

SP140-[PHD-BRD] (His) (Speckled 140 kDa Nuclear Body Protein, PHD-Bromodomain)

CATALOG NO.: RD-11-198

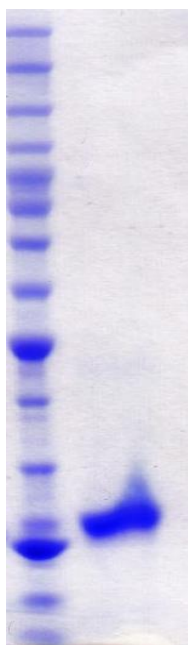
LOT NO.:

DESCRIPTION: Human recombinant SP140 PHD-bromodomain construct (residues 688-862; Genbank Accession # NM_007237; MW = 23.6 kDa) expressed with an N-terminal His-tag in *E. coli*.

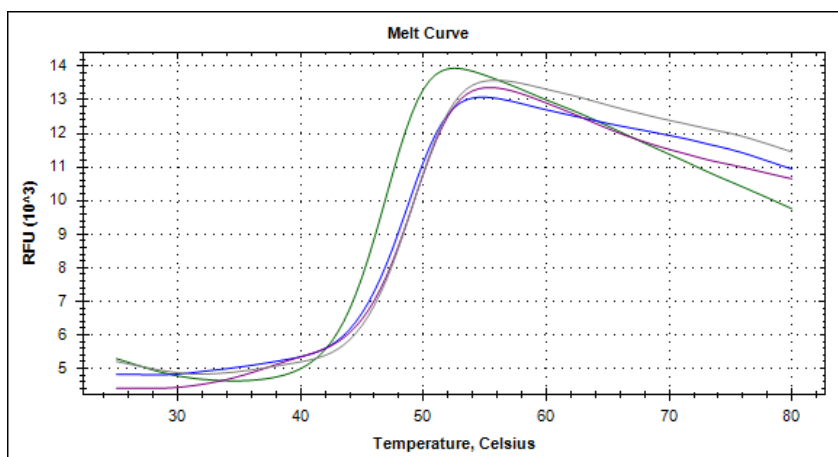
PURITY: >80% by SDS-PAGE

SUPPLIED AS: µg/µL in 50 mM Tris HCl, pH 7.5, 500 mM NaCl, 1 mM TCEP, 10 % glycerol

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 (12% acrylamide) of 5 µg of RBC SP140-[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 SP140-[PHD-BRD] (His) Thermal denaturation of SP140-[PHD-BRD] (His) is detected (CFX384 TMTouch thermal cycler, 'FRET' channel; Bio- Rad) by increased binding and fluorescence of the dye SYPRO®Orange (Life Technologies). Addition of 25 µM PF11 (grey), CBP112 (purple) and Bromosporine (blue) stabilizes the protein folding and shifts the T_m (inflection point) from 47°C to 49°C, 49°C and 48.5°C respectively.

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