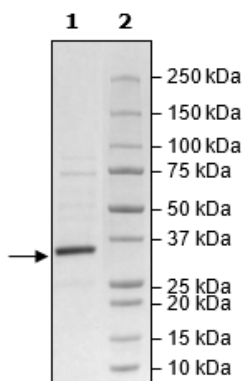


Product Information

Construct:	TYK2 (JH2 domain), (FLAG-575-869)
Concentration:	2.34 mg/ml
Species:	Human
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, and 3 mM DTT
Expression System:	Sf9
Format:	Aqueous buffer solution
Stability:	At least 6 months at -80°C. Avoid freeze/thaw cycles.
Storage:	-80°C
Genbank Accession:	NM_003331
MW:	34 kDa
Purity:	84%
Assay Conditions:	Assay was done according to TYK2 JH2 Pseudokinase Domain Inhibitor Screening Assay Kit (BPS Bioscience #78107) with various amounts of TYK2 (JH2 Domain).
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



TYK2 (JH2 Domain) Activity

