

Product Information

Description:	Recombinant human CD137L, also known as TNF superfamily 9 (TNFSF9), encompassing amino acids 71-254(end). This construct contains an N-terminal His-tag (6xHis). The recombinant protein was affinity purified.
Species	Human
Construct:	CD137L (His-71-254(end))
Concentration:	0.17 mg/ml
Expression System:	<i>E. coli</i>
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
MW:	21 kDa
Genbank Accession:	NM_003811
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

