

Data Sheet

EZH2 A738T /EED/SUZ12/RbAp48/AEBP2

Human, Recombinant, N-terminal FLAG-Tag (EED), N-Terminal His-Tag (Rest)
Catalog #: 51061
Lot #: 161118-1 **Conc.:** 0.83 mg/ml

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

Stability: At least 6 months at -80°C . Avoid freeze/thaw cycles. Protein may be diluted to $\geq 100 \mu\text{g/ml}$ in PBS + glycerol and stored at -80°C .

References:

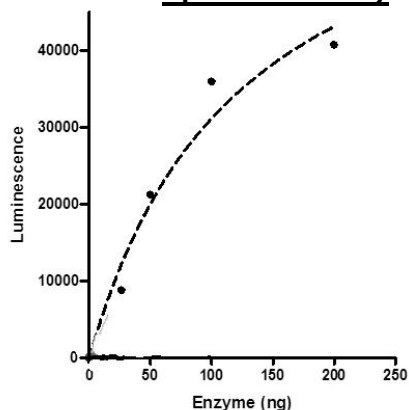
- Varambally, S. et al. Science 322 (5908), 1695-1699 (2008).
- Rakotobe, D. et al., Virol. J. 5, 32 (2008).

Description: Mutant version of our regular EZH2 5-member complex (Cat. # 51004), but with an Ala-to-Thr mutation on a.a. 738 of the EZH2 protein. Complex of human EZH2 (GenBank Accession No. NM_004456), (a-a 2-end ; A738T*) 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: 1.33 pmol/min/ μg
 50 μl reaction mix (20 mM phosphate buffer pH 7.4, 0.05% Tween-20, 10 μM S-adenosylmethionine, and various amount of 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



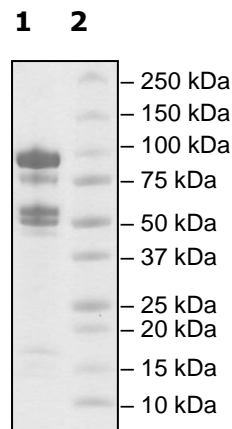
6-20% SDS-PAGE Coomassie

Lane 1:
EZH2(A738T)-5m

Lane 2:
Protein Marker

Purity: $\geq 90\%$

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



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