

MMP8 Inhibitor Screening Assay Kit (Colorimetric) ab139452

[1 References](#) [4 图像](#)

概述

产品名称	MMP8抑制剂Screening Assay试剂盒(Colorimetric)
检测方法	Colorimetric
样品类型	Inhibitor compounds
检测类型	Enzyme activity
产品概述	Abcam MMP8 Inhibitor Screening Assay Kit (Colorimetric) (ab139452) is a complete assay system designed to screen MMP8 inhibitors using a thiopeptide as a chromogenic substrate (Ac-PLG-[2-mercapto-4-methyl-pentanoyl]-LG-OC ₂ H ₅). The MMP cleavage site peptide bond is replaced by a thioester bond in the thiopeptide. Hydrolysis of this bond by an MMP produces a sulfhydryl group, which reacts with DTNB [5,5'-dithiobis(2-nitrobenzoic acid), Ellman's reagent] to form 2-nitro-5-thiobenzoic acid, which can be detected by its absorbance at 412 nm ($\epsilon=13,600 \text{ M}^{-1}\text{cm}^{-1}$ at pH 6.0 and above). The assays are performed in a convenient 96-well microplate format.
说明	This kit is useful to screen inhibitors of MMP8, a potential therapeutic target. The MMP inhibitor NNGH is also included as a prototypic control inhibitor. Thiol inhibitors should not be used with this kit, as they may interfere with the colorimetric assay.
平台	Microplate reader

性能

存放说明 Please refer to protocols.

组件	1 x 96 tests
96-well Clear Microplate (1/2 Volume)	1 unit
Colorimetric Assay Buffer	1 x 20ml
MMP Inhibitor	1 x 50 μ l
MMP Substrate	1 x 50 μ l
MMP8 Enzyme (Human, Recombinant)	1 x 22 μ l

功能 Can degrade fibrillar type I, II, and III collagens.

组织特异性

Neutrophils.

序列相似性

Belongs to the peptidase M10A family.

Contains 4 hemopexin-like domains.

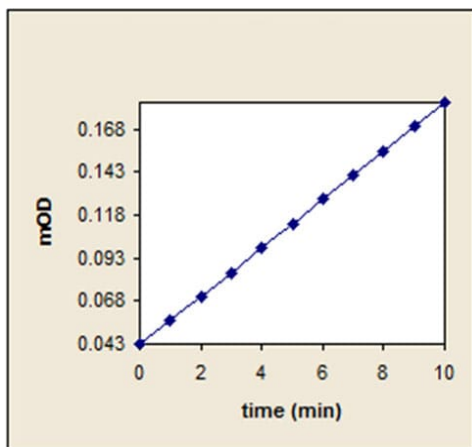
结构域

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

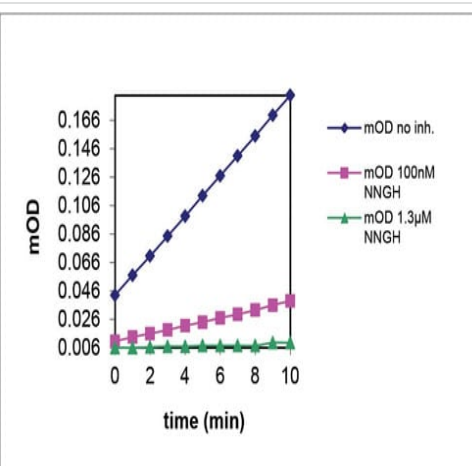
Cytoplasmic granule. Secreted > extracellular space > extracellular matrix. Stored in intracellular granules.

图片



Slope = $V = 1.39E-02$ OD/min

Plot of OD vs. time.

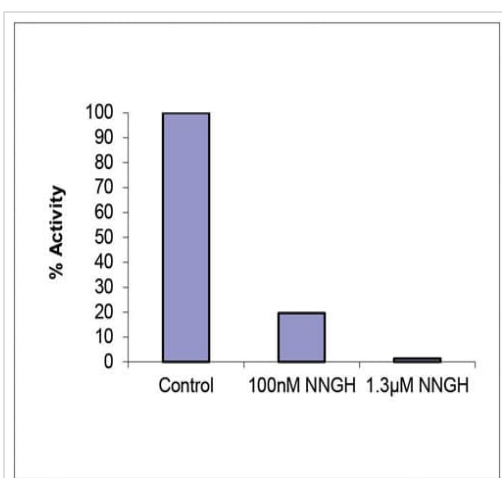


control slope = $1.39E-02$ OD/min

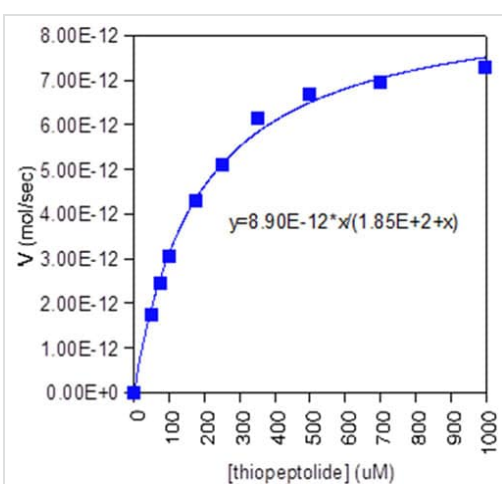
inhibitor slope (100nM) = $2.73E-03$ OD/min

inhibitor % activity remaining = $(2.73E-03/1.39E-02) \times 100 = 19.6\%$

Inhibition of MMP8 by NNGH



Inhibition of MMP8 by NNGH



$K_m = 185 \mu\text{M}$

$V_{max} = 8.90 \text{ pmol/sec}$

Example graph for K_m and V_{max} determination

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