

Product Information

Description:	Recombinant human KEAP1 (kelch like ECH associated protein 1), full length, encompassing amino acids 2-624. This construct contains an N-terminal His-tag (6xHis). This recombinant protein was affinity purified.
Species:	Human
Construct:	KEAP1 (His-2-624(end))
Concentration:	0.25 mg/ml
Expression System:	<i>E. coli</i>
Purity:	80%
Format:	Aqueous buffer solution
Formulated In:	40 mM Tris-HCl pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol
MW:	70 kDa
Genbank Accession:	NM_203500
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

