

SP140L-[PHD-BRD] (His)

(Nuclear body protein SP140-like protein)

CATALOG NO.: RD-11-201

LOT NO.:

DESCRIPTION: Human recombinant SP140L PHD-bromodomain construct (residues 401-580 (C-term.); Genbank Accession # NM_138402; MW = 24.0 kDa) expressed in *E. coli* with an N-terminal His-tag.

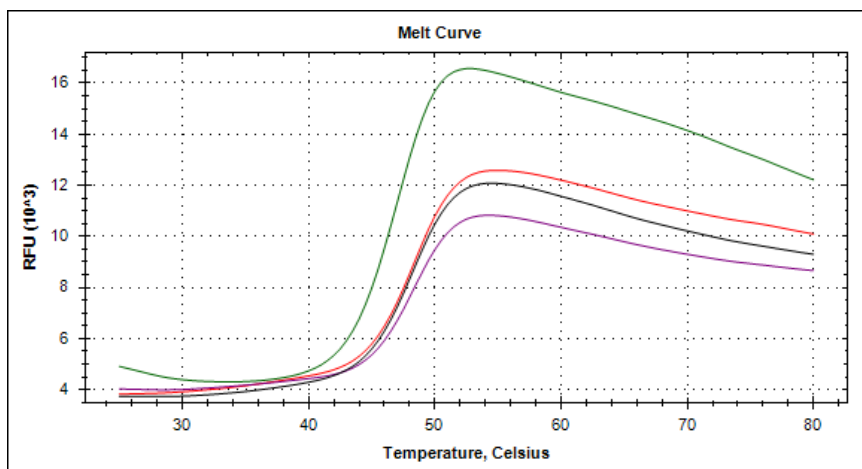
PURITY: >90% by SDS-PAGE

SUPPLIED AS: $_ \mu\text{g}/\mu\text{L}$ in 50 mM Tris HCl, pH 7.5, 500 mM NaCl, 1 mM TCEP, 10% glycerol (v/v) 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 protein is not recommended.



Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 4 μg of RBC SP140L-[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 SP140L-[PHD-BRD] (His) in presence or absence of common bromodomain ligands. Thermal denaturation of SP140L-[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 JQ1 (black), CBP112 (purple) or SGC-CBP30 (red) stabilizes the protein folding and shifts the T_m (inflection point) from 47°C to 48°C, 48.5°C or 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