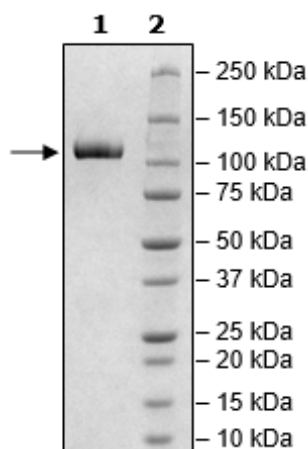


Product Information

Description:	Recombinant human ERAP1 (tumor necrosis factor receptor shedding aminopeptidase regulator), full length, encompassing amino acids 1-941(end). This construct contains a C-terminal Avi-tag™ followed by a His-Tag (10xHis). The recombinant protein was enzymatically biotinylated using the Avi-Tag™ and affinity purified.
Label:	This protein is enzymatically biotinylated using Avi-Tag™ technology. Biotinylation confirmed to be ≥90%.
Species	Human
Construct:	ERAP1 (1-941(end)-Avi-His)-(Biotin)
Concentration:	0.31 mg/ml
Expression System:	HEK293
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	50 mM HEPES, pH 7.4, 500 mM NaCl, 250 mM Imidazole, and 20% glycerol
MW:	110 kDa
Genbank Accession:	EAW96077
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

