

Human HDAC8 ELISA Kit ab282872

重组 SimpleStep ELISA

5 图像

概述

产品名称	人HDAC8 ELISA试剂盒				
检测方法	Colorimetric				
精确度	批次内				
	样品	n	Mean	SD	CV%
	Extract	8			1.9%
	批次间				
	样品	n	Mean	SD	CV%
	Extract	0			0%
样品类型	Cell Lysate				
检测类型	Sandwich (quantitative)				
灵敏度	3.5 pg/ml				
范围	46.88 pg/ml - 3000 pg/ml				
回收率	特定样本回收率				
	样品类型	平均%		范围	
	Cell Lysate	101		98% - 107%	
检测时间	1h 30m				
实验步骤	One step assay				
种属反应性	与反应: Human				
产品概述	Human HDAC8 ELISA kit (ab282872) is a single-wash 90 min sandwich ELISA designed for the				

SimpleStep ELISA® technology employs capture antibodies conjugated to an affinity tag that is

recognized by the monoclonal antibody used to coat our SimpleStep ELISA® plates. This approach to sandwich ELISA allows the formation of the antibody-analyte sandwich complex in a single step, significantly reducing assay time. See the SimpleStep ELISA® protocol summary in the image section for further details. Our SimpleStep ELISA® technology provides several benefits:

- Single-wash protocol reduces assay time to 90 minutes or less
- High sensitivity, specificity and reproducibility from superior antibodies
- Fully validated in biological samples
- 96-wells plate breakable into 12 x 8 wells strips

A 384-well SimpleStep ELISA® microplate (**ab203359**) is available to use as an alternative to the 96-well microplate provided with SimpleStep ELISA® kits.

说明

HDAC8 (Histone deacetylase 8) is a nuclear enzyme responsible for the deacetylation of lysine residues on the N-terminal part of the core histones (H2A, H2B, H3 and H4). Histone deacetylation gives a tag for epigenetic repression and plays an important role in transcriptional regulation, cell cycle progression and developmental events. Histone deacetylases act via the formation of large multiprotein complexes. HDAC8 is also involved in the deacetylation of cohesin complex protein SMC3 regulating release of cohesin complexes from chromatin. Mouse and rat HDAC8 share 96% amino acid sequence identity with human HDAC8 protein.

Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.

It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.

平台

Pre-coated microplate (12 x 8 well strips)

性能

存放说明

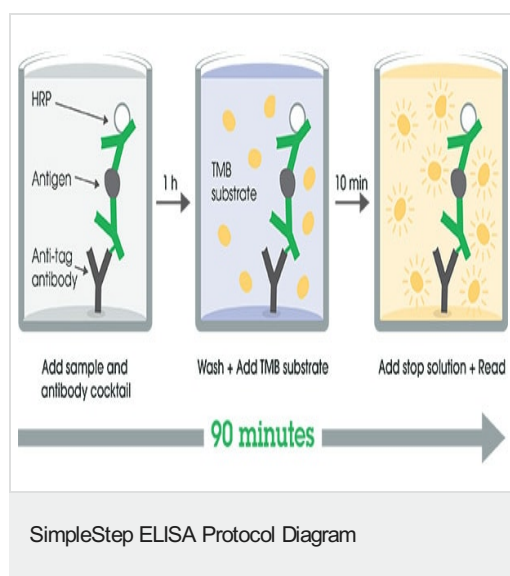
Store at +4°C. Please refer to protocols.

组件	1 x 96 tests
10X Human HDAC8 Capture Antibody	1 x 600µl
10X Human HDAC8 Detector Antibody	1 x 600µl
10X Wash Buffer PT (ab206977)	1 x 20ml
50X Cell Extraction Enhancer Solution (ab193971)	1 x 1ml
5X Cell Extraction Buffer PTR (ab193970)	1 x 10ml
Antibody Diluent CPI2	1 x 6ml
Human HDAC8 Lyophilized Recombinant Protein	2 vials
Plate Seals	1 unit
Sample Diluent NS (ab193972)	1 x 12ml

组件	1 x 96 tests
SimpleStep Pre-Coated 96-Well Microplate (ab206978)	1 unit
Stop Solution	1 x 12ml
TMB Development Solution	1 x 12ml

功能	Responsible for the deacetylation of lysine residues on the N-terminal part of the core histones (H2A, H2B, H3 and H4). Histone deacetylation gives a tag for epigenetic repression and plays an important role in transcriptional regulation, cell cycle progression and developmental events. Histone deacetylases act via the formation of large multiprotein complexes. May play a role in smooth muscle cell contractility.
组织特异性	Weakly expressed in most tissues. Expressed at higher level in heart, brain, kidney and pancreas and also in liver, lung, placenta, prostate and kidney.
序列相似性	Belongs to the histone deacetylase family. HD type 1 subfamily.
翻译后修饰	Phosphorylated by PKA on serine 39. Phosphorylation reduces deacetylase activity observed preferentially on histones H3 and H4.
细胞定位	Nucleus. Cytoplasm. Excluded from the nucleoli. Found in the cytoplasm of cells showing smooth muscle differentiation.

图片



SimpleStep ELISA technology allows the formation of the antibody-antigen complex in one single step, reducing assay time to 90 minutes. Add samples or standards and antibody mix to wells all at once, incubate, wash, and add your final substrate. See protocol for a detailed step-by-step guide.

Powered by
recombinant antibodies



**Research with
confidence**
Consistent and
reproducible results



**Long-term and
scalable supply**
Recombinant
technology



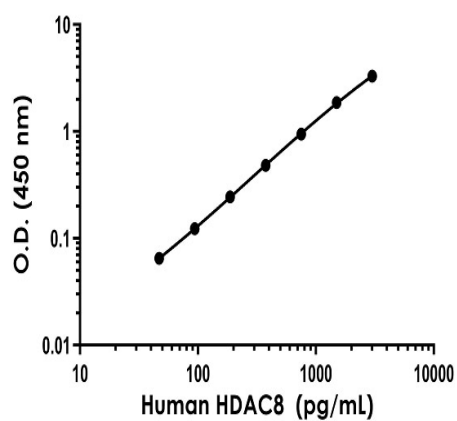
**Success from the
first experiment**
Confirmed
specificity



**Ethical standards
compliant**
Animal-free
production

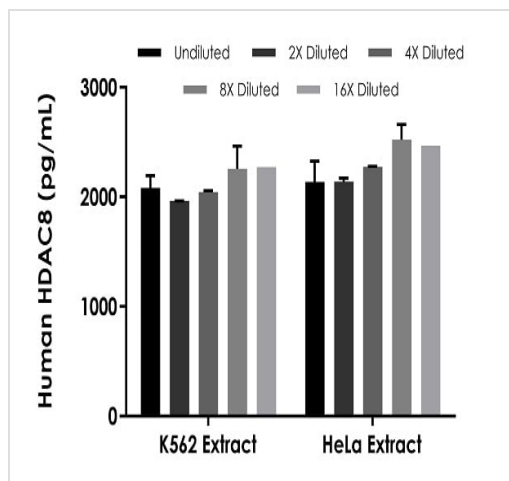
Recombinant Antibody Benefits

To learn more about the advantages of recombinant antibodies see [here](#).



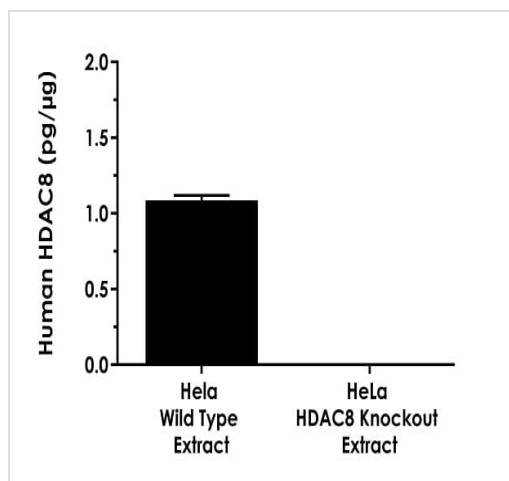
Example of human HDAC8 standard curve in 1X Cell
Extraction Buffer PTR.

The HDAC8 standard curve was prepared as described in Section 10. Raw data values are shown in the table. Background-subtracted data values (mean \pm SD) are graphed.



Interpolated concentrations of native HDAC8 in human K562 cell and HeLa cell based on 250 µg/mL and 500 µg/mL extract loads, respectively.

The concentrations of HDAC8 were measured in duplicate and interpolated from the HDAC8 standard curve and corrected for sample dilution. The interpolated dilution factor corrected values are plotted (mean \pm SD, n=2). The mean HDAC8 concentration was determined to be 2,107 pg/mL in K562 extract and 2,289 pg/mL in HeLa extract.



Interpolated concentrations of native HDAC8 in human control wild type HeLa cell and HDAC8 knockout HeLa cell based on 400 µg/mL extract loads.

The concentrations of HDAC8 were measured in duplicate and interpolated from the HDAC8 standard curve and expressing the interpolated concentrations as pg of protein per µg of extract (mean \pm SD, n=2). The mean HDAC8 concentration was determined to be 1.09 pg/µg in wild type HeLa extract and undetectable in HDAC8 knockout HeLa extract.

Please note: All products are "FOR RESEARCH USE ONLY. NOT FOR USE IN DIAGNOSTIC PROCEDURES"

Our Abpromise to you: Quality guaranteed and expert technical support

- Replacement or refund for products not performing as stated on the datasheet
- Valid for 12 months from date of delivery
- Response to your inquiry within 24 hours
- We provide support in Chinese, English, French, German, Japanese and Spanish
- Extensive multi-media technical resources to help you
- We investigate all quality concerns to ensure our products perform to the highest standards

If the product does not perform as described on this datasheet, we will offer a refund or replacement. For full details of the Abpromise, please visit <https://www.abcam.cn/abpromise> or contact our technical team.

Terms and conditions

- Guarantee only valid for products bought direct from Abcam or one of our authorized distributors