

Data Sheet

Fluorogenic MMP9 (Q279R) Assay Kit

Catalog #79915
Size: 96 reactions

BACKGROUND: MMP9 (matrix metalloproteinase 9), also known as Gelatinase B, is a member of the matrix metalloproteinase (MMP) family involved in the degradation of the extracellular matrix. Abnormal expression of MMP9 is observed in many diseases, including cancer and coronary artery disease. The (Q279R) mutation, found within the fibronectin type II domain, is associated with increased risk for in-transit metastasis.

DESCRIPTION: The *Fluorogenic MMP9 (Q279R) Assay Kit* is designed to measure MMP9 (Q279R) 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 MMP9 (Q279R) enzyme, fluorogenic substrate, and MMP9 assay buffer for 100 enzyme reactions.

COMPONENTS:

Catalog #	Component	Amount	Storage	
80215	Recombinant Human MMP9 (Q279R)*	>1 µg	-80°C	Avoid freeze/ thaw cycles!
79916	1 mM MMP Substrate	50 µl	-80°C	
79917	1X MMP Assay Buffer 1	25 ml	-20°C	
79685	Black, low binding black microtiter plate	1	Room Temperature	

*We have given excess for ease of retrieval.

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

Fluorescent microplate reader capable of reading exc/em=328 nm/393 nm

APPLICATIONS: Great for studying enzyme kinetics and HTS applications.

STABILITY: One year from date of receipt when stored as directed.

REFERENCE(S):

1. Wang C., *et al.*, 2020 Feb 12. *J. Cell Physiol.* doi: 10.1002/jcp.29485.
2. Owyong M., *et al.*, 2019 Dec; *Life Sci Alliance* **2(6)**: e201800226.
3. Qu J., *et al.*, 2019; *Am. J. Cancer Res.* **9(7)**: 1415–1428.

ASSAY PROTOCOL:

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

Please visit our website at: www.bpsbioscience.com

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 μ M solution. Dilute only enough as is required for the assay.
- 2) Prepare the substrate solution: N wells \times (20 μ l 1X assay buffer + 5 μ l diluted (10 μ M) MMP Substrate).
- 3) Add 25 μ l of the substrate solution to each well (Final concentration of the MMP substrate in a 50 μ l reaction is 1 μ M).
- 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 IC₅₀ 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 μ l inhibitor solution to each well designated "Test Sample." Add 5 μ l of 10% DMSO in water (inhibitor buffer) to "Blank" and "Positive Control" wells.
- 6) Thaw MMP9 (Q279R) on ice. Upon first thaw, briefly spin tube containing enzyme to recover the full content of the tube. Aliquot MMP9 into single use aliquots. Store remaining undiluted enzyme in aliquots at -80°C. Note: MMP9 (Q279R) enzyme is sensitive to freeze/thaw cycles. Do not re-use diluted enzyme.
- 7) Dilute MMP9 (Q279R) in 1x assay buffer at 0.1 ng/ μ l (2 ng per reaction).
- 8) Add 20 μ l diluted MMP9 (Q279R) enzyme solution to wells designated as "Positive Control" and "Test Sample." Add 20 μ l 1X assay buffer to the "Blank" 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**

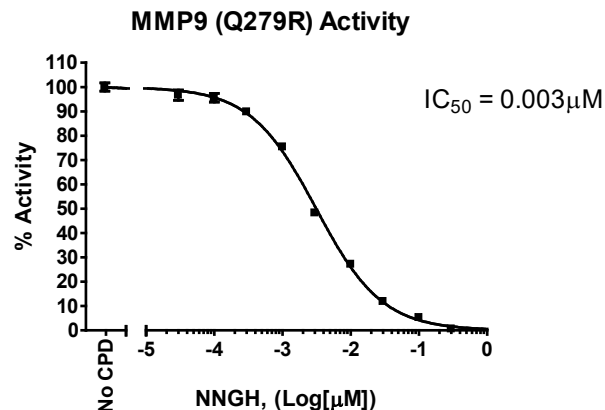
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com

Component	Positive Control	Test Sample	Blank
Substrate solution	25 μ l	25 μ l	25 μ l
Test Inhibitor	-	5 μ l	-
10% DMSO in water (Inhibitor buffer)	5 μ l	-	5 μ l
MMP9 (Q279R, 0.1 ng/ μ l)	20 μ l	20 μ l	-
1X Assay Buffer	-	-	20 μ l
Total	50 μl	50 μl	50 μl

- 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 nm and detection of emission at a wavelength 393 nm. The fluorescence intensity can also be measured kinetically. "Blank" value is subtracted from all other values.

EXAMPLE OF ASSAY RESULTS:



Inhibition of MMP9 (Q279R) enzyme activity by NNGH, measured using the *Fluorogenic MMP9 (Q279R) Assay Kit (BPS Bioscience #79915)*. 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*

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

Please visit our website at: 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

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
MMP3 (K45E) Inhibitor Screening Assay Kit	79907	384 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

Please visit our website at: www.bpsbioscience.com