

**ATAD2 (His)**

**CATALOG NO.:** RD-11-251

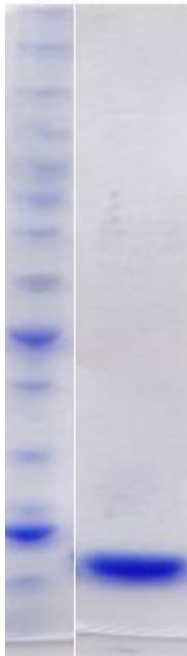
**LOT NO.:**

**DESCRIPTION:** Human recombinant ATAD2 bromodomain (residues 981-1146; Genbank Accession # NM\_014109; MW = 17.95 kDa) expressed as an N-terminal His-fusion protein in *E. coli*.

**PURITY:** >95% 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

**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 (12% acrylamide) of 5  $\mu\text{g}$  of RBC ATAD2 (His).** 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

