

Data Sheet

EZH2 Protein Complex

Human, Recombinant, N-terminal FLAG-Tag (EED), N-Terminal His-Tag (Rest)

Catalog #: 51004

Lot #: 161117

Conc.: 0.26 mg/ml

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

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Storing diluted enzyme is not recommended, if necessary, use carrier protein (BSA 0.1 – 0.5%).

References:

- Morin, R., *et al.* (2010). *Nature Genetics* **42**, 181 - 185
- Varambally, S. *et al.* (2008). *Science* **322 (5908)**, 1695-1699.

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, 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.

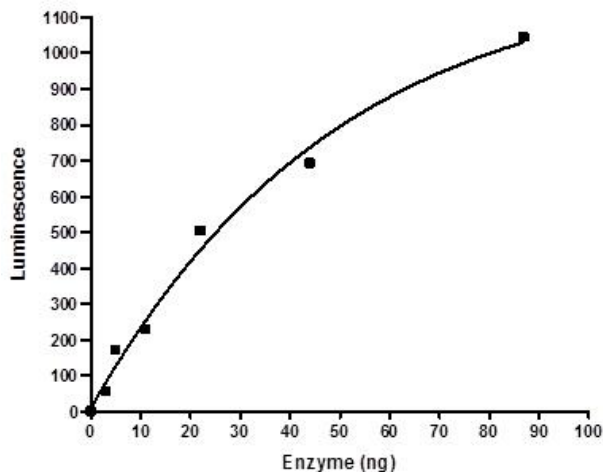
Specific Activity: ≥ 0.08 pmol/min/ μg

Assay conditions: 50 μl reaction mix (20 mM phosphate buffer pH 7.4, 0.05% Tween-20, 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



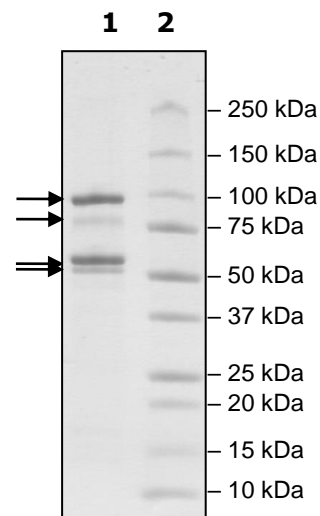
4-20% SDS-PAGE
Coomassie staining

Lane 1:
6 μg EZH2 5-complex

Lane 2:
Protein Marker

MW: EZH2 86 kDa
 SUZ12 87 kDa
 AEBP 53 kDa
 EED 51 kDa
 RbAp48 48 kDa

Purity: $\geq 90\%$



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