

## Data Sheet

### **Fluorogenic MMP13 Assay Kit**

**Catalog #79991**  
**Size: 96 reactions**

**BACKGROUND:** MMP13 (matrix metalloproteinase 13), also known as collagenase 3, is a member of the matrix metalloproteinase (MMP) family involved in the degradation of the extracellular matrix. MMP13 primarily cleaves type II collagen, and it is expressed in the skeleton where it plays a role in restructuring the collagen matrix for bone mineralization and cartilage turnover. MMP13 is also overexpressed in rheumatoid arthritis, osteoarthritis, and many human carcinomas, suggesting its potential as a therapeutic target.

**DESCRIPTION:** The *Fluorogenic MMP13 Assay Kit* is designed to measure MMP13 activity for screening and profiling applications, in a homogeneous assay with no time-consuming washing steps. The kit comes in a convenient 96-well format, with purified MMP13 enzyme, fluorogenic substrate, and MMP assay buffer for 100 enzyme reactions.

**COMPONENTS:**

Catalog #	Component	Amount	Storage	
11345	MMP13, His-tag	4 µg	-80°C	<b>Avoid freeze/ thaw cycles!</b>
79919	Fluorogenic MMP Substrate (1 mM)	10 µl	-80°C	
79917	1X MMP Assay Buffer 1	25 ml	-20°C	
79685	Black, low binding black microtiter plate	1	Room Temperature	

**APPLICATIONS:** Great for studying enzyme kinetics and HTS applications.

**STABILITY:** At least one year from date of receipt when stored as directed.

**REFERENCE(S):**

1. Wang, M., *et al.* MMP13 is a critical target gene during the progression of osteoarthritis. *Arthritis Res Ther* **15**: R5 (2013).
2. Li, H., *et al.* New insights on the MMP-13 regulatory network in the pathogenesis of early osteoarthritis. *Arthritis Res Ther* **19**: 248 (2017).

**MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:**

Fluorescent microplate reader capable of reading  $\lambda_{exc}/\lambda_{em}=328\text{ nm}/393\text{ nm}$

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401**, Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)

**ASSAY PROTOCOL:**

**All samples and controls should be tested in duplicate.**

- 1) Dilute 1 mM MMP substrate 1:100 in 1X assay buffer, to make a 10  $\mu$ M solution. Dilute only enough as is required for the assay. Store remaining 1 mM substrate in aliquots at -80°C.
- 2) Prepare the substrate solution: N wells  $\times$  (20  $\mu$ l 1X assay buffer + 5  $\mu$ l diluted (10  $\mu$ M) MMP Substrate).
- 3) Add 25  $\mu$ l of the substrate solution to each well (Final concentration of the MMP substrate in a 50  $\mu$ l reaction is 1  $\mu$ M).

Component	Positive Control	Test Sample	Blank
Substrate solution	25 $\mu$ l	25 $\mu$ l	25 $\mu$ l
Test Inhibitor	-	5 $\mu$ l	-
Inhibitor buffer (usually 10% DMSO in water)	5 $\mu$ l	-	5 $\mu$ l
MMP13 (1.6 ng/ $\mu$ l)	20 $\mu$ l	20 $\mu$ l	-
1X Assay Buffer	-	-	20 $\mu$ l
<b>Total</b>	<b>50 <math>\mu</math>l</b>	<b>50 <math>\mu</math>l</b>	<b>50 <math>\mu</math>l</b>

- 4) Prepare the inhibitor solution.

The final concentration of DMSO in the assay should not exceed 1%. If the inhibitor compound is dissolved in DMSO, make a 100-fold higher concentration of the compound than the highest concentration you want to test in DMSO. Then make a 10-fold dilution in 1X assay buffer (at this step the compound concentration is 10-fold higher than the final concentration in 10% DMSO). To determine an IC<sub>50</sub> or to test lower concentrations of the compound, prepare a series of further dilutions in 1X assay buffer containing 10% DMSO (the final concentration of the DMSO will be 1% in all samples).

If the inhibitor compound is dissolved in water, make a solution of the compound 10-fold higher than the final concentration in 1X assay buffer.

- 5) Add 5  $\mu$ l inhibitor solution to each well designated "Test Sample." Add 5  $\mu$ l of inhibitor buffer (same solution without inhibitor; usually 10% DMSO in water ) to the "Blank" and "Positive Control" wells.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

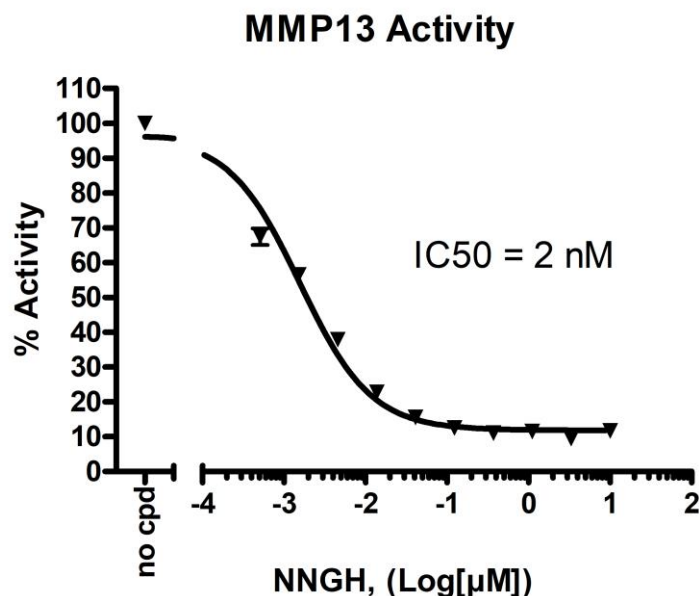
To place your order, please contact us by Phone **1.858.202.1401**, Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)

- 6) Thaw MMP13 on ice. Upon first thaw, briefly spin tube containing enzyme to recover the full content of the tube. Aliquot MMP13 into single use aliquots. Store remaining undiluted enzyme in aliquots at  $-80^{\circ}\text{C}$ . Note: MMP13 enzyme is sensitive to freeze/thaw cycles. Do not re-use diluted enzyme.
- 7) Dilute MMP13 in 1x assay buffer at  $1.6\text{ ng}/\mu\text{l}$  ( $32\text{ ng}$  per reaction).
- 8) Add  $20\ \mu\text{l}$  diluted MMP13 enzyme solution to wells designated as "Positive Control" and "Test Sample." Add  $20\ \mu\text{l}$  1X assay buffer to the "Blank" wells.
- 9) Incubate at room temperature for 30 minutes. Measure the fluorescence intensity in a microtiter plate-reading fluorimeter capable of excitation at a wavelength  $328\text{ nm}$  and detection of emission at a wavelength  $393\text{ nm}$ . The fluorescence intensity can also be measured kinetically. "Blank" value is subtracted from all other values.

#### EXAMPLE OF ASSAY RESULTS:



Inhibition of MMP13 enzyme activity by NNGH, measured using the *Fluorogenic MMP13 Assay Kit* (BPS Bioscience #79991). Fluorescence intensity was measured using a Tecan fluorescent microplate reader. *Data shown is lot-specific. For lot-specific information, please contact BPS Bioscience, Inc. at [info@bpsbioscience.com](mailto:info@bpsbioscience.com)*

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401**, Fax **1.858.481.8694**

Or you can Email us at: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

Please visit our website at: [www.bpsbioscience.com](http://www.bpsbioscience.com)



6042 Cornerstone Court West, Ste. B  
San Diego CA 92121  
Tel: 1.858.202.1401  
Fax: 1.858.481.8694  
Email: [info@bpsbioscience.com](mailto:info@bpsbioscience.com)

## RELATED PRODUCTS

<u>Product</u>	<u>Cat. #</u>	<u>Size</u>
MMP1, His-Tag (Human)	80214	20 µg
MMP2, His-Tag (Human)	80213	20 µg
MMP3(K45E), His-Tag (Human)	11346	100 µg
MMP8, His-Tag (Human)	100552	100 µg
MMP9(Q279R), His-Tag (Human)	80215	20 µg
Fluorogenic MMP1 Assay Kit	79983	96 rxns.
Fluorogenic MMP2 Assay Kit	79918	96 rxns.
Fluorogenic MMP3 (K45E) Assay Kit	79907	384 rxns.
Fluorogenic MMP8 Assay Kit	79929	96 rxns.
Fluorogenic MMP9 (Q279R) Assay Kit	79915	96 rxns.
Fluorogenic MMP10 Assay Kit	79986	96 rxns.

OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401**, Fax **1.858.481.8694**

Or you can Email us at: [\*\*info@bpsbioscience.com\*\*](mailto:info@bpsbioscience.com)

Please visit our website at: [\*\*www.bpsbioscience.com\*\*](http://www.bpsbioscience.com)