

## Product Information

<b>Description:</b>	Recombinant TARC (CCL17) is a monomeric protein consisting of 71 amino acid residues and migrates as an approximately 8 kDa protein under non-reducing conditions and reducing conditions in SDS-PAGE. Optimized DNA sequence encoding Human TARC (CCL17) mature chain was expressed in <i>E. coli</i> .
<b>Species:</b>	Human
<b>Expression System:</b>	<i>E. coli</i>
<b>Purity:</b>	>98% by SDS-PAGE and HPLC
<b>Formulated In:</b>	Lyophilized from a 0.2 µm filtered 20 mM phosphate buffer, 100 mM NaCl solution, pH 7.5.
<b>Endotoxin Level:</b>	<0.1 ng/µg (1 EU/µg), using the LAL gel clot method.
<b>MW:</b>	8 kDa
<b>Storage:</b>	The lyophilized protein is stable for at least years from date of receipt at -20° C. Upon reconstitution, this cytokine can be stored in working aliquots at +4°C for one month, or at -20° C for six months, with a carrier protein without detectable loss of activity. Avoid repeated freeze/thaw cycles.
<b>Reconstitution:</b>	A quick spin of the vial followed by reconstitution in distilled water to a concentration not less than 0.1 mg/ml. This solution can then be diluted into other buffers.
<b>Biological Activity</b>	Determined by its ability to chemoattract human T cells using a concentration range of 2.0-40.0 ng/ml.