

## Product Information

<b>Description:</b>	Recombinant IGFBP-3 is a monomer protein consisting of 269 amino acid residues, migrates as an approximately 28 kDa protein under non-reducing and reducing conditions in SDS-PAGE. Optimized DNA sequence encoding Human IGFBP-3 mature chain fused to an N-terminal His tag was expressed in CHO Cells.
<b>Species:</b>	Human
<b>Expression System:</b>	CHO
<b>Tag:</b>	N-terminal His-Tag
<b>Purity:</b>	>95% by SDS-PAGE and HPLC
<b>Formulated In:</b>	Lyophilized from a 0.2 µm filtered PBS solution.
<b>Endotoxin Level:</b>	<0.1 ng/µg (1 EU/µg), using the LAL gel clot method.
<b>MW:</b>	28 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>	The ED(50) was determined by the ability to inhibit IGF-II induced proliferation of MCF-7 and was found to be ≤ 0.2 µg/ml in the presence of 15 ng/ml human IGF-II.