

Data Sheet

EZH2 Y641F /EED/SUZ12/RbAp48/AEBP2

Human, recombinant

Catalog #: 51017

Lot #: 120227 **Conc.:** 0.25 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 200 mM Imidazole, 20% Glycerol.

Stability: >6 months at -80°C

References:

1. Varambally, S. et al. *Science* 322 (5908), 1695-1699 (2008).
2. Rakotobe, D. et al., *Virology* 397, 32 (2008).

Application Reference:

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

Description:

Mutant version of our regular EZH2 5-member complex (Cat. # 51004), but with a Tyr-to-Phe mutation on a.a. 646 of the EZH2 protein. Complex of human EZH2 (GenBank Accession No. NM_004456), (a-a 2-end ; Y646F*) 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. *This mutation is often cited as Y641F in variant 3 (GenBank Accession No. NM_001203247).

Application:

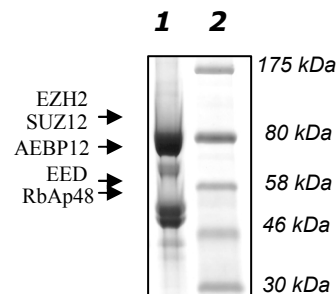
Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Assurance

10% gradient SDS-PAGE
Coomassie staining

Lane 1:
 2.5 µg enzyme complex

Lane 2:
 Protein Marker
 BioLabs (#P7708L)



Purity ≥ 91%

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