

## VHL/CUL2/ELOB/ELOC/RBX1 Complex

Cat. # UB331

### Background

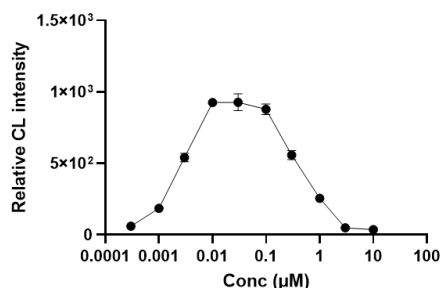
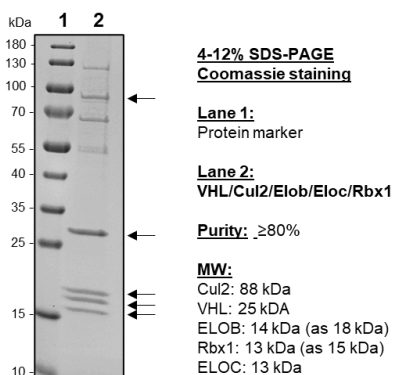
The VHL complex is a multi-subunit ubiquitin ligase constituted with VHL, Elongin B (ELOB), Elongin C (ELOC), Culin-2, and Rbx1. ELOB and ELOC as a heterodimer bind to the BC-box motif present in SOCS- & VHL-box protein families<sup>1</sup>. VHL is the substrate recognition component that brings target specificity. VHL is linked to scaffolding proteins Cullin-2/Rbx1 of the complex via adaptor protein ELOC with ELOB stabilizing the complex<sup>2</sup>.

### Applications

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

### Product Information

<b>Purity</b>	≥ 80% by SDS-PAGE
<b>Molecular Weight</b>	VHL, 25 kDa; CUL2, 88 kDa; ELOB, 14 kDa; ELOC, 13 kDa; Rbx1, 13 kDa
<b>Genbank Accession</b>	VHL, NM_000551; Cul2, NM_003591; ELOB, NM_007108; Rbx1, NM_014248; ELOC, NM_005648
<b>Physical State</b>	Liquid
<b>Quantity</b>	10 µg, 50 µg
<b>Buffer</b>	40 mM Tris-HCL, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 2 mM DTT
<b>Storage</b>	-80°C. Avoid repeated freeze/thaw cycles



**In vitro ubiquitination assay to test the activity of the VHL complex.** In vitro ubiquitination reaction was performed in the presence of various doses of LC2, a VHL degrader of KRAS G12C. Ubiquitinated KRAS G12C was captured on the microtiter plate coated with TUBEs and detected using anti-KRAS antibody. Chemiluminescence intensities were plotted against PROTAC doses to evaluate the extent of ubiquitination.

### 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