

## Product datasheet

# MMP14 Inhibitor Screening Assay Kit (Fluorometric) ab139455

[1 Image](#)

### Overview

<b>Product name</b>	MMP14 Inhibitor Screening Assay Kit (Fluorometric)
<b>Detection method</b>	Fluorescent
<b>Sample type</b>	Inhibitor compounds
<b>Assay type</b>	Enzyme activity
<b>Product overview</b>	Abcam MMP14 Inhibitor Screening Assay Kit (Fluorometric) (ab139455) is a complete assay system designed to screen MMP14 inhibitors using a quenched fluorogenic peptide: MMP Fluorogenic Substrate Mca-Pro-Leu-Gly-Leu-Dpa-Ala-Arg-NH <sub>2</sub> [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.
<b>Notes</b>	This kit is useful to screen inhibitors of MMP14, a potential therapeutic target. The MMP inhibitor NNGH is also included as a prototypic control inhibitor.
<b>Platform</b>	Microplate reader

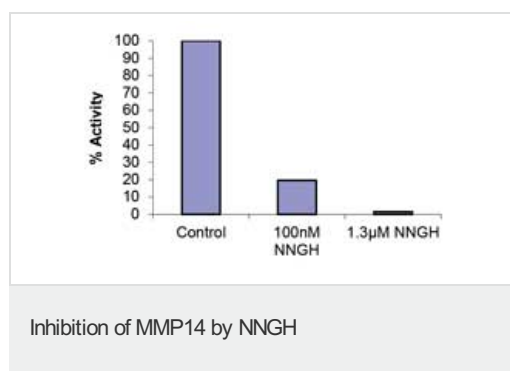
### Properties

**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
MMP14 Enzyme (Human, Recombinant)	1 x 25µl

<b>Function</b>	Seems to specifically activate progelatinase A. May thus trigger invasion by tumor cells by activating progelatinase A on the tumor cell surface. May be involved in actin cytoskeleton reorganization by cleaving PTK7.
<b>Tissue specificity</b>	Expressed in stromal cells of colon, breast, and head and neck. Expressed in lung tumors.
<b>Sequence similarities</b>	Belongs to the peptidase M10A family. Contains 4 hemopexin-like domains.
<b>Domain</b>	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.
<b>Post-translational modifications</b>	The precursor is cleaved by a furin endopeptidase.
<b>Cellular localization</b>	Membrane. Melanosome. Identified by mass spectrometry in melanosome fractions from stage I to stage IV.

## 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