

Product Information

Description:	Recombinant human TYK2 (tyrosine kinase 2), encompassing amino acids 442-1187(end). This construct contains an N-terminal GST-tag followed by a Thrombin Cleavage site to facilitate the removal of the GST. This recombinant protein was affinity purified.
Species:	Human
Construct:	TYK2 (GST-442-1187(end))
Concentration:	0.10 mg/ml
Expression System:	Sf9
Purity:	86%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, and 20% glycerol
MW:	111 kDa
Genbank Accession:	NM_003331
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	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

