

ZMYND8-[PHD-BRD-PWWP] (His)

CATALOG NO.: RD-11-472

LOT NO.:

DESCRIPTION: Human recombinant ZMYND8-[PHD-BRD] bromodomain (residues 103-426; Genbank Accession # NM_183047.2; MW = 41.0 kDa) expressed as an N-terminal 6xHis-tagged protein in *E. coli*.

PURITY: >70% by SDS-PAGE

SUPPLIED AS: _ µg/µL in 50 mM Tris HCl, pH 7.5, 500 mM NaCl, 1 mM TCEP, 10 % glycerol as determined by OD₂₈₀.

STORAGE: -70°C. Thaw quickly and store on ice before use. The remaining, unused, undiluted protein should be snap frozen, for example in a dry/ice ethanol bath or liquid nitrogen. Minimize freeze/thaws if possible, but very low volume aliquots (<5 µl) or storage of diluted enzyme is not recommended.



Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 4 µg of RBC ZMYND8-[PHD-BRD-PWWP] (His). MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, **50**, 40, 30, 25, 20, 15, 10 kDa.

This product is not intended for therapeutic or diagnostic use in animals or in humans.

Reaction Biology

1 Great Valley Parkway, Malvern PA, USA 19355

requests@reactionbiology.com www.reactionbiology.com