

## ASH1L-[BRD] (GST)

(ASH-1 like protein; huASH1; KMT2H)

**CATALOG NO.:** RD-11-240

**LOT NO.:**

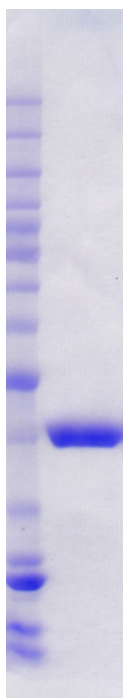
**DESCRIPTION:** Human recombinant ASH1L bromodomain (residues 2428-2559; Genbank Accession # NM\_018489; MW = 42.0 kDa) expressed in *E. coli* with an N-terminal GST-tag. ASH1L is a large (333 kDa), multi-domain protein associated with actively transcribed regions of chromatin. Its bromodomain lies C-terminal to its SET domain, which confers histone H3K36 methyltransferase activity<sup>1-3</sup>. ASH1L is the human homolog of *Drosophila* Ash1, a Trithorax group protein. Like its counterpart in *Drosophila*<sup>4</sup>, ASH1L has been found to play a role in the regulation of Hox gene expression<sup>5,6</sup>. The ASH1L bromodomain displays strong binding to various Lys(Ac) residues in singly acetylated histone peptide microarrays (histones H1.4K74Ac, H2AK36Ac, H2BK85Ac, H3K56Ac, H4K59Ac/K79Ac)<sup>7</sup>.

**PURITY:** >95% by SDS-PAGE

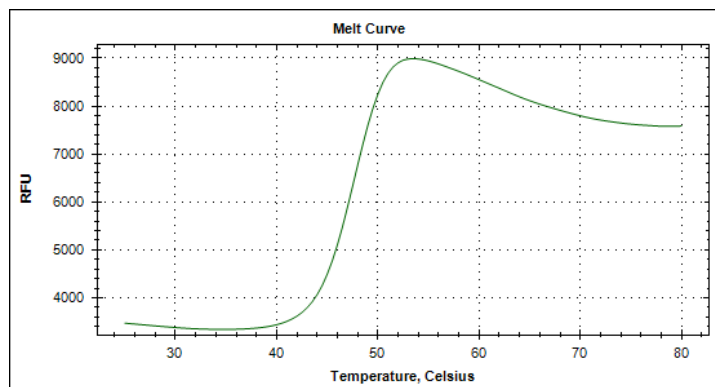
**SUPPLIED AS:**  $\mu\text{g}/\mu\text{L}$  in 50mM Tris HCl, pH 7.5, 500mM NaCl, 1mM TCEP, 10% glycerol (v/v) as determined by OD<sub>280</sub>

**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  $\mu\text{l}$ ) or storage of diluted protein is not recommended.

**REFERENCES:** 1) Y Tanaka *et al. Gene* 2007 **397** 161; 2) S. An *et al. J. Biol. Chem.* 2011 **286** 8369; 3) D.S. Cabianca *et al. Cell* 2012 **149** 819; 4) C. Beisel *et al. Nature* 2002 **419** 857; 5) G. Gregory & *et al. Mol. Cell. Biol.* 2007 **27** 8466; 6) Y Tanaka *et al. PLOS One* 2011 **6** e28171; 7) P. Filippakopoulos *et al. Cell* 2012 **149** 214



**Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 4  $\mu\text{g}$  of RBC ASH1L-[BRD] (GST).** 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 ASH1L-[BRD] (GST) in presence or absence of common bromodomain ligands.**

Thermal denaturation of ASH1L-[BRD](GST) is detected (CFX384 TMTouch thermal cycler, 'FRET' channel; Bio- Rad) by increased binding and fluorescence of the dye SYPRO®Orange (Life Technologies). Apo form of ASH1L-BRD-G displays a T<sub>m</sub> of 47.5°C and is not stabilized in the presence of various known bromodomain ligands (JQ1, PF11, CBP112, Bromosporine, SGC-CBP30, BET151 and RVX-208; all tested at 25  $\mu\text{M}$ ).

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