

MMP Activity Assay Kit (Fluorometric - Red) ab112147

14 References [画像数 4](#)

医薬用外毒物

製品の概要

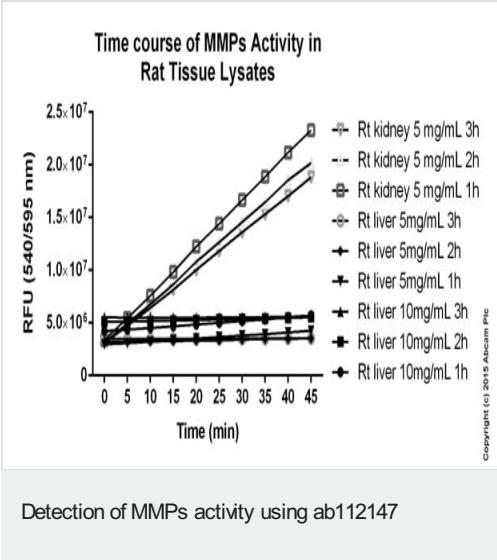
製品名	MMP Activity Assay Kit (Fluorometric - Red)
検出方法	Fluorescent
サンプルの種類	Purified protein, Tissue Lysate
アッセイタイプ	Direct
種交差性	交差種: Mammals, Other species
製品の概要	<p>MMP Activity Assay Kit (Fluorometric Red) ab112147 uses a fluorescence resonance energy transfer (FRET) peptide as a MMP substrate. In the intact FRET peptide, the fluorescence of one part is quenched by the other. Upon cleavage into two separate fragments by MMPs, the fluorescence is recovered.</p> <p>The MMP assay is designed to check the general activity of a MMP enzyme in a tissue sample. It can also be used to screen MMP inhibitors when a purified MMP enzyme is used.</p> <p>With excellent fluorescence quantum yield and longer wavelength, the substrate used in this assay shows less interference from autofluorescence of test compounds and cellular components and is much more sensitive than an EDANS/Dabcyl FRET substrate.</p> <p>The MMP assay signal can be easily read by a fluorescence microplate reader at Ex/Em = 540/590 nm. The pH-independent fluorescence makes the assay reading available for the whole physiological pH range.</p> <p>The high photostability of this FRET peptide provides a useful imaging probe. Many labs have used this kit for the high throughput screening of MMP inhibitors as potential anticancer drug candidates. This assay might be also used for monitoring cancer cells.</p>
特記事項	<p>ab112147 should be stored Desiccated</p> <p>Abcam has not and does not intend to apply for the REACH Authorisation of customers' uses of products that contain European Authorisation list (Annex XIV) substances.</p> <p>It is the responsibility of our customers to check the necessity of application of REACH Authorisation, and any other relevant authorisations, for their intended uses.</p>
試験プラットフォーム	Microplate reader

製品の特性

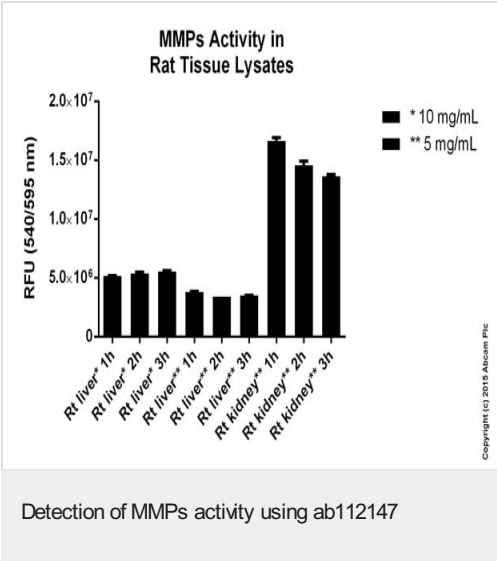
保存方法Store at -20°C. Please refer to protocols.

内容	100 tests
APMA, 4-Aminophenylmercuric Acetate	1 x 20µl
Assay Buffer	1 x 20ml
MMP Red Substrate	1 x 60µl

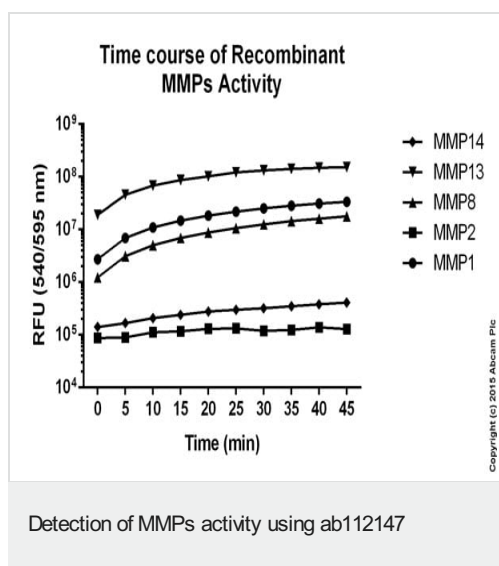
画像



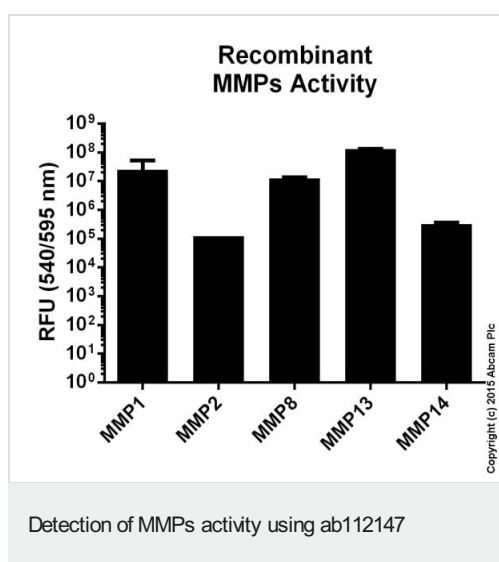
Tissues were lysed with RIPA buffer and activated with 2 mM APMA (1:1) for 1,2 and 3 hours at 37°C. Samples were then diluted to 5 mg/mL and 10 mg/mL with Assay Buffer. 50 µl of MMP containing sample was mixed with MMP Red Substrate. The fluorescence signal was monitored 30 min after the start of the reaction by using a microplate reader with a filter set of Ex/Em = 540/595 nm. The reading from all wells was subtracted with the reading from substrate control, which contains MMP Red Substrate but no MMPs.



Tissues were lysed with RIPA buffer and activated with 2 mM APMA (1:1) for 1,2 and 3 hours at 37°C. Samples were then diluted to 5 mg/mL and 10 mg/mL with Assay Buffer. 50 µl of MMP containing sample was mixed with MMP Red Substrate. The fluorescence signal was monitored 30 min after the start of the reaction by using a microplate reader with a filter set of Ex/Em = 540/595 nm. The reading from all wells was subtracted with the reading from substrate control, which contains MMP Red Substrate but no MMPs.



MMPs were activated with 2 mM APMA (1:1). Samples were then diluted to 0.6 µg/mL (30 ng per well) with Assay Buffer. 50 µl of MMP containing sample was mixed with MMP Red Substrate. The fluorescence signal was monitored 30 min after the start of the reaction by using a microplate reader with a filter set of Ex/Em = 540/595 nm. The reading from all wells was subtracted with the reading from substrate control, which contains MMP Red Substrate but no MMPs.



MMPs were activated with 2 mM APMA (1:1). Samples were then diluted to 0.6 µg/mL (30 ng per well) with Assay Buffer. 50 µl of MMP containing sample was mixed with MMP Red Substrate. The fluorescence signal was monitored 30 min after the start of the reaction by using a microplate reader with a filter set of Ex/Em = 540/595 nm. The reading from all wells was subtracted with the reading from substrate control, which contains MMP Red Substrate but no MMPs.

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