

MLL-[PHD-BRD] (His)

CATALOG NO.: RD-11-318

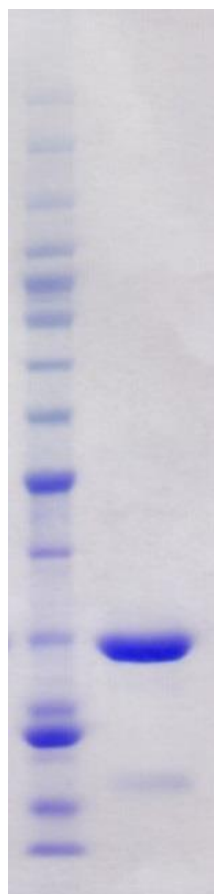
LOT NO.:

DESCRIPTION: Human recombinant MLL PHD-bromodomain construct (residues 1566-1784; Genbank Accession # NM_005933.3; MW = 27.8 kDa) expressed with an N-terminal His-tag in *E. coli*.

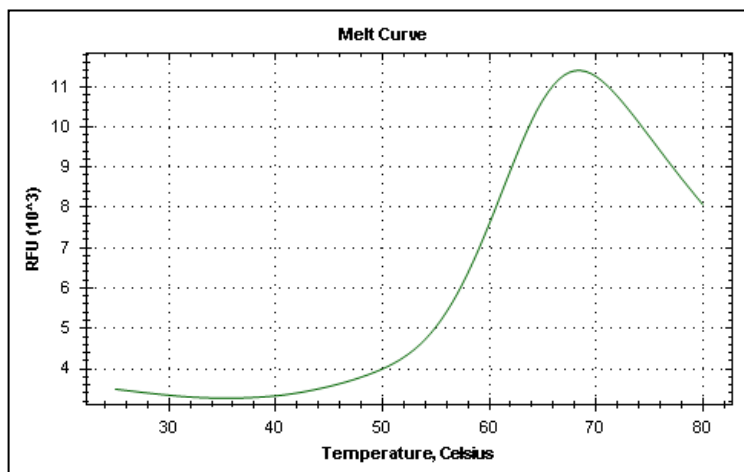
PURITY: >85% by SDS-PAGE

SUPPLIED AS: $_ \mu\text{g}/\mu\text{L}$ in 50 mM HEPES pH 7.5, 500 mM NaCl, 10% glycerol, 1 mM TCEP 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 MLL-[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 MLL-[PHD-BRD] (His). Thermal denaturation of MLL-[PHD-BRD] (His) is detected (CFX384™ Touch thermal cycler, 'FRET' channel; Bio-Rad) by increased binding and fluorescence of the dye SYPRO® Orange (Life Technologies). The apo form of MLL-[PHD-BRD] (His) displays a T_m of 61.0°C and is not stabilized in the presence of various known bromodomain ligands (JQ1, PFI1, CBP112, Bromosporine, SGC-CBP30, BET151, RVX-208, GSK2801 and PFI-3; not shown; all tested at 25 μM).

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