

Product Information

Description:	Complex of Rbx1 (Ring-Box 1), Cul4B (cullin 4B), Dcaf11 (Ddb1- and Cul4-associated factor 11) and Ddb1 (Damage specific DNA binding protein 1. Recombinant Rbx1, encompassing amino acids 2-108(end), mutated at M5I, contains an N-terminal His-tag (6xHis). Recombinant Cul4B, encompassing amino acids 3-913(end), contains an N-terminal His-tag (6xHis). Recombinant Dcaf11, encompassing amino acids 2-546(end), contains an N-terminal FLAG-tag. Recombinant human Ddb1, encompassing amino acids 2-1140(end) contains an N-terminal FLAG-tag. These recombinant proteins were co-expressed, and affinity purified.
Background:	Eukaryotic cells have evolved complex mechanisms to regulate cellular pathways and protein quality. Cullin proteins are a family of hydrophobic proteins that serve as scaffold and can complex with RING and adaptor proteins to form CRLs (E3 ligase-Cullin-RING ligase) complexes, which can interact with different substrates and regulate multiple cellular processes. When the CRL Rbx1-Cul4b-Ddb1 binds to Dcaf11 it can target p21 for ubiquitination and regulate cell cycle progression.
Species:	Human
Construct:	Rbx1 (M5I) (His-2-108(end)) / Cul4B (His-3-913(end)) / Dcaf11 (FLAG-2-546(end)) / Ddb1 (FLAG-2-1140(end))
Mutation:	Rbx1: M5I
Concentration:	0.55 mg/ml
Expression System:	HEK293
Purity:	89%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 0.04% Tween-20, and 20% glycerol
MW:	Rbx1: 13 kDa; Cul4B: 105 kDa; Dcaf11: 63 kDa; Ddb1: 128 kDa
Genbank Accession:	Rbx1: NM_014248; Cul4B: NM_003588.4; Dcaf11: NM_025230.5; Ddb1: NM_001923
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining

