

Data Sheet

CD83, Avi-His-Tag

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

Catalog #: 100224

Lot#: 181011-1 **Conc.:** 1.86 mg/ml

Description: Human CD83, also known as cluster of differentiation 83, (GenBank Accession No. NM_004233), a.a. 20-144 with C-terminal Avi-His-Tag, expressed in a HEK293 expression system. MW = 19 kDa.

Formulated in: 8mM phosphate, pH7.4, 110mM NaCl, 2.2mM KCl, and 20% glycerol.

Stability: At least 6 months at -80°C. Avoid freeze/thaw cycles.

References:

1. Zhou, L., *et al.*, *Journal of Immunology*. 1995; **154(8)**: 3821-3835.
2. Lechmann, M. *et al.*, *Trends in Immunology*. 2011; **23(6)**: 273-275.

Quality Assurance

4-20% SDS-PAGE Coomassie staining

Lane 1:

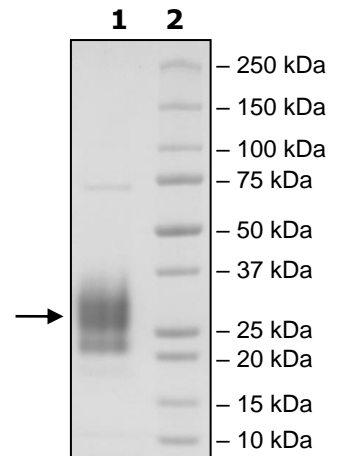
3 µg CD83

Lane 2:

Protein Marker

MW: 19 kDa+glycans

Purity: >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**

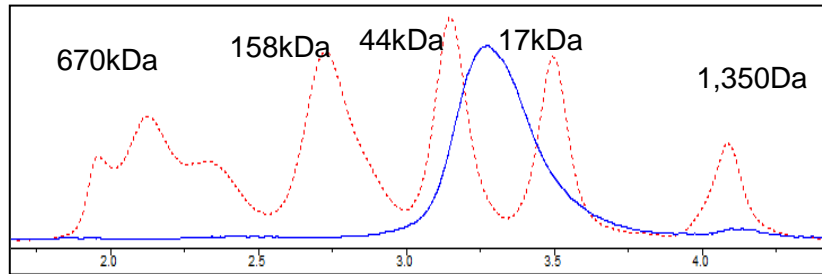
Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com



6042 Cornerstone Court West, Suite 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