

HDAC8 (human), (recombinant)

Citations: 3

[View Online »](#)

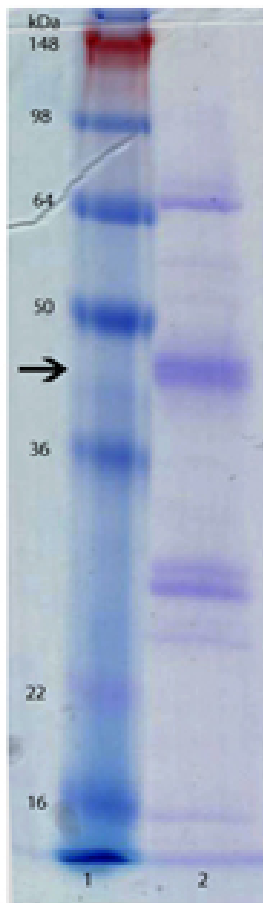
Ordering Information

[Order Online »](#)

BML-SE145-0100	100U
----------------	------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE Analysis: 1. MWM, 2. 2µg HDAC8 (human), (recombinant) (His-tag) (Prod. No. SE145).

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 dilution and/or prolonged storage on ice. It is recommended that thawing and addition of the enzyme to assay wells be done within as short a time as possible before start of the assay and that the thawed enzyme be kept on ice throughout. Best results will be obtained by adding the chilled, undiluted enzyme directly to prewarmed buffer and proceeding immediately to the addition of pre-warmed substrate. Either add a small volume of undiluted enzyme to each well (e.g. 1 U per well) or dilute into pre-warmed buffer (e.g. 1 U per 25 µl) and immediately aliquot to wells (see attached assay conditions). The remaining, unused enzyme should be refrozen quickly by, for example, snap freezing in a dry/ice ethanol bath or liquid nitrogen. The enzyme is stable to at least 4 freeze/thaw cycles. To minimize the number of freeze/thaw cycles, aliquot the HDAC8 into separate tubes and store at -70°C. NOTE: When stored under the above conditions, this enzyme is stable at the concentration supplied, in its current storage buffer. Procedures such as dilution of the enzyme followed by refreezing, could lead to loss of activity.
Long Term Storage	-80°C
Shipping	Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name	Histone deacetylase 8
Application Notes	Useful tool to study HDAC8 kinetics and inhibitor sensitivity and also the effect of the enzyme in transcriptional regulation, cell cycle progression, and oncogenesis.
Biological Activity	Like other class I HDACs, HDAC8 exhibits trichostatin A-inhibitable histone deacetylase activity and can mediate transcription repression.
Formulation	Liquid. In 10mM TRIS, pH 7.5, containing 100mM NaCl, 3mM MgCl ₂ and 10% glycerol.
Gene/Protein Identifier	NM_018486 (RefSeq)
MW	42 kDa
Purity Detail	Partially purified by single-step affinity chromatography and gel filtration.

Source	Produced in <i>E. coli</i> . Recombinant HDAC8 from human cDNA (377 aa)
Specific Activity	≥0.4 U/μg. One unit is defined as the amount of enzyme that deacetylates 1pmol/min of FLUOR DE LYS® deacetylase substrate (Prod. No. BML-KI104) at 37°C /250μM.
UniProt ID	Q9BY41

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)