

Data Sheet

Human CD46-BC, His-Tag

Human recombinant, C-terminal His-tag

Catalog #: 71230

Lot #: 150519-D

Conc.: 1.0 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. Storing diluted protein is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

1. Liszewski, MK., Atkinson, JP., *J. Immunol.* 1996; **156(11)**:4415-4421.
2. Gerlier, D., *et al.*, *Gen. Virol.* 1994. **75(Pt 9)**:2162-2171.

Description: Human CD46-BC, also known as membrane co-factor protein, Genbank Accession No. BC_030594, a.a. 35–328 with C-terminal His-tag, expressed in a HEK293 cell expression system. MW = 38 kDa (precursor), 34 kDa (mature). Note: this protein is expressed with a 34 amino acid signal sequence that is cleaved after secretion and is heavily glycosylated, resulting in a protein that runs at a higher apparent MW.

Application: Useful for studies of protein function and protein-protein interaction.

Quality Assurance

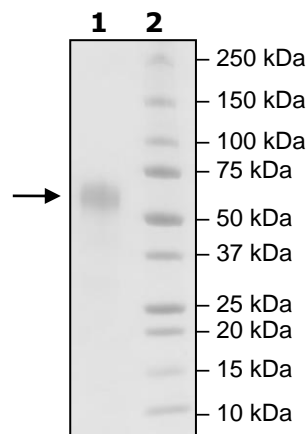
**4-20% SDS-PAGE
Coomassie staining**

Lane 1:
3 μg CD46-BC

Lane 2:
Protein Marker

MW: 34 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 West, Suite E
San Diego, CA 92121
Tel: 1.858.829.3082
Fax: 1.858.481.8694
Email: info@bpsbioscience.com

Sequence

MEPPGRRECPFPSWRFPGLLLAAMVLLLYSFSDACEEPPTFEAMELIGKPKPYEIGER
VDYKCKKGYFYIPPLATHHTICDRNHTWLPVSDDACYRETCPYIRDPLNGQAVPANGTYE
FGYQMHFICNEGYLIGEEILYCELKGSVAIWSGKPPICEKVLCTPPPKIKNGKHTFSEV
EVFEYLDAVTYSCDPAPGPDPSLIGESTIYCGDNSVWSRAAPECKVVKCRFPVVENG
KQISGFGKFFYYKATVMFECDKGFYLDGSDTIVCDSNSTWDPPVPKCLKVSTSSTTKS
PASSASGPRPTYKPPVSNYPGYPKPEEGILDSLDRHHHHHHHHHH

Blue: Signal sequence

Black: CD46-BC (a.a. 35 – 328)

Red: 10x His

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