

## Product Information

<b>Description:</b>	Recombinant human TGIT (T-cell immunoreceptor with Ig and ITIM domain), encompassing amino acids 22-141. This construct contains a C-terminal Fc region. The recombinant protein was affinity purified.
<b>Species</b>	Human
<b>Construct:</b>	TIGIT (22-141-Fc(IgG1))
<b>Concentration:</b>	0.92 mg/ml
<b>Expression System:</b>	HEK293
<b>Purity:</b>	88%
<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>	40 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Genbank Accession:</b>	NM_173799
<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

