

Product Information

Description:	Recombinant human Skd3 (also known as ClpB, caseinolytic peptidase B protein homolog), encompassing amino acids 92-707(end). This construct contains an MBP (maltose binding protein) tag followed by a TEV protease site, and an E455Q mutation. This protein was affinity purified.
Background:	Skd3 is a mitochondrial AAA ATPase protein that acts as a chaperone. It uses ATP hydrolysis to chaperone protein unfolding and disassembly, and to reverse protein aggregation. Sdk3 also plays a role in apoptosis by solubilizing HAX1 (HCLS-associated protein X-1). Mutations lead to metabolic disorders characterized by neutropenia, 3-methylglutaconic aciduria and developmental delays. Interestingly, Skd3 has been linked to Venetoclax resistance in the treatment of acute myeloid leukemia (AML), revealing its importance in AML therapy.
Species:	Human
Construct:	SKD3 (E455Q) (MBP-TEV-92-707(end))
Mutation:	E455Q
Concentration:	3.09 mg/ml
Expression System:	<i>E. coli</i>
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	50 mM Tris, pH 8.0, 300 mM KCl, 10% glycerol, 5 mM ATP, 10 mM MgCl ₂ , 3 mM DTT, and 10 mM Maltose
MW:	111 kDa
Genbank Accession:	NM_030813.6
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

