

# HDAC3 (human), (recombinant) (His- tag)

Highly active

HDAC3, although a class I HDAC and has been assigned to its own phylogenetic subclass and its caspase-dependent cleavage and relocation to the cytoplasm may be critical to the progression of apoptosis, implying an important role in pathways affecting heart disease.

Citations: 6

[View Online »](#)

Ordering Information

[Order Online »](#)

BML-SE507-0050	50µg
----------------	------

Manuals, SDS & CofA

[View Online »](#)

## Handling & Storage

Long Term Storage      -80°C

Shipping      Dry Ice

**Regulatory Status**      RUO - Research Use Only

## Product Details

Alternative Name      Histone deacetylase 3

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

Gene/Protein Identifier      NM\_003883 (RefSeq)

MW      49 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       $\geq 0.7$  U/ $\mu$ g. One U=1 pmol/min at 37°C, 100  $\mu$ M, FLUOR DE LYS<sup>®</sup>-SIRT1 deacetylase substrate (Prod. No. BML-K1177).

UniProt ID      O15379



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)