

BPTF-[PHD-BRD] (His) (Bromodomain and PHD finger-containing Transcription Factor; FALZ; FAC1)

CATALOG NO.: RD-11-204

LOT NO.:

DESCRIPTION: Human recombinant BPTF PHD-bromodomain construct (residues 2722-2920 (C-term.); Genbank Accession # NM_182641 MW = 26.0 kDa) expressed in *E. coli* with an N-terminal His-tag. Full-length BPTF^{1,2} is a DNA and histone-binding component of the NURF nucleosome remodeling complex (see review³). This construct comprises BPTF's second PHD zinc-finger, which primarily binds histone H3K4me3^{4,5}, and the adjacent bromodomain, which has binding affinity for various histone tail acetyllysines^{6,7}, including H4 K5Ac⁷, K12Ac⁶ and K16Ac⁶. BPTF expression is elevated in developing neurons⁸, but also in neuronal tissue under various neurodegenerative conditions^{9,10}. Amplification of the BPTF-coding chromosomal locus is prevalent in various cancers and knockdown of BPTF restricts proliferation in cultured cells with an engineered pre-malignant phenotype¹¹.

PURITY: >95% by SDS-PAGE

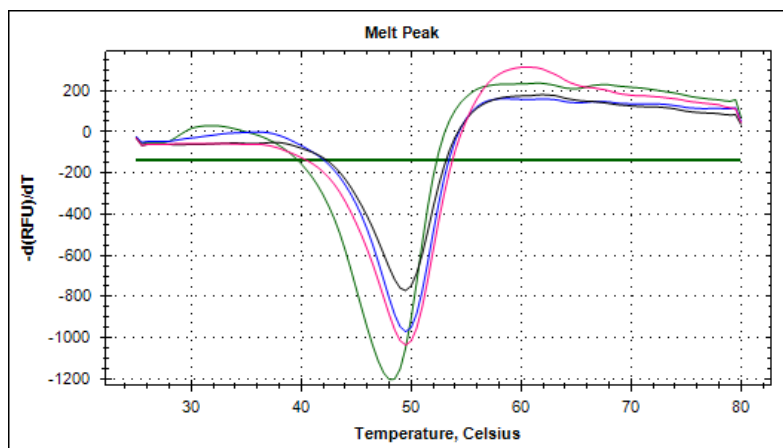
SUPPLIED AS: μ g/ μ L in 50 mM Tris-HCl, pH 7.5, 500 mM NaCl, 1 mM TCEP, 10% glycerol (v/v)

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 μ L) or storage of diluted protein is not recommended.

REFERENCES: 1) K.L. Jordan-Sciutto *et al. J. Biol. Chem.* 1999 **274** 35262; 2) M.H. Jones *et al. Genomics* 2000 **63** 35; 3) S.G. Alkhatib & J.W. Landry *FEBS Lett.* 2011 **585** 3197; 4) H. Li *et al. Nature* 2006 **442** 91; 5) H. Li *et al. Mol. Cell* 2007 **28** 677; 6) A.J. Ruthenburg *et al. Cell* 2011 **145** 692; 7) P. Filippakopoulos *et al. Cell* 2012 **149** 214; 8) K.L. Jordan-Sciutto *et al. Biochem. Biophys. Res.* 1999 **260** 785; 9) S. Schoonover *et al. J. Neuropathol. Exp. Neurol.* 1996 **55** 444; 10) X. Mu *et al. Exp. Neurol.* 1997 **146** 17; 11) Y. Buganim *et al. PLoS One* 2010 **5** e9657



Coomassie blue-stained SDS-PAGE (12% acrylamide) of 5 μ g of RBC BPTF-[PHD-BRD] (His). MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.



Differential Scanning Fluorimetry of RBC BPTF-[PHD-BRD] (His) in the Presence or Absence of Common Bromodomain Ligands. Thermal denaturation of BPTF-[PHD-BRD] (His) (0.1 mg/mL) is detected (CFX384 TMTouch thermal cycler, 'FRET' channel; Bio- Rad) by increased binding and fluorescence of the dye SYPRO[®]Orange (Life Technologies). Shown are the first derivatives of fluorescence vs. temperature plots in which the "peaks" indicate the T_m's (inflection points) of the original plots. Addition of either PF11 (pink) or CBP112 (black) or Bromosporine (blue) (all 25 μ M) stabilizes the protein folding and shifts the T_m (inflection point) from 48°C to 49.5°C.

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