

## Rabbit Anti-Human/Rat Kelch-like 1 Protein

### Instruction Manual

<b>Catalog Number</b>	PK-AB718-3015
<b>Synonyms</b>	KLHL1 Antibody: Kelch-like 1 protein
<b>Description</b>	The mammalian Kelch-like 1 (KLHL1) was initially discovered as a homolog to the <i>Drosophila</i> Kelch gene that is highly expressed in several brain tissues. The predicted protein domain structure of KLHL1 is characteristic of a number of proteins that bind actin, form dimers, and often act as actin-organizing proteins. Based on the presence of anti-sense RNA that spans the transcription and translation start sites as well as the first splice site of KLHL1 in brain tissue of individuals suffering from the neurodegenerative disorder spinocerebellar ataxia type 8 (SCA8), it has been suggested that KLHL1 is involved in this disease and that regulation of KLHL1 protein may be affected by antisense RNA expression.
<b>Quantity</b>	100 µg
<b>Source / Host</b>	Rabbit
<b>Immunogen</b>	Rabbit polyclonal KLHL1 antibody was raised against a 14 amino acid peptide from near the amino-terminus of human KLHL1 (Genbank accession No. AAF81719).
<b>Purification Method</b>	Immunoaffinity chromatography purified IgG
<b>Clone / IgG Subtype</b>	Polyclonal antibody
<b>Species Reactivity</b>	Human, Rat
<b>Specificity</b>	
<b>Formulation</b>	Antibody is supplied in PBS containing 0.02% sodium azide.
<b>Reconstitution</b>	During shipment, small volumes of antibody will occasionally become entrapped in the seal of the product vial. For products with volumes of 200 µl or less, we recommend gently tapping the vial on a hard surface or briefly centrifuging the vial in a tabletop centrifuge to dislodge any liquid in the container's cap.
<b>Storage &amp; Stability</b>	Antibody can be stored at 4°C for three months and at -20°C for up to one year. As with all antibodies care should be taken to avoid repeated freeze thaw cycles. Antibodies should not be exposed to prolonged high temperatures.
<b>Applications</b>	E, WB, IHC INote: Antibody might be suitable for other applications not tested so far. Optimal concentrations for each application have to be determined individually. Application have to be determined individually. KLHL1 antibody can be used for the detection of KLHL1 by Western blot at 2 - 4 µg/mL. Antibody can also be used for immunohistochemistry starting at 10 µg/mL.
<b>Images</b>	Available upon request.
<b>References</b>	Nemes JP, Benzo KA, and Koob MD. The SCA8 transcript is an antisense RNA to a brain-specific transcript encoding a novel actin-binding protein (KLHL1). <i>Human Mol. Gen.</i> 2000; 9:1543-51. Adams J, Kelso R, and Cooley L. The kelch repeat superfamily of proteins: propellers of cell function. <i>Trends Cell Biol.</i> 2000; 10:17-24. Robinson DN and Cooley L. <i>Drosophila</i> kelch is an oligomeric ring canal actin organizer. <i>J. Cell Biol.</i> 1997; 138:799-810.
<b>Images</b>	Available upon request.
<b>Related Products</b>	Blocking Peptide, Cat. No. PK-AB718-3013P; EL4 Cell Lysate, Cat. No. PK-AB718-1287 CARD8 Antibody, Cat. No. PK-AB718-3187; CARD9 Antibody, Cat. No. PK-AB718-2515 CARMA1 Antibody, Cat. No. PK-AB718-3189

FOR IN VITRO RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC PROCEDURES.