

HDAC1 (human), (recombinant) (His- tag)

Highly active

Human HDAC1 (HD1) was the first protein to be linked to histone deacetylase activity. It is homologous to the yeast protein Rpd31, a relationship which has since come to define the “class I HDACs”. HDAC1 promotes transcriptional repression by deacetylating lysine ϵ -amino groups in histone N-terminal tails, a function frequently carried out in association with multi-protein transcription repression complexes such as NuRD3, Sin34 and CoREST6. Ubiquitously expressed in human tissues HDAC1-containing complexes appear to contribute the greater part of (at least class I) deacetylase activity in HeLa nuclear extracts. Aside from its interaction with co-repressors, HDAC1 activity may be regulated by post-translation modifications such as phosphorylation⁹ and sumoylation or binding to the inhibitor maspin, a tumor-suppressive serpin homolog. Although originally described as a “histone deacetylase”, HDAC1 has been shown to catalyze the regulatory deacetylation of non-histone proteins, including p53. Overexpression of HDAC1 has been found in various cancer types. HDAC inhibitors (HDACi) have shown considerable promise as anti-cancer agents and HDACi compounds from multiple chemical classes are in stages of drug development ranging from preclinical to phase III trials.

Citations: 30

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE456-0050	50µg
----------------	------

Manuals, SDS & CofA

[View Online »](#)

Handling & Storage

Use/Stability	The enzyme is stable on ice for the time typically required to set up an experiment (30-60 min.), but may lose activity with prolonged storage on ice. It is recommended that thawing and dilution of the enzyme be done within as short a time as possible before start of the assay. The remaining, unused, undiluted enzyme should be refrozen quickly by, for example, snap freezing in a dry/ice ethanol bath or liquid nitrogen. Freezing and storage of diluted enzyme is not recommended.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Histone deacetylase 1
Formulation	Liquid. In 50mM TRIS, pH 8.0, 138mM sodium chloride and 10% glycerol.
Gene/Protein Identifier	NM_004964 (RefSeq)
MW	55 kDa
Purity Detail	Partially purified by single-step affinity chromatography and gel filtration.
Source	Produced in insect cells. HDAC1 from human cDNA (482 aa). Produced in a baculovirus expression system.
Specific Activity	≥5 U/μg. One U=1 pmol/min at 37°C, 100μM, FLUOR DE LYS [®] -SIRT1 deacetylase substrate (Prod. No. BML-KI177).
UniProt ID	Q13547

Last modified: May 29, 2024



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

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

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

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