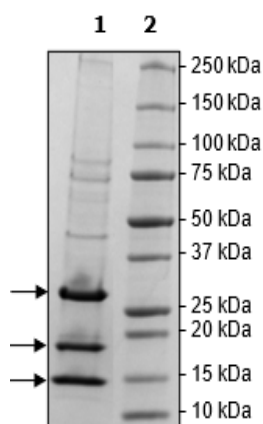


Product Information

Description:	Complex of ELOB (Elongin-B), ELOC (Elongin-C) and VHL (von Hippel-Lindau disease tumor suppressor), expressed individually and affinity purified as a complex. Recombinant human ELOB, full length encompassing amino acids 2-118 (end). This construct contains an N-terminal FLAG-tag. Recombinant human ELOC, full length encompassing amino acids 2-112 (end). This construct contains an N-terminal His-tag (6xHis). Recombinant human VHL, full length encompassing amino acids 2-213 (end). This construct contains an N-terminal FLAG-tag.
Species:	Human
Construct:	ELOB (FLAG-2-118(end)) / ELOC (His-2-112(end)) / VHL (FLAG-2-213(end))
Concentration:	0.14 mg/ml
Expression System:	HEK293
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, and 3 mM DTT
MW:	VHL: 25 kDa; ELOB: 14 kDa; ELOC: 13 kDa
Genbank Accession:	VHL: NM_000551; ELOB: NM_007108; ELOC: NM_005648
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.
Assay Conditions:	Assay run according to VHL Binding Assay Kit (BPS Bioscience #78805) with ELOB/ELOC/VHL Complex titrated at various concentrations. This assay uses fluorescence polarization to determine the binding of the complex to a fluorescent probe.
Applications:	Binding studies and screening of binding inhibitors.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



VHL Binding to probe

