

ab139455

MMP14 Inhibitor

Screening Assay Kit

(Fluorometric)

Instructions for Use

For the screening of MMP14 inhibitors

This product is for research use only and is not intended for diagnostic use.

Table of Contents

1.	Background	3
2.	Principle of the Assay	3
3.	Protocol Summary	4
4.	Materials Supplied	5
5.	Storage and Stability	6
6.	Materials Required, Not Supplied	7
7.	Assay Protocol	8
8.	Data Analysis	12

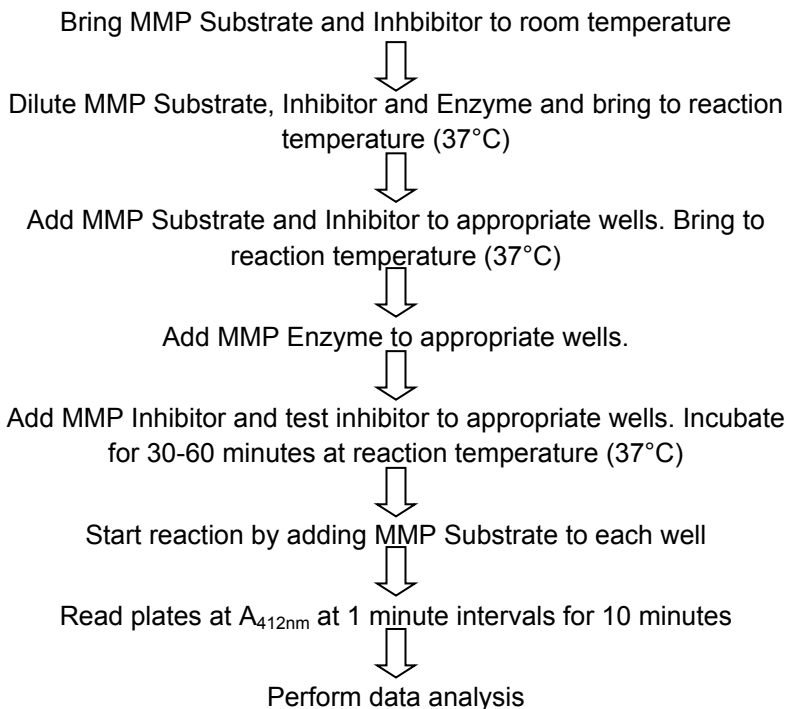
1. Background

Matrix metalloproteinase-14 (MMP14, membrane-type MMP1, MT1-MMP) is a member of the MMP family of extracellular proteases. These enzymes play a role in many normal and disease states by virtue of their broad substrate specificities. Targets of MMP14 include collagen, gelatin, aggrecan, CD44, and pro- α 1-3. MMP14 is secreted as a 63 kDa proenzyme (as measured by SDS PAGE), and activated by cleavage to 60 kDa and below. MMP14 is an important target for inhibitor screening due to its involvement in cancer and ocular pathology.

2. Principle of the Assay

Abcam MMP14 Inhibitor Screening Assay Kit (Fluorometric) (ab139455) is a complete assay system designed to screen MMP14 inhibitors using a quenched fluorogenic peptide: fluorogenic substrate Mca-Pro-Leu-Gly-Leu-Dpa-Ala-Arg-NH₂ [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. The kit is useful to screen inhibitors of MMP14, a potential therapeutic target.

3. Protocol Summary



4. Materials Supplied

Item	Quantity	Storage
96-well White Microplate ½ Volume	1 unit	RT
Fluorometric Assay Buffer	1 x 20 mL	RT
MMP Calibration Standard (40 µM (17.8 µg/mL) in DMSO)	1 x 50 µL	-80°C
MMP Fluorogenic Substrate (400 µM (437 µg/mL) in DMSO)	1 x 200 µL	-80°C
MMP Inhibitor (1.3mM NNGH in DMSO)	1 x 50 µL	-80°C
MMP14 Enzyme (Human, Recombinant) (12 U/µL)	1 x 25 µL	-80°C

5. Storage and Stability

- Store components as stated in table for the highest stability.
- The MMP14 enzyme should be handled carefully in order to retain maximal enzymatic activity. It is stable, in diluted or concentrated form, for several hours on ice.
- As supplied, MMP14 enzyme is stable for 5 freeze/thaw cycles. To minimize the number of freeze/thaw cycles, aliquot the MMP14 into separate tubes and store at -80°C.
- When setting up the assay, do not maintain diluted components at reaction temperature (e.g. 37°C) for an extended period of time prior to running the assay.

6. Materials Required, Not Supplied

- Microplate reader capable of excitation at 328 nm and emission at 420 nm. The following Ex/Em have also been used: 320, 340/393, 400, 405.
- Pipettes or multi-channel pipettes capable of pipetting 1-100 μL accurately.
(Note: reagents can be diluted to increase the minimal pipetting volume to $>10 \mu\text{L}$).
- Ice bucket to keep reagents cold until use.
- Water bath or incubator for component temperature equilibration.

7. Assay Protocol

1. Briefly warm kit components MMP Fluorogenic Substrate, MMP Calibration Standard and MMP Inhibitor to RT to thaw DMSO.
2. Dilute MMP inhibitor 1/200 in Assay Buffer as follows. Add 1 μL inhibitor into 200 μL Assay Buffer, in a separate tube. Warm to reaction temperature (e.g. 37°C).
3. Dilute sufficient volume of MMP Fluorogenic Substrate to 40 μM in assay buffer to required total volume (10 μL are needed per well). Warm to reaction temperature (e.g. 37°C).
4. Dilute MMP14 Enzyme 1/100 in assay buffer to required total volume (20 μL are needed per well). Warm to reaction temperature (e.g. 37°C) shortly before assay.
5. Pipette assay buffer into each desired well of the $\frac{1}{2}$ volume microplate as follows:
 - Calibration = 80 μL in 3 wells (see step 11)
 - Control (no inhibitor) = 70 μL
 - MMP Inhibitor = 50 μL
 - Test inhibitor = varies (see Table 1)
6. Allow microplate to equilibrate to assay temperature (e.g. 37°C).

7. Add 20 μL MMP14 Enzyme (diluted in step 4) to the control, inhibitor, and test inhibitor wells. Final amount of MMP14 will be 2.4 U per well (24.0 mU/ μL). Remember to not add MMP14 to the calibration wells!
8. Add 20 μL MMP inhibitor (diluted in step 2) to the inhibitor wells only! Final inhibitor concentration=1.3 μM .
9. Add desired volume of test inhibitor to appropriate wells. See Table 1.
10. Incubate plate for 30-60 minutes at reaction temperature (e.g. 37°C) to allow inhibitor/enzyme interaction.
11. In the meantime, calibrate the fluorescent microplate reader, using Ex/Em=328/420: Prewarm assay buffer to reaction temperature in 3 wells in the microplate, then to each add 10 μL MMP Fluorogenic Substrate to give the concentration to be used in the assay (e.g., for 4 μM final add 10 μL 40 μM) and mix. When the fluorescent signal is constant, use this reading as the zero (Blank) value in arbitrary fluorescence units (RFUs). Using the same wells, with their mixtures of MMP Substrate and Assay Buffer, add 10 μL MMP Calibration standard to give 3 different final molar concentrations ranging between 2 and 10% of the substrate peptide molar concentration (e.g., 80, 200, and 400 nM) and measure their fluorescence. Use these values to build a standard curve relating micromolar MMP Calibration Standard concentration (x axis) to RFUs (y axis). The slope of the line is the conversion factor (CF). If multiple

concentrations of MMP Substrate peptide are used, such as in kinetic determinations, Section 7, step 11 must be performed for each concentration, due to absorptive quenching by the substrate peptide. Note: this calibration can be done at any time.

12. Start reaction by the addition of 10 μ L MMP14 Fluorogenic Substrate (diluted and equilibrated to reaction temperature in step 3). Final substrate concentration = 4 μ M.
13. Continuously read plates fluorescent microplate reader, using Ex/Em=328/420. For example, record data at 1 minute time intervals for 10 minutes.
14. Perform data analysis (see next section).

NOTE: Retain microplate for future use of unused wells.

Table 1. Example of Samples

Sample	Assay Buffer	MMP14 (120 mU/μL)	Inhibitor (6.5 μM)	Substrate (40 μM)	Total Volume
Control	70 μ L	20 μ L	0 μ L	10 μ L	100 μ L
MMP Inhibitor	50 μ L	20 μ L	20 μ L	10 μ L	100 μ L
Test Inhibitor*	X μ L	20 μ L	Y μ L	10 μ L	100 μ L

*Test inhibitor is the experimental inhibitor. Dissolve/dilute inhibitor into assay buffer and add to appropriate wells at desired volume "Y". Adjust volume "X" to bring the total volume to 100 μ L.

Example of plate:	well#	sample
	A1	Calibration
	B1	Calibration
	C1	Calibration
	D1	Control
	E1	Control
	F1	MMP Inhibitor
	G1	MMP Inhibitor
	H1	Test inhibitor
	A2 ...	Test inhibitor...

8. Data Analysis

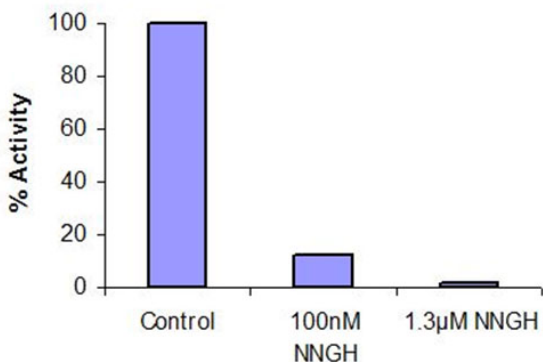
1. Plot data as RFUs (minus Blank RFU value determined during calibration, step 11) versus time for each sample.
2. Determine the range of time points during which the reaction is linear.
3. Obtain the initial reaction velocity (**V**) in RFUs/min: determine the slope of a line fit to the linear portion of the data plot using an appropriate routine.
4. It is best to use a range of inhibitor concentrations, each in duplicate. Average the slopes of duplicate samples.

A. To determine inhibitor % remaining activity:

Inhibitor % activity remaining = $(V_{\text{inhibitor}}/V_{\text{control}}) \times 100$.

See Figure 1 for example of results.

Figure 1. Inhibition of MMP14 by NNGH



B. To find the activity of the samples expressed as picomoles substrate hydrolyzed per minute:

Employ the following equation:

$$X \text{ pmoles substrate/min} = 1/\text{CF} \times V \times \text{vol}$$

Where:

CF is the conversion factor (micromolar concentration/RFUs, from step 11, Assay Protocol)

V is initial reaction velocity in (RFUs/min)

vol. is the reaction volume in microliters (100)

UK, EU and ROW

Email: technical@abcam.com | Tel: +44-(0)1223-696000

Austria

Email: wissenschaftlicherdienst@abcam.com | Tel: 019-288-259

France

Email: supportscientifique@abcam.com | Tel: 01-46-94-62-96

Germany

Email: wissenschaftlicherdienst@abcam.com | Tel: 030-896-779-154

Spain

Email: soportecientifico@abcam.com | Tel: 911-146-554

Switzerland

Email: technical@abcam.com

Tel (Deutsch): 0435-016-424 | Tel (Français): 0615-000-530

US and Latin America

Email: us.technical@abcam.com | Tel: 888-77-ABCAM (22226)

Canada

Email: ca.technical@abcam.com | Tel: 877-749-8807

China and Asia Pacific

Email: hk.technical@abcam.com | Tel: 108008523689 (中國聯通)

Japan

Email: technical@abcam.co.jp | Tel: +81-(0)3-6231-0940

www.abcam.com | www.abcam.cn | www.abcam.co.jp