

## Data Sheet

### MMP3(K45E), His-tag

Human, recombinant, C-terminal His-tag

**Catalog #:** 11346

**Lot #:** 130730

**Conc.:** 1.20 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 200 mM Imidazole, and 20% glycerol.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles.

**References:**

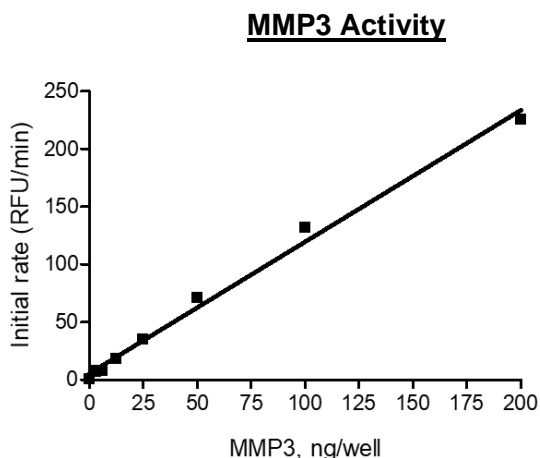
1. Sternlicht, M.D., *et al.*, *Cell*, 1999; **98(2)**:137-146.
2. Buttice, G., *et al.*, *Oncogene*. 1996; **13(11)**:2297-2306.

**Description:** Human matrix metalloproteinase-3 (MMP3), also known as Stromelysin-1 (STMY1), GenBank Accession No. NM\_002422, a.a. 18-477(end) with a Lys to Glu mutation at a.a. 45 and a C-terminal His-tag, expressed in a HEK293 cell expression system. MW=54 kDa.

**Assay Conditions:** Assay buffer: 25 mM Tris pH 9.0, 25  $\mu\text{M}$   $\text{ZnCl}_2$ , 0.05% Tween-20 Enzyme was diluted to 100 ng/ $\mu\text{l}$  in a solution containing 1 mM AMPA (amino-phenyl mercuric acetate) in assay buffer activated for 1 hour at  $37^{\circ}\text{C}$ . Then, 25  $\mu\text{l}$  of 2X 2-fold serial dilution enzyme titration in assay buffer was mixed with 25  $\mu\text{l}$  of 20  $\mu\text{M}$  substrate solution (Mca-RPKPVE-Nval-WRK(Dnp)-NH<sub>2</sub> Fluorogenic MMP Substrate, R&D Systems, ES002, 10 mg/ml = 6 mM stock in DMSO, 10  $\mu\text{M}$  final assay concentration) in a Corning 384-well HV black microplate. Fluorescence was kinetically recorded at 320 nm excitation/405 nm emission for 30 minutes.

**Applications:** Useful for the study of enzyme kinetics, screening inhibitors, and selectivity

### Quality Assurance

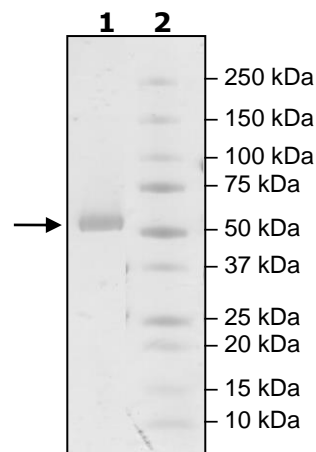


**4-20% SDS-PAGE**  
**Coomassie staining**

**Lane 1:**  
2.4 $\mu\text{g}$  MMP3(K45E)

**Lane 2:**  
Protein Marker

**MW:** 54 kDa  
**Purity:**  $\geq 90\%$



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)