

**BRD8-1 (GST)**

(Bromodomain containing protein 8, bromodomain 1)

**CATALOG NO.:** RD-11-305

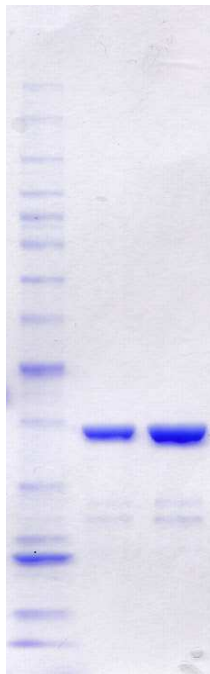
**LOT NO.:**

**DESCRIPTION:** Human recombinant BRD8, bromodomain 1 (residues 701-819; Genbank Accession # NM\_139199.1; MW = 40.9 kDa) expressed as an N-terminal GST-fusion protein in *E. coli*.

**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 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 enzyme is not recommended.



**Coomassie blue-stained SDS-PAGE (4-12% acrylamide) of 1 and 2  $\mu\text{g}$  of RBC BRD8-1 (GST).** MW markers (left) are, from top, 220, 160, 120, 100, 90, 80, 70, 60, 50, 40, 30, 25, 20, 15, 10 kDa.

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