

MMP7 Inhibitor Screening Assay Kit (Fluorometric)
ab139445

1 Image

Overview

Product name	MMP7 Inhibitor Screening Assay Kit (Fluorometric)
Detection method	Fluorescent
Sample type	Inhibitor compounds
Assay type	Enzyme activity
Product overview	Abcam MMP7 Inhibitor Screening Assay Kit (Fluorometric) (ab139445) is a complete assay system designed to screen MMP7 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.
Notes	This kit is useful to screen inhibitors of MMP7, a potential therapeutic target. The MMP inhibitor NNGH is also included as a prototypic control inhibitor.
Platform	Microplate reader

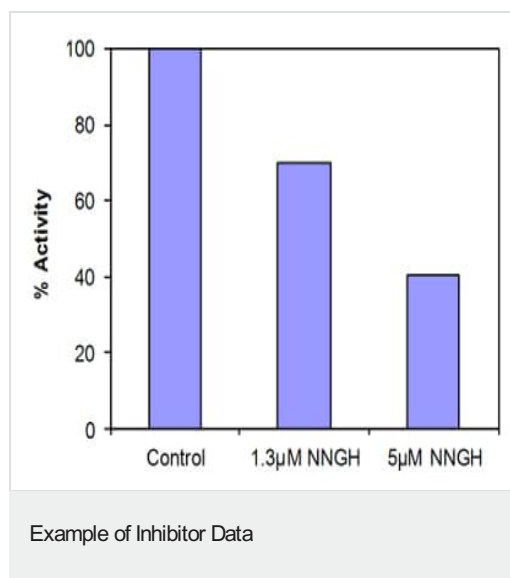
Properties

Storage instructions Please refer to protocols.

Components	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
MMP7 Enzyme (Human, Recombinant)	1 x 33.3µl

Function	Degrades casein, gelatins of types I, III, IV, and V, and fibronectin. Activates procollagenase.
Sequence similarities	Belongs to the peptidase M10A family.
Domain	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.
Cellular localization	Secreted > extracellular space > extracellular matrix.

Images



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.com/abpromise> or contact our technical team.

Terms and conditions

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