

Human Vimentin Profiling ELISA Kit ab173190

1 References 3 图像

概述

产品名称

检测方法

精确度

人Vimentin Profiling ELISA试剂盒

Colorimetric

批次内

样品	n	Mean	SD	CV%
Overall	3			7.5%

批次间

样品	n	Mean	SD	CV%
Overall	3			12%

样品类型

检测类型

灵敏度

回收率

Adherent cells, Suspension cells, Tissue Homogenate, Tissue Lysate

Quantitative

8 µg/ml

> 80 %

特定样本回收率		
样品类型	平均%	范围
Cell culture media	80	71% - 87%
Fetal Bovine Serum	89	81% - 93%
Bovine Serum Albumin	89	83% - 84%

实验步骤	Multiple steps standard assay
种属反应性	与反应: Human
产品概述	Abcam's Vimentin Human Profiling <i>in vitro</i> ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for the accurate qualitative measurement of total vimentin protein in Human samples.

The assay employs an antibody specific to Vimentin coated onto well plate strips. Samples are pipetted into the wells and Vimentin present in the sample is bound to the wells by the

说明

平台

性能

存放说明

功能

组织特异性

疾病相关

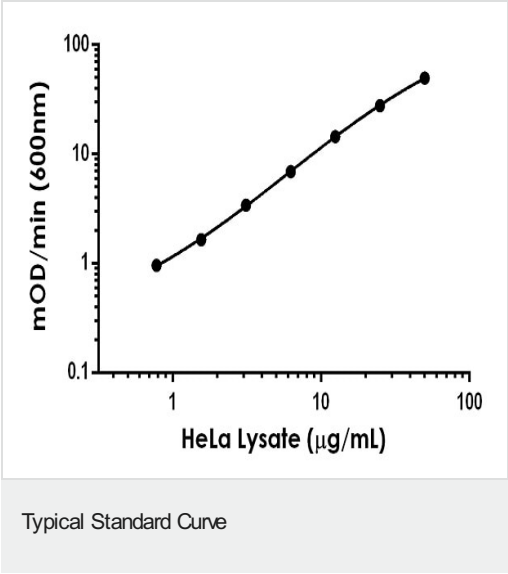
序列相似性

结构域

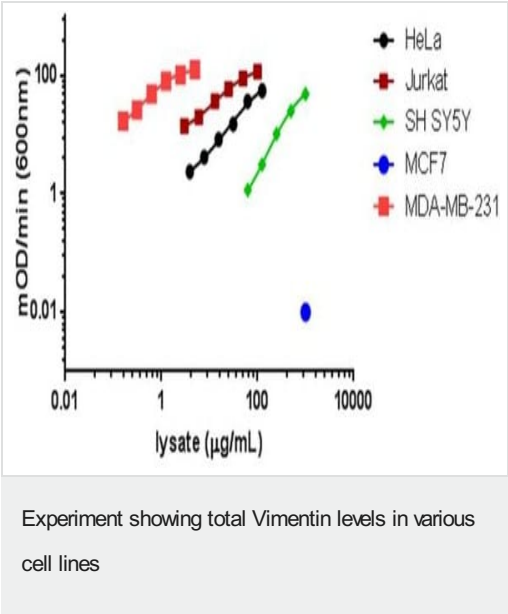
The [IL]-x-C-x-x-[DE] motif is a proposed target motif for cysteine S-nitrosylation mediated by the

	iNOS-S100A8/A9 transnitrosylase complex.
翻译后修饰	<p>Filament disassembly during mitosis is promoted by phosphorylation at Ser-55 as well as by nestin (By similarity). One of the most prominent phosphoproteins in various cells of mesenchymal origin. Phosphorylation is enhanced during cell division, at which time vimentin filaments are significantly reorganized. Phosphorylation by PKN1 inhibits the formation of filaments. Phosphorylated at Ser-56 by CDK5 during neutrophil secretion in the cytoplasm. Phosphorylated by STK33.</p> <p>O-glycosylated during cytokinesis at sites identical or close to phosphorylation sites, this interferes with the phosphorylation status.</p> <p>S-nitrosylation is induced by interferon-gamma and oxidatively-modified low-density lipoprotein (LDL(ox)) possibly implicating the iNOS-S100A8/9 transnitrosylase complex.</p>
细胞定位	Cytoplasm.

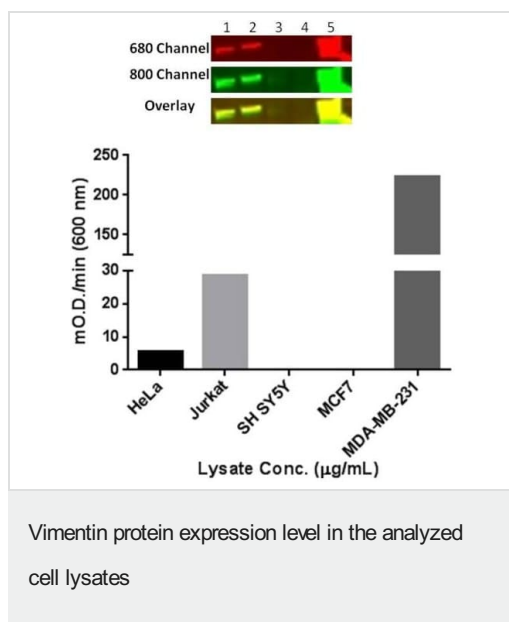
图片



Typical standard curve generated when using the Vimentin Human Profiling ELISA kit.



HeLa cell lysate, Jurkat cell lysate, SH SY5Y cell lysate, MCF7 cell lysate (not shown) and MDA-MB-231 cell lysate were tested using this kit. HeLa cell lysate shown as standard control sample, varying concentration other cell Lysates were analyzed within working range of the assay.



Western blot of cell lysates (top panel): 1) HeLa, 2) Jurkat, 3) Sy5y, 4) MCF7, 5) MDA-MB-231 (10 µg/lane) using primary antibodies: rabbit anti-vimentin (1/1000 dilution) and mouse anti-vimentin (**ab8978**, 1/1000 dilution); Secondary antibodies used were goat anti-rabbit 680-RD (Red, 1/10,000) and goat anti-mouse 800 (Green, 1/10,000). Blot was scanned using a LI-COR® Odyssey® imager and overlay shows target specificity for total vimentin (yellow). Bar graph (bottom panel) represents relative mOD/min (600 nm) observed for 10 µg/mL cell lysate, except for MDA-MB-231 cell lysate which reports mOD/min (600 nm) for 1 µg/mL.

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