

PRODUCT INFORMATION



HDAC8 (human, recombinant)

Item No. 19380

Overview and Properties

Synonym:	Histone Deacetylase 8
Source:	Active recombinant C-terminal hexahistidine-tagged protein expressed in <i>E. coli</i>
Amino Acids:	2-377 (full length)
Uniprot No.:	Q9BY41
Molecular Weight:	45.3 kDa
Storage:	-80°C (as supplied); avoid freeze/thaw cycles by aliquoting protein
Stability:	≥6 months
Purity:	≥80% estimated by SDS-PAGE
Supplied in:	10 mM Tris, pH 7.5, containing 100 mM NaCl, 3 mM MgCl ₂ , and 20% glycerol
Protein Concentration:	<i>batch specific</i> mg/ml
Activity:	<i>batch specific</i> U/ml
Specific Activity:	<i>batch specific</i> U/mg
Unit Definition:	One unit is the amount of enzyme required to release 1 nmol/min acetate from 100 μM acetylated p53 peptide (Item No. 10010995) at 37°C in 25 mM Tris, pH 8.0, 137 mM NaCl, 2.7 mM KCl, and 1 mM MgCl ₂ .

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

Images

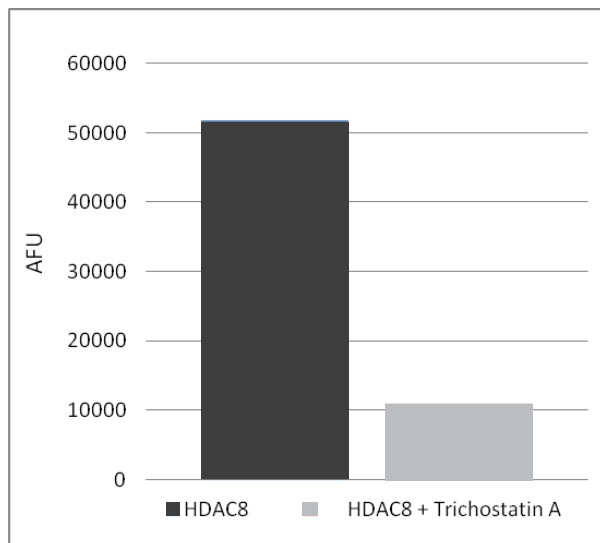
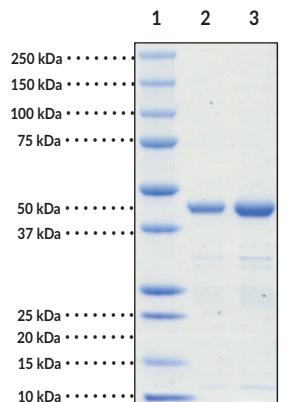


Figure 1: HDAC8 activity was inhibited by a known inhibitor Trichostatin A.



Lane 1: MW Markers
Lane 2: HDAC8 (2 μg)
Lane 3: HDAC8 (4 μg)

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, 06/21/2023

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 8 (HDAC8) is a class I HDAC that catalyzes the zinc-dependent deacetylation of core histones.^{1,2} It is composed of a single α/β domain with an N-terminal deacetylase active site, which can be disrupted by phosphorylation of a proximal serine.³ HDAC8 is ubiquitously expressed in human tissues and, unlike most other HDACs, is localized to both the cytosol and nucleus.⁴ It associates with other proteins and transcription factors to regulate gene expression. HDAC8 has a flexible active site that can deacetylate lysines from both histone tails and cytosolic proteins.³ It deacetylates pyruvate kinase M2 (PKM2), inducing its relocalization to the nucleus where it binds β -catenin and promotes cyclin D1 expression and cell cycle progression in HepG2 hepatocellular carcinoma (HCC) cells.⁵ HDAC8 associates with RUNX family transcription factor 1 (RUNX1) and suppresses the expression of RUNX1-regulated genes, which code for proteins involved in cytokine expression, cell differentiation, cell cycle progression, and tumor suppression, in COS-7 cells.⁶ Inhibition of HDAC8 decreases tumor volume and increases intratumoral apoptosis without reducing body weight in a neuroblastoma mouse xenograft model.⁷ Cayman's HDAC8 (human, recombinant) protein can be used for enzyme activity assay and Western blot applications.

References

1. Lin, H.-Y., Chen, C.-S., Lin, S.-P., *et al.* Targeting histone deacetylase in cancer therapy. *Medicinal Research Reviews* **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).
3. Somoza, J.R., Skene, R.J., Katz, B.A., *et al.* Structural snapshots of human HDAC8 provide insights into the class I histone deacetylases. *Structure* **12(7)**, 1325-1334 (2004).
4. Xia, J.-K., Qin, X.-Q., Zhang, L., *et al.* Roles and regulation of histone acetylation in hepatocellular carcinoma. *Front. Genet.* **13**, 982222 (2022).
5. Zhang, R., Shen, M., Wu, C., *et al.* HDAC8-dependent deacetylation of PKM2 directs nuclear localization and glycolysis to promote proliferation in hepatocellular carcinoma. *Cell Death Dis.* **11(12)**, 1036 (2020).
6. Durst, K.L., Lutterbach, B., Kummalue, T., *et al.* The *inv(16)* fusion protein associates with corepressors via a smooth muscle myosin heavy-chain domain. *Mol. Cell Biol.* **23(2)**, 607-619 (2003).
7. Rettig, I., Koeneke, E., Trippel, F., *et al.* Selective inhibition of HDAC8 decreases neuroblastoma growth *in vitro* and *in vivo* and enhances retinoic acid-mediated differentiation. *Cell Death Dis.* **6(2)**, e1657 (2015).

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