

Data Sheet

VHL/CUL2/ELOB/ELOC/RBX1 Complex

Human, Recombinant

Catalog #: 100373

Lot #: 190503

Conc.: 0.31 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 20% glycerol, 3 mM DTT.

Stability: At least 6 months at -80°C .
 Avoid freeze/thaw cycles.

References:

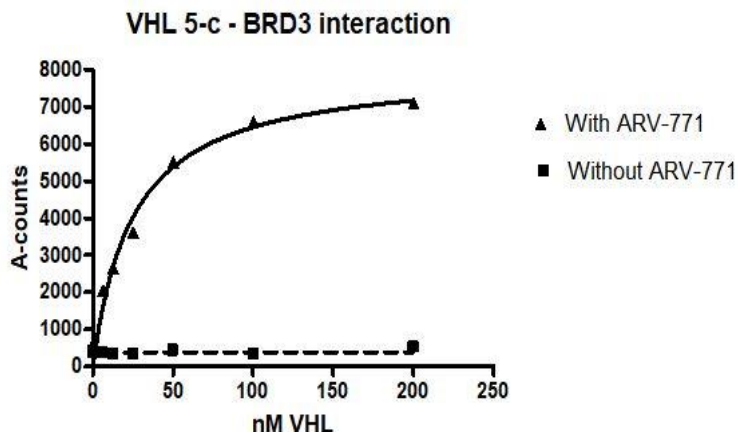
1. Kamura, T., *et al.*, *Science*. 1999; **284(5414)**: 657-661.
2. Stebbins, C.E., *et al.*, *Science*. 1999; **284(5413)**: 455-461.

Description: VHL/CUL2/ELOB/ELOC/RBX1 complex of human **VHL** (GenBank Accession No. NM_000551) a.a. 2-212(end) with N-terminal FLAG-tag and MW= 25 kDa, human **CUL2** (GenBank Accession No. NM_003591), a.a. 2-745(end) with N-terminal FLAG-tag and MW= 88 kDa, 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 kDa human **Rbx1** (GenBank Accession No. NM_014248), a.a. 2-108(end) with N-terminal His-tag and MW= 13 kDa, individually expressed in a HEK293 expression system.

Assay Conditions: ARV-771-mediated binding of VHL BPS #100373 to BET bromodomain was assessed using A-screen technology. Various amounts of VHL were incubated in 10- μl reaction with ARV-771 and GST-BRD3(BD2) BPS #31033 at room temperature for one hour. After this, FLAG-acceptor beads and GSH-donor beads were added followed by detection of A-counts.

Applications: Useful for the study of protein degradation and PROTAC design, and selectivity profiling.

Quality Assurance



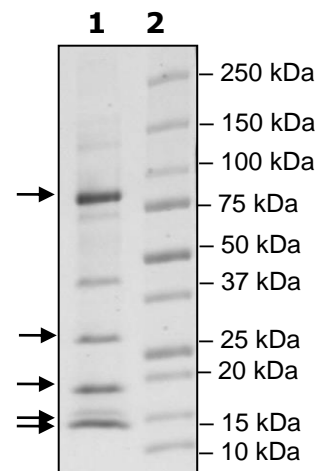
4-20% SDS-PAGE Coomassie staining

Lane 1:
VHL/Cul2/Elob/Eloc/Rbx1

Lane 2:
Protein Marker

Purity: $\geq 80\%$

MW:
 CUL2: 88 kDa
 VHL: 25 kDa
 ELOB: 14 kDa (as 18 kDa)
 Rbx1: 13 kDa (as 15 kDa)
 ELOC: 13 kDa



OUR PRODUCTS ARE FOR RESEARCH USE ONLY. NOT FOR DIAGNOSTIC OR THERAPEUTIC USE.

To place your order, please contact us by Phone **1.858.202.1401** Fax **1.858.481.8694**

Or you can Email us at: info@bpsbioscience.com

Please visit our website at: www.bpsbioscience.com