

## Data Sheet

### EZH2 (E740K) /EED/SUZ12/RbAp48/AEBP2

Human, recombinant  
**Catalog #:** 51032  
**Lot #:** 120514      **Conc.:** 0.38 mg/ml

**Formulated in:** 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 200 mM imidazole, and 20% glycerol.

**Stability:** >6 months at -70°C. Avoid freeze/thaw cycles.

**References:**

1. Varambally, S. *et al. Science* **322 (5908):** 1695-1699 (2008).
2. Rakotobe, D. *et al., Virol. J.* **5 : 32** (2008).

**Application Reference:**

Morin, R.D., *et al. Nat Genet.* 2010 February; **42(2):** 181-185.

**Description:** Mutant version of BPS's regular EZH2 5-member complex (Cat. # 51004), but with a Glu-to-Lys mutation on a.a. 745 of the EZH2 protein.\* Complex of human EZH2 (GenBank Accession No. NM\_004456), (a.a. 2-end ; E745K) with N-terminal His-tag, MW= 86 kDa, human EED (NM\_003797) (a.a. 2-end) with N-terminal FLAG-tag, MW= 51 kDa, human SUZ12 (NM\_015355) (a.a. 2-end) with N-terminal His-tag, MW = 87 kDa, Human AEBP2 (NM\_153207) (a.a. 2-end) with N-terminal His-tag, MW= 53 kDa, and human RbAp48 (NM\_005610) (a.a. 2-end) with N-terminal His-tag, MW = 48 kDa, co-expressed in a baculovirus expression system. \*This mutation is equivalent to E740K in variant 3 (GenBank Accession No. NM\_001203247).

**Application:** Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

### **Quality Assurance**

**4-20% gradient SDS-PAGE**  
**Coomassie staining**

**Lane 1:**  
 6.4 µg enzyme complex

**Lane 2:**  
 Protein Marker

**Purity:** ≥81%



