

## Product Information

<b>Description:</b>	Recombinant human CD34 (also known as Hematopoietic progenitor cell antigen CD34), encompassing amino acids 32-290. The C-terminus of this construct is fused with the Fc fragment of a human IgG1 followed by an Avi-Tag™. The recombinant protein was affinity purified.
<b>Background:</b>	CD34 (Cluster of Differentiation-34) is a single-pass transmembrane 1phosphor-glycoprotein and a stem cell antigen. It is a receptor for L-Selectin (CD62L) and CrkL ligands, as well as E- and P-Selectin in the bone marrow extracellular matrix. It can act both as a cell adhesion molecule and as an adhesion blocking molecule against mast cells, dendritic precursors and eosinophils. CD34 is often expressed on the surface of hematopoietic stem cells (HSC) and is commonly used for positive selection and identification of CD34 <sup>+</sup> cells for stem cell therapy. More recently CD34 has been identified as a potential predictor of neoadjuvant chemotherapy efficacy in cervical cancer patients. Infusion of CD34 <sup>+</sup> cells has been applied in the treatment of spinal cord injury, liver cirrhosis and vascular disease, in addition to hematological malignancies.
<b>Species:</b>	Human
<b>Construct:</b>	CD34 (32-290-Fc(IgG1)-Avi)
<b>Concentration:</b>	1.10 mg/ml
<b>Expression System:</b>	HEK293
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
<b>MW:</b>	56 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Genbank Accession:</b>	NM_001773
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	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

