 and U-937 cells. Elevates the expression of extracellular matrix proteins fibronectin and gelsolin. Induces apoptosis in P-glyoprotein (Pgp) positive and Pgp negative multidrug resistant cells.

Citations: 7

View Online »

Ordering Information

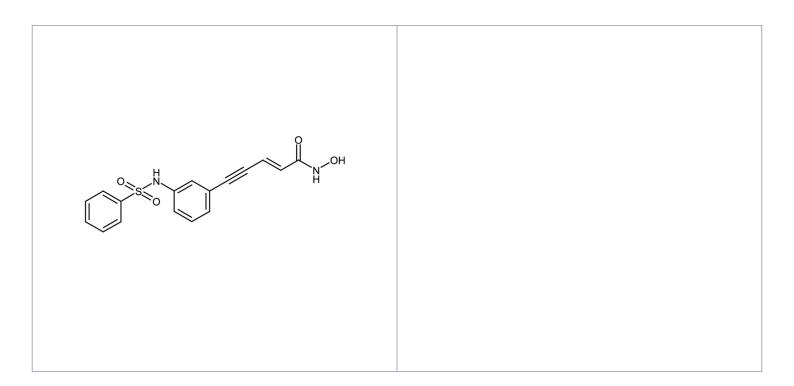
Order Online »

ALX-270-379-M005

5mg

Manuals, SDS & CofA

View Online »



Handling & Storage

Use/Stability As indicated on product label or CoA when stored as recommended.

Handling Protect from light. Packaged under inert gas. After reconstitution, prepare aliquots and

store at -20°C.

Long Term Storage +4°C

Blue Ice **Shipping**

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name (2E)-5-[3-(Phenylsulfonylamino)phenyl]pent-2-en-4-

ynohydroxamic acid

Appearance Tan to yellow solid.

CAS 151720-43-3

Couple Target HDAC

Couple Type Inhibitor

Formula $C_{17}H_{14}N_2O_4S$

342.4 MW

Purity ≥95% (HPLC)

Solubility Soluble in DMSO, methanol or acetonitrile.

Last modified: May 29, 2024

info-

