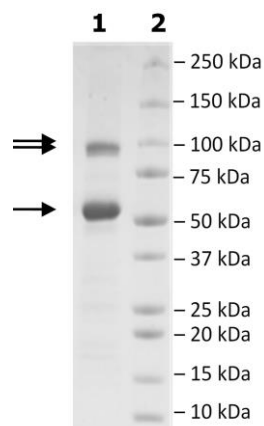


Product Information

Description:	Recombinant human EZH2 (Histone-lysine N-methyltransferase) in complex with EED (Embryonic ectoderm development protein) and SUZ (Suppressor of zeste). EZH2, full length encompassing amino acids 2-751 (end), contains an N-terminal His-tag (6xHis). Polycomb protein EED, full length encompassing amino acids 2-441(end), contains an N-terminal FLAG-tag. Polycomb protein SUZ12 full length encompassing amino acids 2-739 (end), also contains an N-terminal His-tag (6xHis). The complex was affinity purified using the FLAG-tag of EED.
Species	Human
Construct:	EZH2 (His-2-751(end)) / EED (FLAG-2-441(end)) / SUZ12 (His-1-739(end))
Concentration:	5.97 mg/ml
Expression System:	Sf9
Purity:	≥90%
Format:	Aqueous buffer solution.
Formulated In:	40 mM Tris-HCl, pH 8.0, 110 mM NaCl, 2.2 mM KCl, and 20% glycerol
MW:	EZH2: 87 kDa; EED: 51 kDa; SUZ12: 88 kDa
Genbank Accession:	EZH2: NM_004456; EED: NM_003797; SUZ12: NM_015355
Stability:	At least 6 months at -80°C.
Storage:	-80°C
Instructions for Use:	Thaw on ice and gently mix prior to use. DO NOT VORTEX. Perform a quick spin before opening. Aliquot into small volumes and flash freeze for long term storage. Avoid multiple freeze/thaw cycles.
Assay Conditions:	Assay was performed using EZH2 Chemiluminescent Assay Kit (BPS Bioscience #52009L) with EZH2 titrated from 1000 ng/rxn – 0.97 ng/rxn.
Applications:	Useful for the study of enzyme kinetics, screening inhibitors, and selectivity profiling.

Quality Control Data

4-20% SDS-PAGE Coomassie Staining



EZH2/EED/SUZ12 Activity

