

Data Sheet

EZH2/EED/SUZ12/RbAp48 Complex

Human, recombinant

Catalog #: 51099

Lot#: 110901

Conc.: 0.6 mg/ml

Formulated in: 45 mM Tris-HCl, pH 8.0, 124 mM NaCl, 2.4 mM KCl, 3 mM DTT, 225 mM imidazole, and 10% glycerol.

Stability: >6 months at -80 °C

References:

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

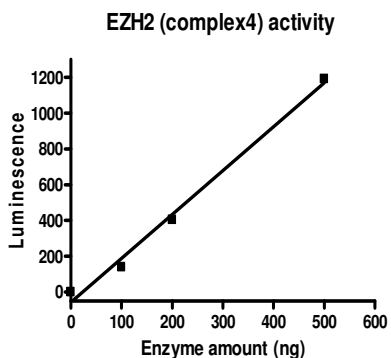
Description: Complex of human EZH2 (GenBank Accession No. NM_004456), (a.a. 2-end) with N-terminal His tag, MW= 86 kDa, human EED (NM_003797) (a.a. 2-end) with N-terminal Flag tag, MW= 51 kDa, human SUZ12 (NM_015355) (a.a. 2-end) with N-terminal His tag, MW = 87 kDa, and human RbAp48 (NM_005610) (a.a. 2-end) with N-terminal His tag, MW = 48 kDa, co-expressed in baculovirus expression system.

Specific Activity: 0.004 pmol/min/μg. Assay conditions: 50 μl reaction mix (50 mM TRIS pH 8.8, 5 mM MgCl₂, 4 mM DTT, 20 μM S-adenosylmethionine, and 50-200 ng enzyme complex) add to the wells coated with the substrate. 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

Specific Activity

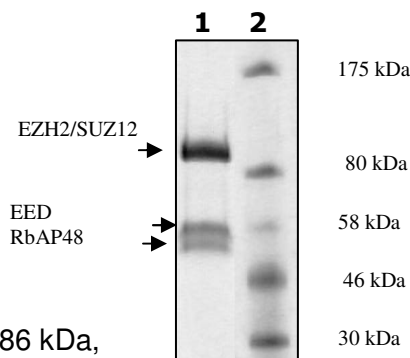


4-20% gradient SDS-PAGE Coomassie staining

Lane 1:
 4 μg enzyme complex

Lane 2:
 Protein Marker
 BioLabs (#P7708L)

MW: 87 kDa, 86 kDa,
 51 kDa, 48 kDa
Purity: >95%



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

To place your order, please contact us by Phone **1.858.829.3082** Fax **1.858.481.8694**

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

Please visit our website at: www.bpsbioscience.com