

Data Sheet

PDPN, Fc-Fusion, Avi-Tag

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

Catalog #: 79341

Lot#: 171025

Conc.: 0.84 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. Wicki, A., et al., Br. J. Cancer. 2007; 96(1): 1-5.
2. Robertson, N., et al., Genomics. 1995; 23(1):42-50.

Description:

Human PDPN, known as Podoplanin, GenBank Accession No. NM_006474, a.a. 23-131, with C-terminal Avi-Tag fused to the Fc portion of Human IgG1 and expressed in a HEK293 cell expression system. This protein runs at a higher MW by SDS-PAGE due to glycosylation. MW = 40 kDa.

Quality Assurance

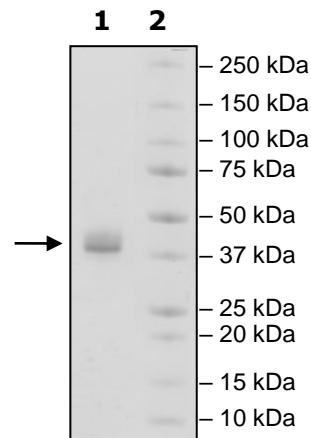
4-20% SDS-PAGE Coomassie staining

Lane 1:
4 μg PDPN

Lane 2:
Protein Marker

MW: 40 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 **1.858.829.3082** Fax **1.858.481.8694**

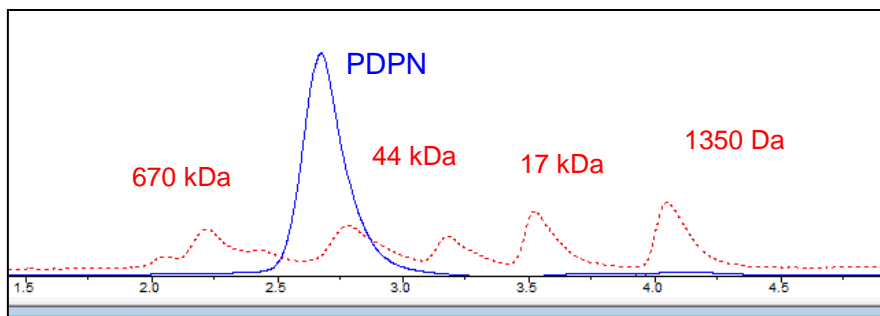
Or you can Email us at: info@bpsbioscience.com

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