

Data Sheet

Mouse CD155, His-tag

Mouse, recombinant, C-terminal, His-tag

Catalog #: 71167

Lot #: 150520

Conc.: 0.54 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. Bottino, C. *et al.*, *J. Exp. Med.* 2003; **198(4)**: 557-567.
2. Pende, D. *et al.*, *Blood*. 2005; **105(5)**: 2066-2073.

Description:

Mouse CD155, also known as Poliovirus receptor and Nectin-like protein 5, GenBank Accession No. NM_027514, a.a. 29-348 with a C-terminal His-tag, expressed in a HEK293 cell expression system. MW = 35.6 kDa. This protein runs at a higher MW by SDS-PAGE due to glycosylation.

Application:

Useful for studying protein binding and for screening small molecules and antibodies.

Quality Assurance

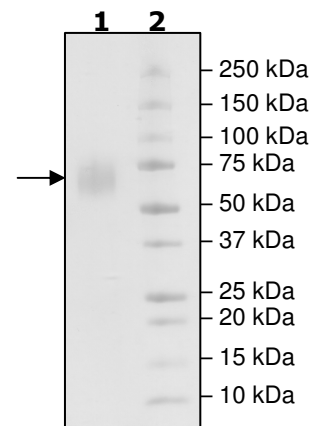
4-20% SDS-PAGE
Coomassie staining

Lane 1:
3 μg CD155

Lane 2:
Protein Marker

MW: 35.6 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**

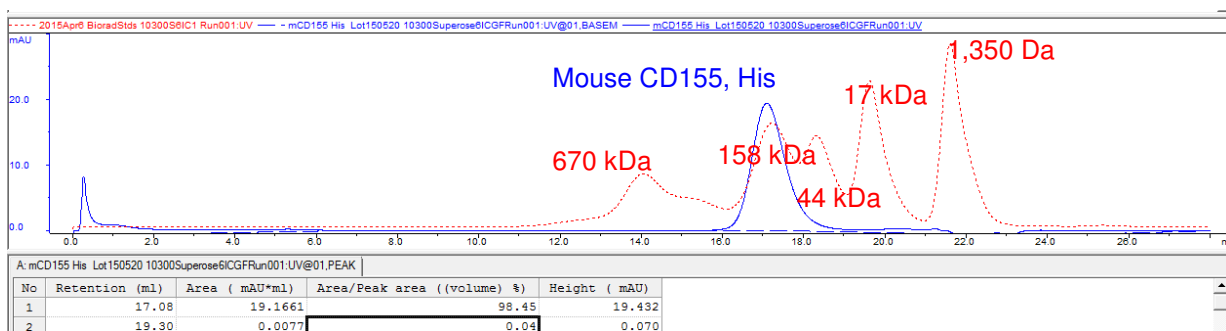
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6044 Cornerstone Court W, Ste E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Gel Filtration



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