

Data Sheet

LILRA1, Avi-His Tag, Biotin-labeled

Human, Recombinant, C-terminal Avi-His-Tag,
Biotin-labeled

Catalog #: 100246

Lot #: 190405-1

Conc.: 0.80 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°C . *Avoid
freeze/thaw cycles.*

Description: Human LILRA1, also known as CD85l, LIR-6, and leukocyte immunoglobulin like receptor A1. GenBank Accession No. NM_006863, a.a. 17-461, with a C-terminal Avi-His-Tag, expressed in an HEK293 cell expression system. This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation is confirmed to be $\geq 90\%$. MW = 51 kDa. This protein runs at a higher molecular weight due to glycosylation.

References:

1. Brown, D. *et al.*, *HLA*. 2004; **64(3)**: 215-225.
2. Burshtyn, D. N., *et al.*, *The Journal of Immunology*. 2016; **196(3)**: 947-955.

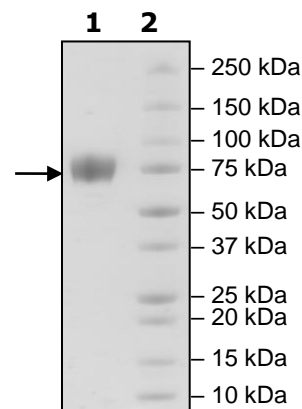
Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:
4 μg LILRA1

Lane 2:
Protein Marker

MW:
51 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