

Human Caspase-9 ELISA Kit ab119508

4 References 1 图像

概述

产品名称	人Caspase-9 ELISA试剂盒			
检测方法	Colorimetric			
精确度	批次内			
	样品	n	Mean	SD
	Overall	8		6.6%
	批次间			
	样品	n	Mean	SD
	Overall	8		9%
样品类型	Cell culture supernatant, Serum, Cell Lysate			
检测类型	Sandwich (quantitative)			
灵敏度	0.4 ng/ml			
范围	1.6 ng/ml - 100 ng/ml			
回收率	103 %			
实验步骤	Multiple steps standard assay			
种属反应性	与反应: Human			
产品概述	Abcam's Caspase-9 Human in vitro ELISA (Enzyme-Linked Immunosorbent Assay) kit is designed for accurate quantitative measurement of Human Caspase-9 concentrations in cell lysate, cell culture supernatant and serum.			
	Caspase-9 specific antibodies have been precoated onto 96-well plates. Standards and test samples are added to the wells and along with a Caspase-9 detection antibody and the microplate is then incubated at room temperature. Following washing a HRP conjugated anti-rabbit-IgG antibody is added to each well, incubated at room temperature then again washed. TMB is added and then catalyzed by HRP to produce a blue color product that changes into yellow after adding acidic stop solution. The density of yellow coloration is directly proportional to the Caspase-9 amount of sample captured in plate.			
平台	Microplate			

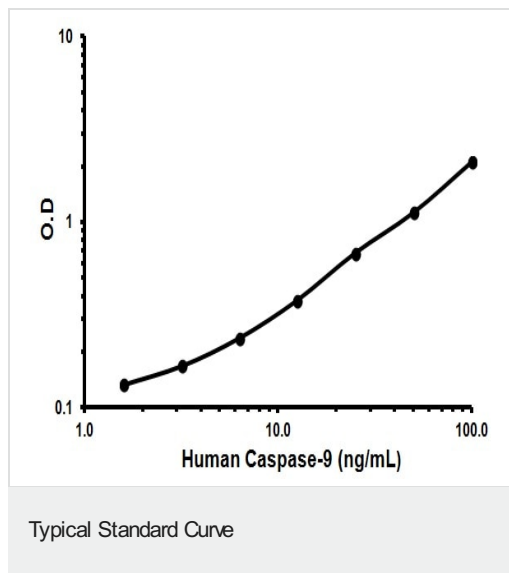
性能

存放说明 Store at +4°C. Please refer to protocols.

组件	1 x 96 tests
20X Assay Buffer Concentrate	1 x 5ml
20X Wash Buffer Concentrate	1 x 50ml
Adhesive Films	4 units
Anti-rabbit-IgG HRP conjugated antibody	1 x 10µl
Human Caspase-9 Standard	2 vials
10X Lysis buffer	1 x 15ml
Microplate coated with monoclonal antibody to Human Caspase-9 (12 x 8 wells)	1 unit
Polyclonal anti-Human Caspase-9 Detection Antibody (rabbit)	1 x 70µl
Sample Diluent	1 x 12ml
Stop Solution (1M Phosphoric acid)	1 x 15ml
TMB Substrate Solution	1 x 15ml

功能	Involved in the activation cascade of caspases responsible for apoptosis execution. Binding of caspase-9 to Apaf-1 leads to activation of the protease which then cleaves and activates caspase-3. Proteolytically cleaves poly(ADP-ribose) polymerase (PARP). Isoform 2 lacks activity is an dominant-negative inhibitor of caspase-9.
组织特异性	Ubiquitous, with highest expression in the heart, moderate expression in liver, skeletal muscle, and pancreas. Low levels in all other tissues. Within the heart, specifically expressed in myocytes.
序列相似性	Belongs to the peptidase C14A family. Contains 1 CARD domain.
发展阶段	Expressed at low levels in fetal heart, at moderate levels in neonate heart, and at high levels in adult heart.
翻译后修饰	Cleavages at Asp-315 by granzyme B and at Asp-330 by caspase-3 generate the two active subunits. Caspase-8 and -10 can also be involved in these processing events.

图片



Representative Standard Curve using ab119508.

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