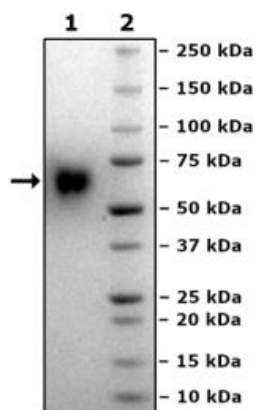


Product Information

Description:	Cynomolgus monkey (<i>Macaca fascicularis</i>) CD46-BC, also known as membrane co-factor protein, encompassing amino acids 35–328. This construct contains a C-terminal His-Tag (10xHis). 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. The recombinant protein was affinity purified. HiP™ indicates a high purity protein (≥90% pure) and less than 10% aggregation as measured by gel filtration.
Species:	Monkey
Construct:	CD46-BC (35-328-His) (Monkey)
Concentration:	1.00 mg/ml
Expression System:	HEK293
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
MW:	34 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Aggregation:	<10%
Genbank Accession:	XM_005540692.2
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



Quality Control Data

Gel Filtration Curve

