

VHL/ELOB/ELOC Complex

Cat. #UB332

Description:

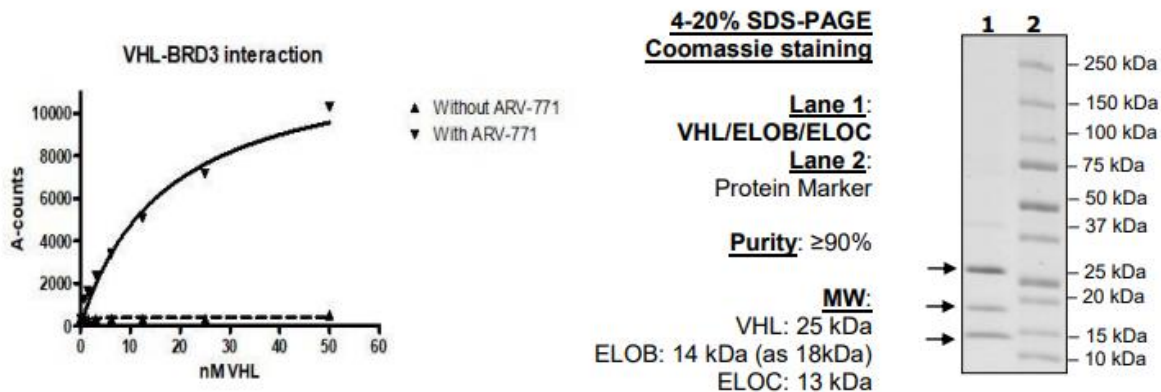
VHL/ELOB/ELOC complex of human ELOB (GenBank Accession No. NM_007108), a.a. 2-118 with N-terminal FLAG-tag and MW= 14 kDa; ELOC (GenBank Accession No. NM_005648), a.a. 2-112(end) with N-terminal His-tag and MW= 13; human VHL (GenBank Accession No. NM_000551) a.a. 2-212(end) with N-terminal FLAG-tag and MW= 25 kDa, individually expressed in a HEK293 expression system.

Application:

- Protein degradation
- PROTAC design
- Selectivity Profiling

Product Information

Purity:	≥ 95% by SDS-PAGE
Molecular Weight:	VHL: 25 kDa, ELOB: 14 kDa, ELOC: 13 kDa
Physical State:	Liquid
Quantity:	10 µg, 50 µg
Buffer:	40 mM Tris-HCL, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 3 mM DTT
Storage:	-80° C. Avoid repeated freeze/thaw cycles



References

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

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