

MMP9 Inhibitor Screening Assay Kit (Fluorometric) ab139449

画像数 1

製品の概要

製品名	MMP9 Inhibitor Screening Assay Kit (Fluorometric)
検出方法	Fluorescent
サンプルの種類	Inhibitor compounds
アッセイタイプ	Enzyme activity
製品の概要	Abcam MMP9 Inhibitor Screening Assay Kit (Fluorometric) (ab139449) is a complete assay system designed to screen MMP9 inhibitors using a quenched fluorogenic peptide: MMP Fluorogenic Substrate Mca-Pro-Leu-Gly-Leu-Dpa-Ala-Arg-NH ₂ [Mca=(7-methoxycoumarin-4-yl)-acetyl; Dpa=N-3-(2,4-dinitrophenyl)-L-α-β-diaminopropionyl]. Mca fluorescence is quenched by the Dpa group until cleavage by MMPs at the Gly-Leu bond separates the two moieties. The assays are performed in a convenient 96-well microplate format.
特記事項	This kit is useful to screen inhibitors of MMP9, a potential therapeutic target. The MMP inhibitor NNGH is also included as a prototypic control inhibitor.
試験プラットフォーム	Microplate reader

製品の特性

保存方法 Please refer to protocols.

内容	1 x 96 tests
96-well White Microplate 1/2 Volume	1 unit
Fluorometric Assay Buffer	1 x 20ml
MMP Calibration Standard	1 x 50μl
MMP Fluorogenic Substrate	1 x 200μl
MMP Inhibitor	1 x 50μl
MMP9 Enzyme (Human, Recombinant)	1 x 48.5μl

機能 May play an essential role in local proteolysis of the extracellular matrix and in leukocyte migration. Could play a role in bone osteoclastic resorption. Cleaves KiSS1 at a Gly-

-Leu bond. Cleaves type IV and type V collagen into large C-terminal three quarter fragments and shorter N-terminal one quarter fragments. Degrades fibronectin but not laminin or Pz-peptide.

組織特異性

Produced by normal alveolar macrophages and granulocytes.

関連疾患

Intervertebral disc disease
Metaphyseal anadysplasia 2

配列類似性

Belongs to the peptidase M10A family.
Contains 3 fibronectin type-II domains.
Contains 4 hemopexin repeats.

ドメイン

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.

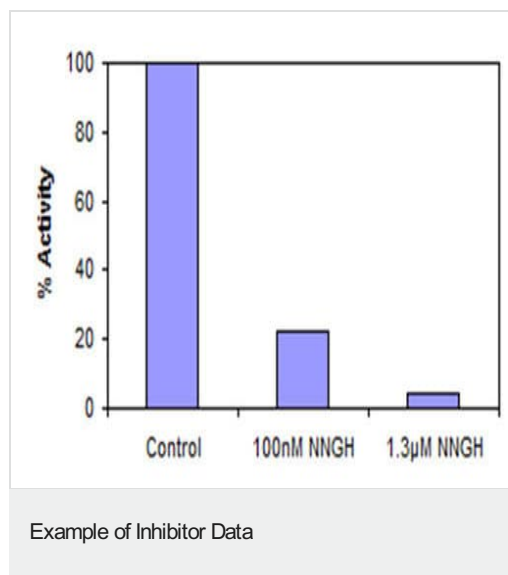
翻訳後修飾

Processing of the precursor yields different active forms of 64, 67 and 82 kDa. Sequentially processing by MMP3 yields the 82 kDa matrix metalloproteinase-9.
N- and O-glycosylated.

細胞内局在

Secreted, extracellular space, extracellular matrix.

画像



Inhibitor of MMP9 by NNGH.

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.co.jp/abpromise> or contact our technical team.

Terms and conditions

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