

## Data Sheet

### **EZH1/EED/SUZ12** **/RbAp48/AEBP2**

Human, recombinant  
**Catalog #:** 51007  
**Lot#:** 130201-EN Conc.: 0.3 mg/ml

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

**Stability:** >6 months at -80°C

**References:**

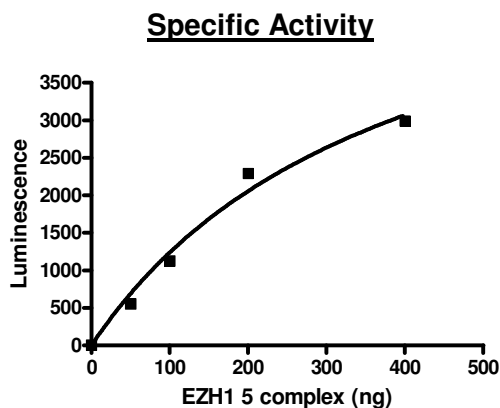
1. Margueron R. *et al. Mol. Cell* **32** (4), 503-518 (2008).
2. Shen X. *et al. Mol. Cell* **32** (4), 491-502 (2008).

**Description:** Complex of human EZH1 (GenBank Accession No. NM\_001991), (a.a. 2-end) 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 baculovirus expression system.

**Assay conditions:** 50 µl reaction mix (20 mM phosphate pH 7.4, 0.05% Tween-20, 40 µM S-adenosylmethionine, and enzyme complex) add to the wells coated with the substrate on a GBioscience Neutravidin white plate. Incubate for 1 hr. Add antibody against methylated K27 residue of histone H3, incubate 1 hr. Then, add secondary HRP-labeled antibody and incubate 30 min. Finally, add HRP chemiluminescent substrates and read luminescence.

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

### Quality Assurance



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

**Lane 1:**  
 3.1 µg enzyme complex

**Lane 2:**  
 Protein Marker

**Purity:** ≥79%

