

Product datasheet

MMP-3 Activity Assay Kit (Fluorometric) ab118972

2 References 2 图像

概述

产品名称	MMP-3 Activity Assay试剂盒(Fluorometric)
检测方法	Fluorescent
样品类型	Cell culture supernatant, Milk, Urine, Tissue Extracts, Cell culture media
检测类型	Enzyme activity
灵敏度	< 0.05 mU/well
检测时间	1h 00m
种属反应性	与反应: Mammals, Other species
产品概述	In Abcam's MMP-3 Activity Assay Kit, MMP-3 hydrolyzes a specific FRET substrate to release the quenched fluorescent group Mca, which can be detected fluorometrically at Ex/Em=325/393 nm. The kit provides a rapid, simple, sensitive and reliable test which can also be used as a high throughput assay of MMP-3 activity. The assay sensitivity is < 50 μ U. Visit our FAQs page for tips and troubleshooting.

说明	The matrix metalloproteinase-3 (MMP-3, stromelysin-1) exhibits a number of activities that would make it a particularly good tumor promoter. Like several other MMPs, MMP-3 was first cloned and later re-cloned as a cancer-specific gene. In addition to degrading numerous extracellular matrix components, MMP-3 can activate gelatinase B, the collagenases and several serpin-type serine proteinase inhibitors. Moreover, it can release a number of cell surface molecules, including E-cadherin, a known contributor to cancer development.
----	--

性能

存放说明 Store at -20°C. Please refer to protocols.

组件	标识符	100 tests
Mca Standard (1mM)	Yellow	1 x 20 μ l
MMP-3 Assay Buffer	WM	1 x 25ml
MMP-3 Positive Control (lyophilized)	Green	1 vial
MMP-3 Substrate	Red	1 x 200 μ l

功能 Can degrade fibronectin, laminin, gelatins of type I, III, IV, and V; collagens III, IV, X, and IX, and

cartilage proteoglycans. Activates procollagenase.

序列相似性

Belongs to the peptidase M10A family.

Contains 4 hemopexin-like domains.

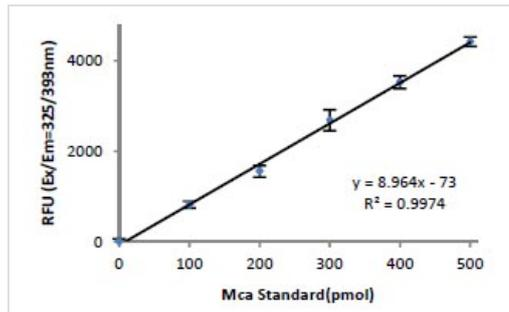
结构域

The conserved cysteine present in the cysteine-switch motif binds the catalytic zinc ion, thus inhibiting the enzyme. The dissociation of the cysteine from the zinc ion upon the activation-peptide release activates the enzyme.

细胞定位

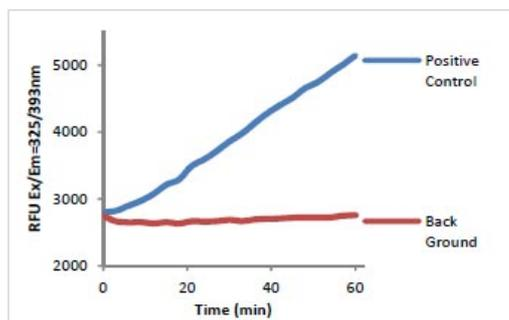
Secreted > extracellular space > extracellular matrix.

图片



MMP-3 Mca Standard

Functional Studies - MMP-3 Activity Assay Kit
(ab118972)



MMP-3 Positive Control

Functional Studies - MMP-3 Activity Assay Kit
(ab118972)

Please note: All products are "FOR RESEARCH USE ONLY AND ARE NOT INTENDED FOR DIAGNOSTIC OR THERAPEUTIC USE"

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