

BRD9 (GST)

(Bromodomain-containing protein 9)

CATALOG NO.: RD-11-187

LOT NO.:

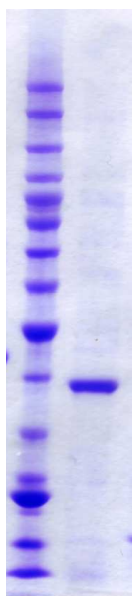
DESCRIPTION: Human recombinant BRD9 bromodomain (residues 135-248; Genbank Accession # NM_023924; MW = 40.1 kDa) expressed as an N-terminal GST-fusion protein in *E. coli*. BRD9 has been identified as a subunit of the BAF-subtype of SWI/SNF nucleosome remodeling complexes^{1,2}. The SWI/SNF complexes are the most highly mutated chromatin regulator in human cancers (19.6% in 44 published sequencing studies), although BRD9 subunit is among the least mutated subunits². However, increased copy numbers of the BRD9 gene occur with high frequency in early stage non-small cell lung cancers³ and in invasive cervical cancers⁴. Triazolophthalazine inhibitors exhibiting low micromolar affinity and some partial selectivity for the BRD9 bromodomain (along with CREBBP, CECR2 and/or BRD4 bromodomains) have been synthesized⁵ and crystal structures of the BRD9 bromodomain determined both with⁵ and without⁶ one such ligand. The BRD9 bromodomain was found to bind, *in vitro*, to various microarrayed histone H3, singly lysine-acetylated peptides⁶.

PURITY: >90% by SDS-PAGE

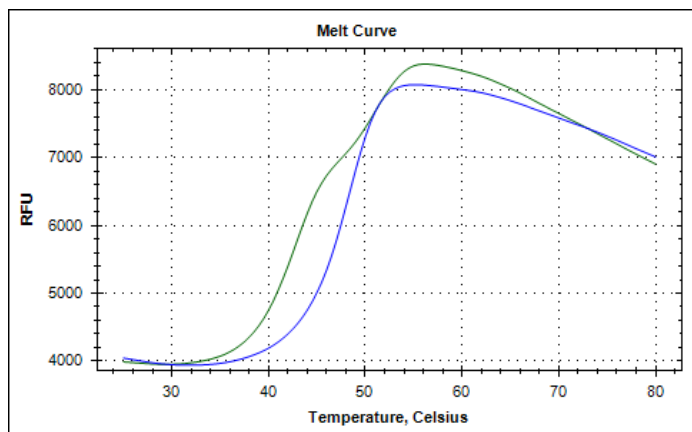
SUPPLIED AS: _ $\mu\text{g}/\mu\text{L}$ in 50 mM HEPES/NaOH, pH 7.0, 500 mM NaCl, 1 mM TCEP, 10 % glycerol as determined by OD₂₈₀

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.

REFERENCES: 1) E. Middeljans *et al. PLoS One* 2012 **7** e33834; 2) C. Kadoch *et al. Nature Genet.* 2013 **45** 592; 3) J.U. Kang *et al. Cancer Genet. Cytogenet.* 2008 **182** 1; 4) L. Scotto *et al. Mol. Cancer* 2012 **7** 58; 5) O. Fedorov *et al. J. Med. Chem.* 2014 **57** 462; 6) P. Filippakopoulos *et al. Cell* 2012 **149** 214



Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 3.6 μg of RBC BRD9 (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 BRD9 (GST) Thermal denaturation of BRD9 (GST) is detected (CFX384™ Touch thermal cycler, 'FRET' channel; Bio-Rad) by increased binding and fluorescence of the dye SYPRO® Orange (Life Technologies). Addition of 25 μM bromosporine stabilizes the protein folding and shifts the T_m (inflection point) from 43°C to 48.5°C.

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