

Product Information

<b>Construct:</b>	IL-12B(23-328)_IL-12A(23-219-Fc(IgG1)-Avi)
<b>Concentration:</b>	0.84 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>	IL12A(p35): NM_000882; IL12B(p40): NM_002187
<b>MW:</b>	87 kDa + glycans
<b>Glycosylation:</b>	This protein runs at a higher MW by SDS-PAGE due to glycosylation.
<b>Purity:</b>	≥90%

Quality Control Data

4-20% SDS-Page Coomassie Staining

