

Data Sheet

EZH2/EED

Human, recombinant

Catalog #: 51002

Lot#: 110627 Conc.: 0.35 mg/ml

Formulated in: 40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, 3 mM DTT, 200 mM imidazole, and 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., *Virol. J.* **5**, 32 (2008).

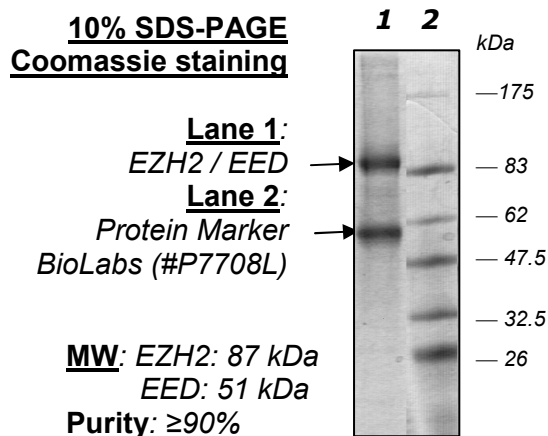
Description:

Complex of human EZH2 (GenBank Accession No. NM_004456), (amino acids 2-end) with N-terminal His tag, MW= 87 kDa, and human EED (GenBank Accession No. NM_003797) (amino acid 2-end) with N-terminal Flag tag, MW= 51 kDa, co-expressed in baculovirus expression system.

Application:

Useful for Western blot analysis and to investigate protein-protein interactions in the EZH2 complex.

Quality Assurance



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