

HDAC2 (full-length) (human), (recombinant) (His- tag)

Highly active HDAC2, like the highly homologous HDAC1, is a class I HDAC first identified as a human homolog of the yeast histone deacetylase Rpd3. HDAC2 found, along with HDAC1 in the Sin3, NuRD. In contrast to class II HDACs, which repress cardiac hypertrophy, HDAC2 activity is required for the cardiac hypertrophic response.

Citations: 2

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE533-0050	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 2
Formulation	Liquid. In 50mM TRIS, pH 8.0, 138 mM NaCl and 10% glycerol.
MW	~66 kDa
Purity Detail	Partially purified by single-step affinity chromatography and gel filtration.
Sequence	HDAC2 from human cDNA. Sequence is identical to residues 1-582 at Genbank accession# NM_00157.2.
Source	Produced in insect cells. Produced in a baculovirus expression system.
Specific Activity	≥0.5 U/μg. One U=1 pmol/min at 37°C, 100 μM, FLUOR DE LYS [®] -SIRT1 substrate (p53 K(Ac)-382 fluorogenic substrate, Prod. No. BML-KI177).
UniProt ID	Q92769

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)