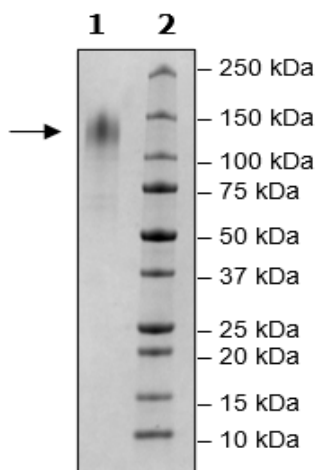


Product Information

| | |
|------------------------------|--|
| Description: | Recombinant human CD34 (CD34 molecule), encompassing amino acids 32-290. This construct contains a C-terminal Avi-Tag™ followed by a His-Tag (6xHis). The recombinant protein was enzymatically biotinylated using Avi-Tag™ and affinity purified. |
| Species | Human |
| Construct: | CD34 (32-290-Avi-His)-(Biotin) |
| Concentration: | 1.02 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: | 31 kDa + glycans |
| Glycosylation: | This protein runs at a higher MW by SDS-PAGE due to glycosylation. |
| Genbank Accession: | NM_001773 |
| Label: | This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation confirmed to be ≥90%. |
| 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



Biotin-Avidin Pulldown

