

Data Sheet

LAIR1, Avi-His-Tag

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

Catalog #: 79483

Lot #: 170217-G **Conc.:** 0.43 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. Protein may be diluted to $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at -80°C .

References:

1. Meyaard, L., *et al.*, *Immunity*. 1997; **7(2)**: 283-290.
2. Lebbink, R., *et al.*, *JEM*. 2006; **203(6)**: 1419.

Description: Human LAIR1 also known as Leukocyte-associated immunoglobulin-like receptor 1, CD305, and LAIR-1. GenBank Accession No. NM_002287, a.a. 22-165, with C-terminal AviTM-His-Tag and expressed in a HEK293 cell expression system. MW = 19 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Quality Assurance

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

Lane 1:

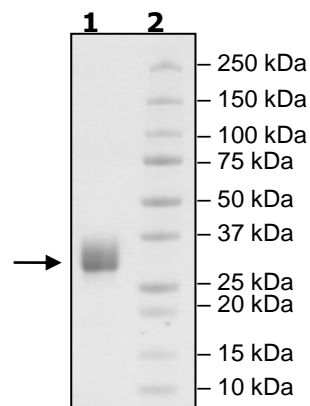
4 μg LAIR1

Lane 2:

Protein Marker

MW: 19kDa+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 **1.858.829.3082** Fax **1.858.481.8694**

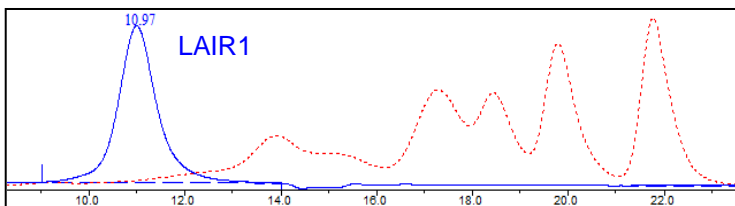
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court W, Ste B
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Gel Filtration Curve



<10% Aggregation

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
Please visit our website at: www.bpsbioscience.com