

## Data Sheet

### LIGHT, His-Tag (Mouse)

Mouse, Recombinant, N-terminal His-tag

**Catalog #:** 79068

**Lot #:** 170809      **Conc:** 0.66 mg/ml

**Formulated in:** 8 mM Phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol.

**Stability:** At least 6 months at  $-80^{\circ}\text{C}$ . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

**References:**

1. Garrido, V., *et al.*, *British Journal of Maematology*. 2012; **158(6)**: 788-797.
2. Ogawa, T., *et al.*, *Journal of Urology*. 2010; **183(3)**: 1206-1212.

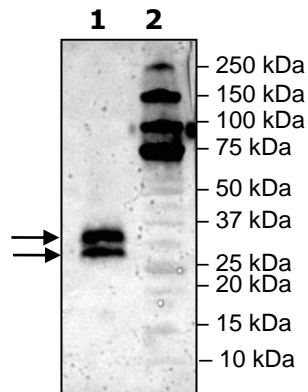
**Description:** Mouse LIGHT protein, also known as TNFSF14, GenBank Accession No. NM\_019418, a.a. 59-239 with N-terminal His-tag. MW = 22 kDa, expressed in a HEK293 cell expression system. This protein runs at a higher M.W. by SDS-PAGE due to glycosylation.

**Application:** Useful as a positive control for western blotting

### Quality Assurance

**Western Blot**

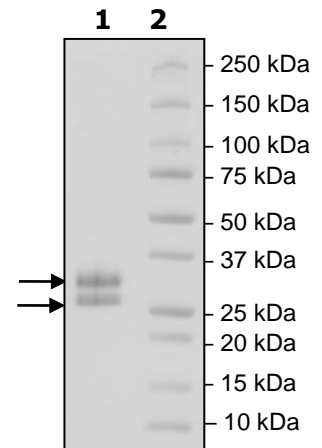
**Lane 1:**  
Light(Mouse)  
**Lane 2:**  
Protein Marker



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

**Lane 1:**  
3µg Light(Mouse)  
**Lane 2:**  
Protein Marker

**MW:** 22 kDa+glycans  
**Purity:** ≥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)