

Data Sheet

SLAMF7, Avi-His-Tag

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

Catalog #: 100271

Lot#: 190205-3

Conc.: 2.58 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.

References:

1. J. van de Donk, N. W. C. *et al.*, *Blood*. 2016; **127**: 681-695.
2. Huaijian, G. *et al.*, *MCB*. 2015; **35(1)**: 41-51.

Description: Human SLAMF7, also known as CD319, CRACC, CS1, and SLAM family member 7, GenBank Accession No. NM_021181.5, a.a. 23-226 with C-terminal Avi-His-Tag™, expressed in a HEK293 expression system. MW = 25 kDa.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:

4 μg SLAMF7

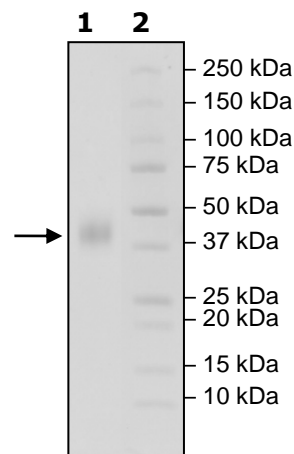
Lane 2:

Protein Marker

MW:

25 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**

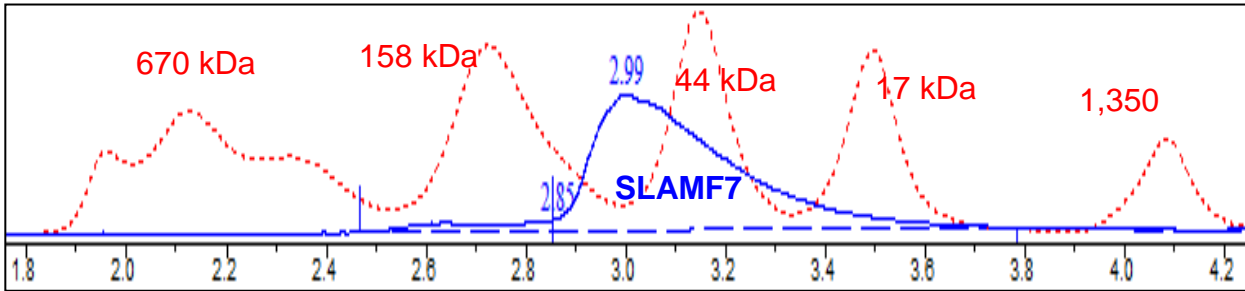
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court West, Suite B
San Diego, CA 92121
Tel: 1.858.202.1401
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 **1.858.202.1401** Fax **1.858.481.8694**
Or you can Email us at: info@bpsbioscience.com
Please visit our website at: www.bpsbioscience.com