



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Data Sheet
Fluorogenic MALT1 Assay Kit
Catalog #79809
Size: 96 reactions

BACKGROUND: MALT1 (Mucosa-associated lymphoid tissue lymphoma translocation protein 1) is a cysteine protease that structurally resembles caspases. MALT1 is a member of the CARMA1-BCL10-MALT1 (CBM) complex that mediates NF- κ B activation downstream of the B cell receptor, T cell receptor and ITAM-coupled natural killer cell receptors. MALT1 cleavage activity is essential to T cell and B cell activation and is linked to the pathogenesis of activated B cell-like diffuse large B cell lymphoma (ABC-DLBCL), a chemoresistant form of DLBCL.

DESCRIPTION: The *Fluorogenic MALT1 Assay Kit* is designed to measure MALT1 activity for screening and profiling applications, in a homogeneous assay with no time-consuming washing steps. The MALT1 assay kit comes in a convenient 96-well format, with purified MALT1 enzyme, fluorogenic substrate, and MALT1 assay buffer for 100 enzyme reactions.

COMPONENTS:

Catalog #	Component	Amount	Storage	
100360	Recombinant Human MALT1	15 μ g	-80°C	Avoid freeze/ thaw cycles!
79810	0.5 mM MALT1 Substrate (in DMSO)	100 μ l	-80°C	
79811	1X MALT1 Assay Buffer*	10 ml	-20°C	
	0.5 M DTT	20 μ l	-20°C	
79685	Black, low binding black microtiter plate	1	Room Temperature	

*Add 0.5 M DTT before use

MATERIALS OR INSTRUMENTS REQUIRED BUT NOT SUPPLIED:

Fluorescent microplate reader capable of reading exc/em=360nm/460nm

APPLICATIONS: Great for studying enzyme kinetics and HTS applications.

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

REFERENCE(S):

Fontan, L. *et al.*, *Cancer Cell*. 2012 Dec 11; **22(6)**: 812–824.

Nagel, D. *et al.*, *Cancer Cell*. 2012 Dec 11; **22(6)**: 825-837.

Rebeaud, F. *et al.*, *Nat Immunol*. 2008; **9**: 272-281

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 W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

ASSAY PROTOCOL:

All samples and controls should be tested in duplicate.

- 1) Add 0.5 M DTT to 1X assay buffer so final DTT concentration is 1 mM. For example, add 10 μ l of 0.5M DTT to 5 ml 1X assay buffer. (DTT should be added freshly. Prepare only enough DTT-containing buffer as required for the assay. Store the remaining 1X assay buffer at -20°C).
- 2) Thaw MALT1 on ice. Upon first thaw, briefly spin tube containing enzyme to recover the full content of the tube. Aliquot MALT1 into single use aliquots. Store remaining undiluted enzyme in aliquots at -80°C. Note: MALT1 enzyme is sensitive to freeze/thaw cycles. Do not re-use thawed aliquots or diluted enzyme.
- 3) Dilute MALT1 in 1x assay buffer (with 1 mM DTT) at 7.5 ng/ μ l (150 ng per reaction).
- 4) Add 20 μ l diluted MALT1 enzyme solution to wells designated as "Positive Control" and "Test Sample". Add 20 μ l 1X assay buffer to the "Blank" wells.
- 5) 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.

- 6) Add 5 μ l inhibitor solution to each well designated "Test Sample". Add 5 μ l 1X assay buffer or 10% DMSO (depending on which inhibitor solution is used) to "Blank" and "Positive Control" wells.
- 7) Incubate the enzyme with the compound solution at room temperature for 30 minutes.

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 W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Component	Positive Control	Test Sample	Blank
MALT1 (7.5 ng/μl)	20 μl	20 μl	–
1X Assay Buffer	–	–	20 μl
Test Inhibitor	–	5 μl	–
Inhibitor Buffer (no inhibitor)*	5 μl	–	5 μl
Incubate at room temperature for 30 minutes			
Substrate solution	25 μl	25 μl	25 μl
Total	50 μl	50 μl	50 μl

- 8) Prepare the substrate solution: N wells × (24 μl 1X assay buffer (with DTT) + 1 μl 0.5 mM MALT1 Substrate).
- 9) Add 25 μl of the substrate solution to each well (Final concentration of the MALT1 substrate in a 50 μl reaction is 10 μM).
- 10) Incubate at room temperature for 30 minutes. Measure the fluorescence intensity in a microtiter plate-reading fluorimeter capable of excitation at a wavelength 360 nm and detection of emission at a wavelength 460 nm. The fluorescence intensity can also be measured kinetically. "Blank" value is subtracted from all other values.

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**

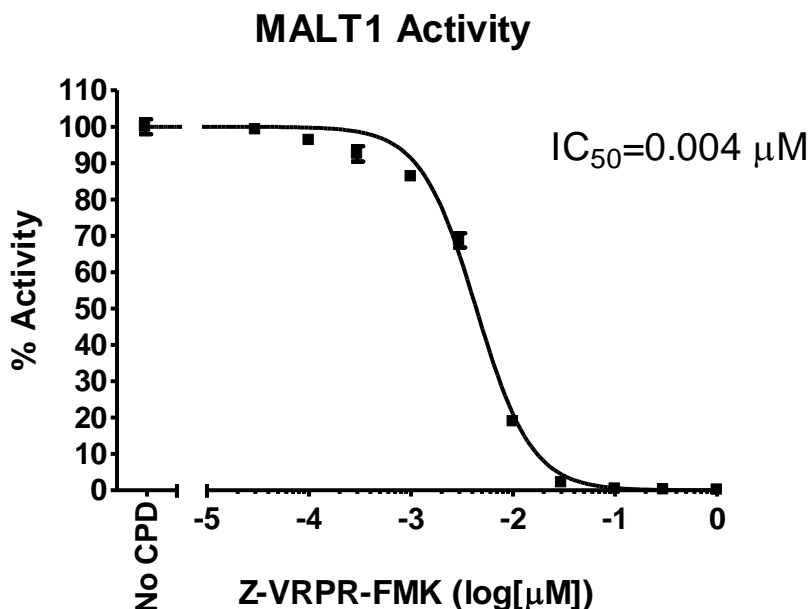
Or you can Email us at: [**info@bpsbioscience.com**](mailto:info@bpsbioscience.com)

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



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.202.1401
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

EXAMPLE OF ASSAY RESULTS:



MALT1 enzyme activity, measured using the *Fluorogenic MALT1 Assay Kit* (BPS Bioscience #79809). 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*

RELATED PRODUCTS

<u>Product</u>	<u>Cat. #</u>	<u>Size</u>
MALT1, FLAG-Tag (Human)	100360	20 µg
MALT1, FLAG-Tag (Mouse)	100054	20 µg
NF-κB reporter (Luc) – NIH/3T3 Cell line	79469	1 vials
A20, His-tag	80393	100 µg

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