

## Product Information

<b>Description:</b>	Recombinant human CD277 (cluster of differentiation 277), encompassing amino acids 31-254. This construct contains a C-terminal Avi-Tag™, followed by an His-Tag (6xHis). This protein was affinity purified.
<b>Background:</b>	CD277 (cluster of differentiation 277), also known as BTN3A (butyrophilin subfamily 3), belongs to the BTN family of transmembrane receptors. CD277 is expressed in T and B cells, monocytes, dendritic and NK cells. It can bind to pAGs (phosphate antigen), stimulating Vγ9Vδ2 T cells, a T cell type mainly found in the peripheral blood which can target cancer cells. It is also found in many cancer types and can be used as a prognosis marker in ovarian, breast and bladder cancer, and pancreatic ductal adenocarcinoma. Its expression levels vary with the type of cancer, with breast cancer showing lower levels, while advanced ovarian cancer can present high protein expression. A deeper understanding of the role of CD277 in each cancer type may open new targeted therapeutic opportunities.
<b>Species:</b>	Human
<b>Construct:</b>	CD277 (31-254-Avi-His)
<b>Concentration:</b>	0.77 mg/ml
<b>Expression System:</b>	HEK293
<b>Purity:</b>	≥90%
<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>	27 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Genbank Accession:</b>	NM_007048.6
<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.
<b>Applications:</b>	Useful for SDS-PAGE.

## Quality Control Data

### 4-20% SDS-PAGE Coomassie Staining

