

Background

The VHL complex is a multi-subunit E3 ubiquitin ligase that mediates substrate-specific ubiquitination and proteasomal degradation. The VHL complex consists of VHL, Elongin B (ELOB), Elongin C (ELOC), Cullin-2 (CUL2), and RBX1. ELOB and ELOC form a heterodimer that binds to the BC-box motif present in SOCS- and VHL-box protein families. 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 ligases.

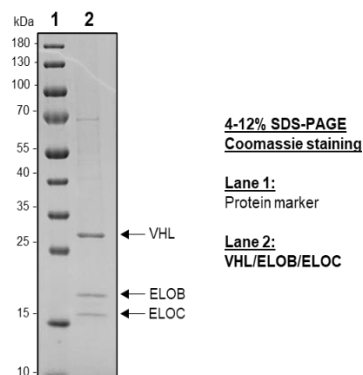
Application(s)

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

Product Specifications

Tag	N-terminal FLAG-tag on ELOB/VHL, N-terminal His-tag (6xHis) on ELOC
Purity	≥ 90% by SDS-PAGE
Molecular Weight	VHL, 25 kDa; ELOB, 14 kDa; ELOC, 13 kDa
Quantity	10 µg, 50 µg
Species	Human. Genbank Accession No: VHL, NM_000551; ELOB, NM_007108; ELOC, NM_005648
Expression System	HEK293
Physical State	Liquid
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

Product QC



References

1. Meyers M, et al., *ACS Chem Biol*. 2024;19(1):58-68.
2. Barankiewicz J, et al., *Cancers (Basel)*. 2022;14(18):4492.
3. Gang, Lu., et al., *Science*. 2014; 343(6168): 305-309.
4. Zhu, Y.X., et al., *Blood*. 2011; 118: 4771-4779.