

HDAC6 (human), (recombinant) (His- tag)

Highly active

HDAC6 is a class II HDAC, a group defined by its member's homology to the yeast HDAC HDA1. It is unique among human HDACs in that it contains two full deacetylase domains. A tubulin deacetylase, HDAC6 may act to promote autophagic clearing of huntingtin aggregates and retard HIV1 infection, but also to promote invasive motility in tumor cells.

Citations: 12

[View Online »](#)

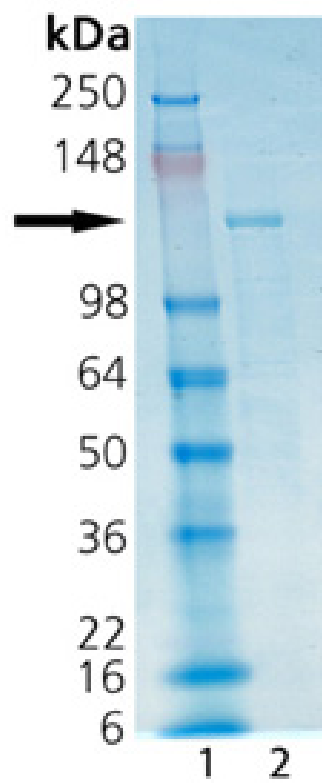
Ordering Information

[Order Online »](#)

BML-SE508-0050	50µg
----------------	------

Manuals, SDS & CofA

[View Online »](#)



SDS-PAGE analysis: Lane 1: MW marker, Lane 2: 2 μ g HDAC6.

Handling & Storage

Long Term Storage -80°C

Shipping Dry Ice

Regulatory Status

RUO - Research Use Only

Product Details

Alternative Name Histone deacetylase 6

Biological Activity Active with various fluorogenic peptide substrates, including FLUOR DE LYS[®] deacetylase substrate (Prod. No. BML-KI104) and FLUOR DE LYS[®]-SIRT1 deacetylase substrate (Prod. No. BML-KI177).

Formulation Liquid. In 50mM TRIS, pH 8.0, 138 mM NaCl and 10% glycerol.

Gene/Protein Identifier NM_006044 (RefSeq)

MW 131 kDa

Purity Detail Partially purified by single-step affinity chromatography and gel filtration.

Source Produced in insect cells. Produced in a baculovirus expression system.

Specific Activity ≥1 U/μg. One U=1 pmol/min at 37°C, 100 μM, FLUOR DE LYS[®]-SIRT1 deacetylase substrate (Prod. No. BML-KI177).

UniProt ID Q9UBN7

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-be@enzolifesciences.com)

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