

PRODUCT INFORMATION



HDAC9 (human, recombinant)

Item No. 10009466

Overview and Properties

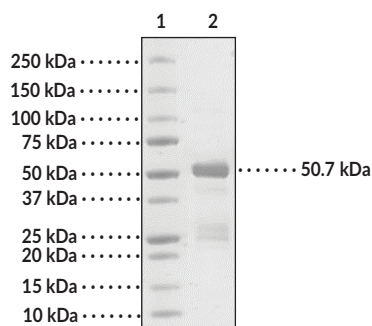
Synonyms: Histone Deacetylase 9
Source: 10 µg of recombinant C-terminal His-tagged protein expressed in baculovirus expression system
Amino Acids: 604-1066
Molecular Weight: 50.7 kDa
Storage: -80°C (as supplied)
Stability: ≥6 months
Purity: ≥75% estimated by SDS-PAGE
Supplied in: 40 mM Tris-HCl, pH 8.0, with 110 mM sodium chloride, 2.2 mM potassium chloride, 20% glycerol, and 200 mM imidazole

Protein

Concentration: *batch specific* mg/ml
Specific Activity: *batch specific* U/mg
Unit Definition: One unit is the amount of enzyme required to release 1 pmol of acetate per minute at 37°C in 25 mM Tris-HCl, pH 8.0, 137 mM sodium chloride, 2.7 mM potassium chloride, 1 mM magnesium chloride, 0.1 mg/ml BSA, and 20 µM fluorogenic HDAC class 2a substrate.

Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

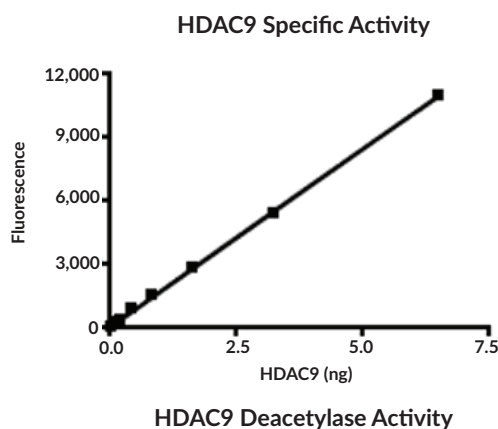
Images



Lane 1: MW Markers
Lane 2: HDAC9 (3.15 µg)

SDS-PAGE Analysis of HDAC9.

Representative gel image shown; actual purity may vary between each batch.



WARNING
THIS PRODUCT IS FOR RESEARCH ONLY - NOT FOR HUMAN OR VETERINARY DIAGNOSTIC OR THERAPEUTIC USE.

SAFETY DATA
This material should be considered hazardous until further information becomes available. Do not ingest, inhale, get in eyes, on skin, or on clothing. Wash thoroughly after handling. Before use, the user must review the [complete](#) Safety Data Sheet, which has been sent via email to your institution.

WARRANTY AND LIMITATION OF REMEDY
Buyer agrees to purchase the material subject to Cayman's Terms and Conditions. Complete Terms and Conditions including Warranty and Limitation of Liability information can be found on our website.

Copyright Cayman Chemical Company, 05/10/2022

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM

PRODUCT INFORMATION



Description

HDAC9 is a Class IIa HDAC which is homologous to the yeast HDAC1 and is larger in size than the other classes of HDACs.^{1,2} Class IIa HDACs contain a highly conserved C-terminal deacetylase catalytic domain (~420 amino acids) and N-terminal domain with no similarity to HDACs in other classes. Class II HDACs can shuttle between the nucleus and cytoplasm, suggesting possible extranuclear functions including regulating the acetylation status of non-histone substrates. By modifying chromatin structure and other non-histone proteins, HDACs play an important role in controlling complex biological events, including cell development, differentiation, programmed cell death, angiogenesis, and inflammation.^{1,2} Considering these major roles, it is conceivable that dysregulation of HDACs and subsequent imbalance of acetylation and deacetylation may be involved in the pathogenesis of various diseases, including cancer and inflammatory diseases.²

References

1. Lin, H.Y., Chen, C.S., Lin, S.P., *et al.* Targeting histone deacetylase in cancer therapy. *Med. Res. Rev.* **26(4)**, 397-413 (2006).
2. Huang, L. Targeting histone deacetylases for the treatment of cancer and inflammatory diseases. *J. Cell. Physiol.* **209(3)**, 611-616 (2006).

CAYMAN CHEMICAL
1180 EAST ELLSWORTH RD
ANN ARBOR, MI 48108 · USA
PHONE: [800] 364-9897
[734] 971-3335
FAX: [734] 971-3640
CUSTSERV@CAYMANCHEM.COM
WWW.CAYMANCHEM.COM