

Oxamflatin

HDAC inhibitor

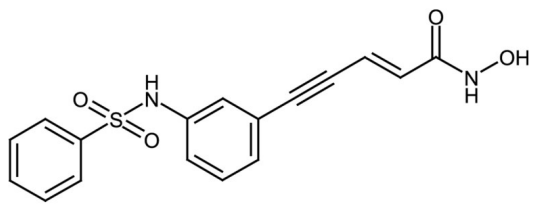
Potent inhibitor of mammalian HDACs (histone deacetylases) (IC_{50} =15.7nM). Acts as a ligand for the enzyme active site metal ion. Upregulates plasminogen activator inhibitor type 2 (PAI-2) expression with concomitant inhibition of urokinase (u-PA) gene and protein expression in HT-1080 and U-937 cells. Elevates the expression of extracellular matrix proteins fibronectin and gelsolin. Induces apoptosis in P-glycoprotein (Pgp) positive and Pgp negative multidrug resistant cells.

Citations: 7

[View Online »](#)

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
Shipping	Blue Ice

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	$C_{17}H_{14}N_2O_4S$
MW	342.4
Purity	≥95% (HPLC)
Solubility	Soluble in DMSO, methanol or acetonitrile.

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
info-usa@enzolifesciences.com

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
info-eu@enzolifesciences.com

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
info-fr@enzolifesciences.com

France
Phone: +33 472 440 655
info-de@enzolifesciences.com

Germany
Phone: +49 7621 5500 526
info-fr@enzolifesciences.com

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
info-uk@enzolifesciences.com