

VHL/ELOB/ELOC Complex

Cat. #UB332

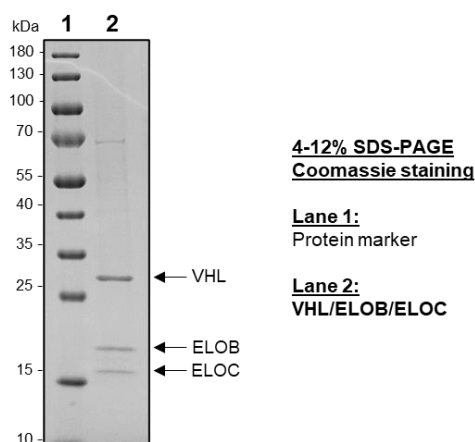
Description: The VHL complex is a multi-subunit ubiquitin ligase constituted with VHL, Elongin B, Elongin C, Culin-2, and Rbx1. Elongin B (ELOB) and Elongin C (ELOC) are heterodimers that bind to the BC-box motif present in SOCS- & VHL-box protein families¹. VHL is the substrate recognition component that brings target specificity. In this specific E3 ligase complex, the ELOB/ELOC works as an adapter between target proteins and Cullin-2/Rbx1 in VHL-box E3 ubiquitin ligases².

Application:

- Protein degradation
- PROTAC and Molecular glue discovery
- Selectivity Profiling

Product Information

Purity	≥ 90% by SDS-PAGE
Molecular Weight	VHL, 25 kDa; ELOB, 14 kDa; ELOC, 13 kDa
Genbank Accession	VHL, NM_000551; ELOB, NM_007108; ELOC, NM_005648
Physical State	Liquid
Tags	N-terminal FLAG-tag on ELOB/VHL, N-terminal His-tag (6xHis) on ELOC
Quantity	10 µg, 50 µg
Buffer	40 mM Tris-HCL, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 2 mM DTT
Storage	-80°C. Avoid repeated freeze/thaw cycles



References

1. Stebbins, C.E., *et al.*, *Science*. 1999; **284(5413)**: 455-461.71-4779.
2. Kamura, T., *et al.*, *Science*. 1999; **284(5414)**: 657-661.

All products are for research use only • not intended for human or animal diagnostic or therapeutic uses
Copyright © 2009 LifeSensors, Inc. All Rights Reserved