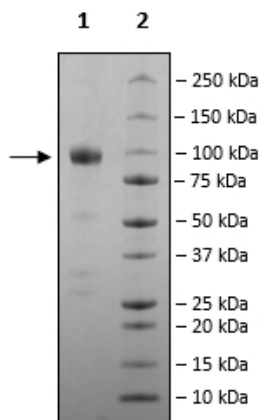


Product Information

Description:	Recombinant human CD66b, encompassing amino acids 35-320. This construct contains a C-terminal human IgG1 Fc domain followed by an Avi-Tag™. This protein was affinity purified.
Background:	CD66b (cluster of differentiation 66b), also known as CEACAM8 (Carcinoembryonic antigen-related cell adhesion molecule 8), belongs to the CEA family of proteins and the immunoglobulin superfamily, and it is expressed in neutrophils, eosinophils, granulocytes, and monocytes. It was initially described as a granulocyte marker, but recent studies have shown that it is also involved in cell adhesion and is a pro-inflammatory mediator. It is highly glycosylated and binds to GPI (glycosylphosphatidylinositol). CD66b ⁺ monocytes can be found in several cancer types and represent a population of cells that do not seem to be involved in immunosuppression but display high phagocytic activity and co-stimulate T cell proliferation and IFN- γ secretion. CD66b ⁺ neutrophils are also present in the tumor microenvironment and it is believed these TIN (tumor infiltrating neutrophils) link to a poor prognosis and can be used to better classify patients for treatment. Anti-CD66b antibodies have been used to generate ADCs (antibody drug conjugates).
Species:	Human
Construct:	CD66b (35-320-Fc(IgG1)-Avi)
Concentration:	1.64 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:	60 kDa + glycans
Glycosylation:	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
Genbank Accession:	NM_001816.4
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.
Assay Conditions:	The protein was validated in ELISA. CD66b (#102029) was coated onto a 96-well plate overnight at 4°C (50 μ l/well at a concentration of 2 μ g/ml in PBS). The plate was washed 3 times with Immuno Buffer 1 (#79311) and blocked with 100 μ l/well of Blocking Buffer 2 (#79728) for 1 hour at Room Temperature (RT). After removing the blocking buffer, 50 μ l/well of serially diluted Anti-CD66b Antibody, Biotin-Labeled (#102078), in Blocking Buffer 2, was added for 30 minutes at RT. The plate was washed 3 times and incubated with Streptavidin-HRP, followed by washing, and incubation with Colorimetric HRP substrate. The reaction was stopped, and absorbance was read at λ =450 nm. The Blank value was subtracted from all values.
Applications:	Useful for binding studies using ELISA.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



CD66:Anti CD66b-Biotin Binding Assay

