

## Data Sheet

### Cathepsin S, His-Tag

Human, Recombinant, C-terminal His-tag  
**Catalog #:** 80008  
**Lot#:** 170214      **Conc.:** 0.20 mg/ml

**Formulated in:** 45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 225 mM imidazole 3 mM DTT and 10% glycerol % Tween-20 and 10% glycerol.

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

**References:**

1. Seo H.R. *et al.*, *Int. J. Cancer* **124** (8), 1794-1801 (2009).
2. Lutzner N. and Kalbacher H. *J. Biol. Chem.* **283** (52), 36185-36194 (2008).

**Description:**

Human Cathepsin S (GenBank Accession No. NM\_004079.3), Cathepsin S (a.a. 17-331), with C-terminal His-tag, MW=37 kDa, expressed in FreeStyle 293-F cells.

**Specific Activity:**  $\geq 186 \text{ pmole/min}/\mu\text{g}$

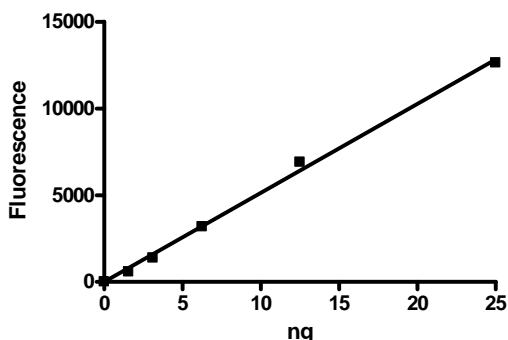
**Assay Condition:** The enzymes were diluted in 25 mM MES buffer pH5, 0.1 M NaCl, 5 mM DTT, and activated by low pH at room temperature for 2 hours. After addition of 10  $\mu\text{M}$  substrate (Z-L-R-AMC) a kinetic reaction was followed for 30 minutes at room temperature and the fluorescence intensity was measured at exc360/em460.

**Application:**

Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Assurance

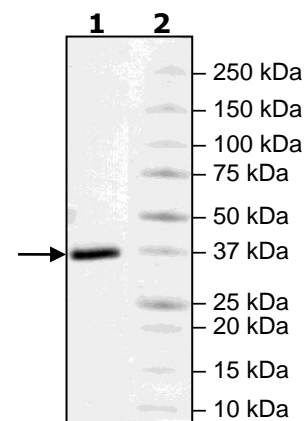
**Specific Activity**



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

**Lane 1:**  
2  $\mu\text{g}$  Cathepsin S  
**Lane 2:**  
Protein Marker

**MW:** 37 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 **1.858.829.3082** 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)