

Data Sheet

MALT1, FLAG-Tag (Mouse)

Mouse, Recombinant, N-terminal FLAG-Tag
Catalog #: 100054
Lot #: 180502-2 **Conc.:** 0.13 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0,
 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-
 20, 20% glycerol 100ng/ul flag peptide

Stability: At least 6 months at -80°C . Avoid
 freeze/thaw cycles. Protein may be diluted to
 $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at
 -80°C .

References:

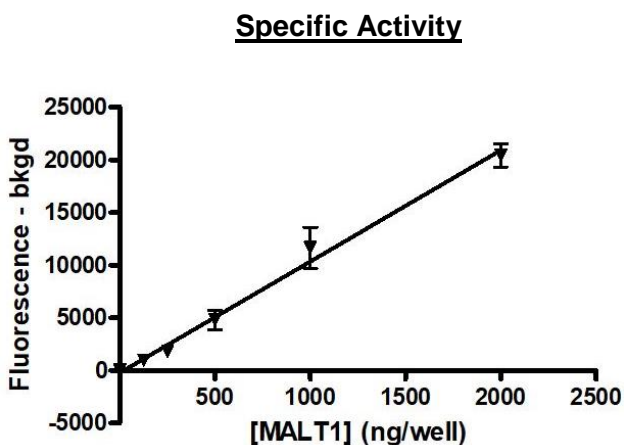
1. Lijun, S., *et al.*, Molecular Cell. 2004. May 14;**3**: 289-301.
2. Lucas, P., *et al.*, JBC. 2001. June 276: 19012-19019.

Description: Mouse MALT1 also known as
 IMD12, MLT, MLT1, PCASP1, and MALT1
 paracaspase, GenBank Accession No.
 NM_172833, a.a. 2-832 (end), and expressed in a
 Baculovirus infected Sf9 cell expression system.
 MW = 93 kDa.

Assay Conditions: The assay was performed by
 fluorescence using a recombinant mMALT1 and
 Ac-LRSR-AMC as substrate. The fluorescence at
 (ex 365 em 460 nm) is correlated with the amount
 of product formed. All of the reactions were
 conducted at 30°C . The reaction mixture is in
 either 96 well black nunc plates or 384 well black
 Greiner plates as indicated. The reaction in
 MALT1 Assay Buffer (50 mM MES, 150 mM NaCl,
 0.1% CHAPS, 1 M Na.citrate, 10 mM DTT added
 fresh, pH 6.8) contains 0-500 μM Ac-LRSR-AMC,
 and 0- 2 μg MALT1. The reaction mixture
 incubated for up to 3h while reading the
 fluorescence signal. For the negative control
 (blank), assay buffer was added instead of the
 enzyme. Absorption signals were measured using
 a Tecan Infinite M1000 plate reader.

Application: Useful as a positive control for
 western blot studies.

Quality Assurance



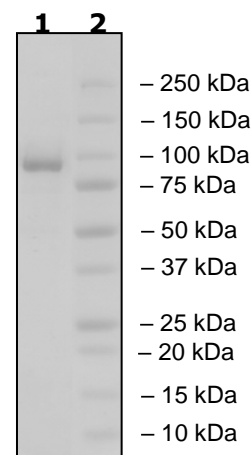
**4-20% SDS-PAGE
 Coomassie staining**

Lane 1:
 MALT1 (3ug)

Lane 2:
 Protein Marker

Purity:
 $>84\%$

MW:
 93 kD



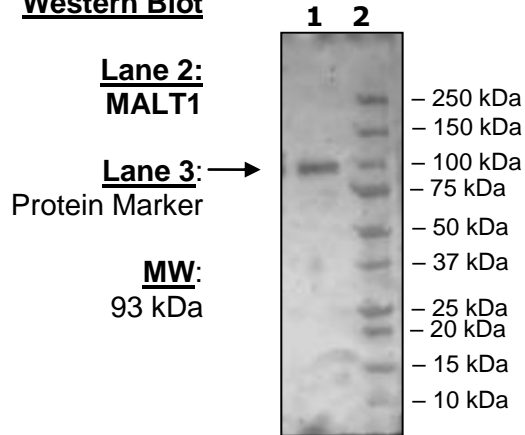
OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.829.3082** 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: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Western Blot



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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com