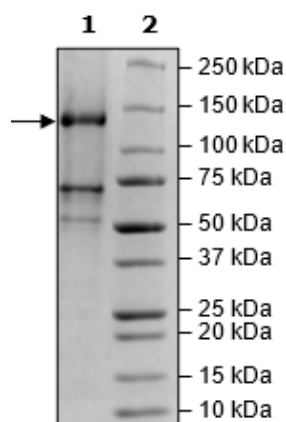


Product Information

Description:	Recombinant human ULK1, also known as ATG1, UNC51, and hATG1 , encompassing amino acids 2–1050 (end). This construct contains an N-terminal FLAG-tag. This protein was affinity purified.
Species:	Human
Construct:	ULK1 (FLAG-2-1050(end))
Concentration:	0.09 mg/ml
Expression System:	HEK293
Purity:	43%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, 20% glycerol, 3 mM DTT, and 80 µg/ml FLAG peptide
MW:	114 kDa
Genbank Accession:	NM_003565
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.
Specific Activity:	≥85 pmol/min/µg
Assay Conditions:	Assay was done according to ULK1 Kinase Assay Kit (BPS Bioscience #78362) with various amounts of ULK1.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



ULK1 Activity

