

## Product Information

<b>Description:</b>	Human secreted interleukin 17 A (IL-17A) also known as cytotoxic T lymphocyte antigen 8 (CTLA8), encompassing amino acids 20-155(end). The construct contains a C-terminal Avi-Tag™ and the recombinant protein was purified by ion exchange chromatography. This protein exists naturally as a homodimer.
<b>Species:</b>	Human
<b>Construct:</b>	IL-17A (20-155(end)-Avi)
<b>Concentration:</b>	0.36 mg/ml
<b>Expression System:</b>	<i>E. coli</i>
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	10 mM Na-phosphate, pH 6.0, and 220 mM NaCl
<b>MW:</b>	Monomer: 18 kDa; Dimer: 35 kDa
<b>Endotoxin Level:</b>	13.9 EU/mg
<b>Genbank Accession:</b>	NM_002190
<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

