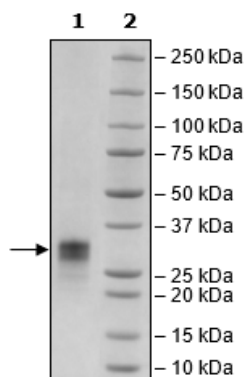


## Product Information

<b>Construct:</b>	OX40L (His-51-183)
<b>Concentration:</b>	1.83 mg/ml
<b>Species:</b>	Human
<b>Formulated In:</b>	8 mM phosphate, pH 7.4, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
<b>Expression System:</b>	HEK293
<b>Format:</b>	Aqueous buffer solution
<b>Stability:</b>	At least 6 months at -80°C. Avoid freeze/thaw cycles.
<b>Storage:</b>	-80°C
<b>Genbank Accession:</b>	NM_003326
<b>MW:</b>	18 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Purity:</b>	≥90%
<b>Assay Conditions:</b>	Assay was performed according to OX40[Biotinylated]:OX40L Inhibitor Screening Assay Kit (BPS Bioscience #72045) with OX40 [Biotin] enzyme titrated at various concentrations.
<b>Applications:</b>	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

## Quality Control Data

## 4-20% SDS-Page Coomassie Staining



## OX40L-OX40[Biotin] Activity

