

# PRODUCT INFORMATION



## HDAC4 (human, recombinant)

Item No. 10009652

### Overview and Properties

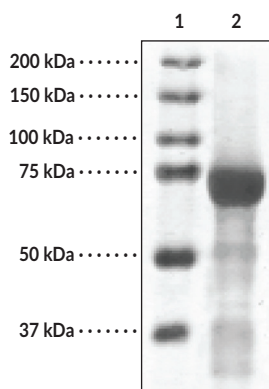
**Synonym:** Histone Deacetylase 4  
**Source:** Active recombinant human N-terminal GST-tagged HDAC4 catalytic domain expressed in insect cells  
**Amino Acids:** 627-1,085  
**Uniprot No.:** P56524  
**Molecular Weight:** 75.2 kDa  
**Storage:** -80°C (as supplied)  
**Stability:** ≥6 months  
**Purity:** ≥60% estimated by SDS-PAGE  
**Supplied in:** 45 mM Tris-HCl, pH 8.0, with 124 mM sodium chloride, 2.4 mM potassium chloride, 18 mM glutathione, and 10% glycerol

### Protein

**Concentration:** *batch specific* mg/ml  
**Activity:** *batch specific* U/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 MgCl<sub>2</sub>, 0.1 mg/ml BSA, and 20 μM fluorogenic HDAC class 2a substrate.

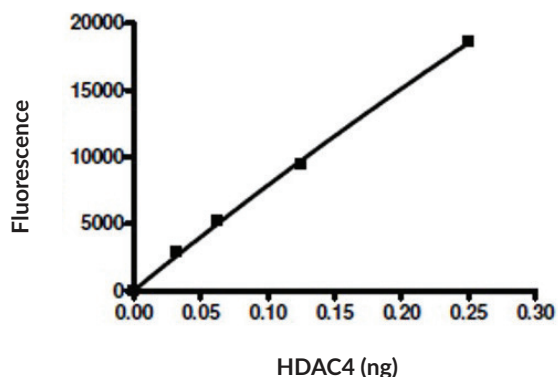
Information represents the product specifications. Batch specific analytical results are provided on each certificate of analysis.

### Images



Lane 1: MW Markers  
Lane 2: HDAC4 (5 μg)

SDS-PAGE Analysis of HDAC4



HDAC4 Deacetylase Activity

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, 04/04/2023

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

---

Histone deacetylase 4 (HDAC4) is a zinc-dependent metalloenzyme and class IIa HDAC.<sup>1</sup> It is composed of an N-terminal regulatory domain, which contains a myocyte-specific enhancer factor 2 (MEF2) binding site, three 14-3-3 binding sites, a nuclear localization signal, and a caspase-3 cleavage site, a catalytic domain, and a C-terminal domain that contains a nuclear export signal. HDAC4 shuttles between the cytoplasm and nucleus and is mainly expressed in skeletal muscle, brain, ovaries, and colon but is also found in the small intestine, heart, kidney, testis, thymus, and leukocytes.<sup>2,3</sup> It acts as a transcriptional corepressor and has many binding partners, including the transcription factors MEF2 and RUNX family transcription factor 2 (RUNX2).<sup>1,4</sup> Knockout of *Hdac4* induces premature ossification and chondrocyte hypertrophy in mice.<sup>4</sup> Mutations in *HDAC4* are associated with brachydactyly-mental retardation syndrome, a disorder characterized by brachydactyly type E, intellectual disability, and facial dysmorphisms.<sup>5</sup> Cayman's HDAC4 (human, recombinant) protein can be used for enzyme activity assays.

## References

---

1. Yang, X.J. and Grégoire, S. Class II histone deacetylases: From sequence to function, regulation, and clinical implication. *Mol. Cell. Biol.* **25(8)**, 2873-2884 (2005).
2. Wang, A.H., Kruhlak, M.J., Wu, J., *et al.* Regulation of histone deacetylase 4 by binding of 14-3-3 proteins. *Mol. Cell. Biol.* **20(18)**, 6904-6912 (2000).
3. Wang, A.H., Bertos, N.R., Vezmar, M., *et al.* HDAC4, a human histone deacetylase related to yeast HDA1, is a transcriptional corepressor. *Mol. Cell. Biol.* **19(11)**, 7816-7827 (1999).
4. Vega, R.B., Matsuda, K., Oh, J., *et al.* Histone deacetylase 4 controls chondrocyte hypertrophy during skeletogenesis. *Cell* **119(4)**, 555-566 (2004).
5. Wakeling, E., McEntagart, M., Bruccoleri, M., *et al.* Missense substitutions at a conserved 14-3-3 binding site in HDAC4 cause a novel intellectual disability syndrome. *HGG Adv.* **2(1)**, 100015 (2021).

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