

Data Sheet

LIGHT His-Avi-Tag, Biotin-labeled

Human, Recombinant, N-terminal His-tag,
 C-terminal Avi-tag, Biotin-labeled

Catalog #: 100032

Lot#: 190829 **Conc.:** 0.53 mg/ml

Formulated in: 8 mM phosphate, 110 mM NaCl, 2.2 mM KCl, 20% glycerol.

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles.

References:

1. Edwards, J.R.. *et al.*, *Arthritis & Rheumatology*. 2006; **54(5)**: 1451-1462.
2. Ogawa, T. *et al.*, *The J. Urology*. 2010; **183(3)**: 1206-1212.

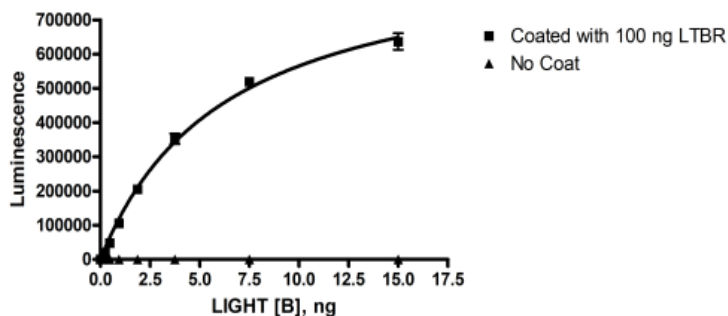
Description: Human LIGHT, also known as TNFSF14, HVEM-L, and LT-y, GenBank Accession No. NM_003807, a.a. 72-240(end) with an N-terminal His-tag and C-terminal Avi-tag, expressed in a HEK293 cell expression system and enzymatically biotin-labeled using Avi-tag™ technology. Biotinylation is confirmed to be $\geq 90\%$. MW=23 kDa. This protein runs at a higher molecular weight due to glycosylation.

Assay Conditions: Reaction done according to LTBR:LIGHT[B] assay kit (BPS Bioscience catalog # 79808) protocol. Buffer is PBS + 0.1% BSA. Blocking buffer is Superblock PBS + 0.05% Tween-20. Coat protein (LTBR) was added overnight at 4 deg 100 ng/well. Various amounts of biotinylated ligand (LIGHT[B]) (0-240 ng) added for 1h at RT. Strep-HRP added for 1h at RT. Luminescence developed with ELISA ECL substrate (BPS Bioscience catalog #79670).

Applications: Useful for studying enzyme kinetics, substrate specificity, and screening inhibitors.

Quality Assurance

LTBR:LIGHT[B] Binding

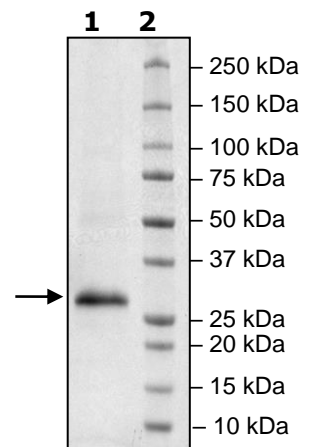


4-20% SDS-PAGE Coomassie staining

Lane 1:
4 μg LIGHT
Lane 2:
Protein Marker

MW: 23 kDa+glycans

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
 Please visit our website at: www.bpsbioscience.com