

## Product Information

<b>Description:</b>	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 corresponds to the wild-type protein. This protein was affinity purified.
<b>Background:</b>	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.
<b>Species:</b>	Human
<b>Construct:</b>	Skd3 (MBP-TEV-92-707(end))
<b>Concentration:</b>	0.62 mg/ml
<b>Expression System:</b>	<i>E. coli</i>
<b>Purity:</b>	≥90%
<b>Format:</b>	Aqueous buffer solution.
<b>Formulated In:</b>	50 mM Tris, pH 8.0, 300 mM KCl, 10% glycerol, 5 mM ATP, 10 mM MgCl <sub>2</sub> , 3 mM DTT, and 10 mM Maltose
<b>MW:</b>	111 kDa
<b>Genbank Accession:</b>	NM_030813.6
<b>Stability:</b>	At least 6 months at -80°C.
<b>Storage:</b>	-80°C
<b>Instructions for Use:</b>	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

