

**BRPF3 (GST)**

(Bromodomain and PHD-finger containing protein 3)

**CATALOG NO.:** RD-11-188

**LOT NO.:**

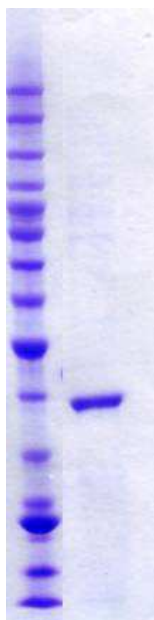
**DESCRIPTION:** Human recombinant BRPF3 bromodomain (residues 588-703; Genbank Accession # NM\_015695; MW = 40.7 kDa) expressed as an N-terminal GST-fusion protein in *E. coli*. A scaffolding component of the MOZ/MORF histone acetyltransferase complex<sup>1,2</sup>, native BRPF3 contains, in addition to its acetyllysine-binding bromodomain, two other presumed histone/chromatin binding domains, a PWWP domain<sup>3</sup> and a PHD zinc-finger.

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

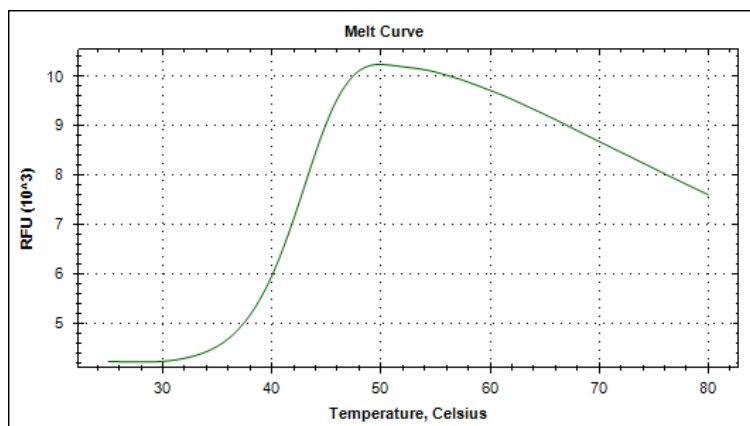
**SUPPLIED AS:** \_ µg/µL in 50 mM HEPES, pH 7.5, 500 mM NaCl, 1 mM TCEP, 10 % glycerol as determined by OD<sub>280</sub>.

**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 µl) or storage of diluted enzyme is not recommended.

**REFERENCES:** 1) Y. Doyon *et al. Mol. Cell* 2006 **21** 51; 2) M. Ullah *et al. Mol. Cell. Biol.* 2008 **28** 6828; 3) H. Wu *et al. PLoS One* 2011 **6** e18919



**Coomassie blue stained SDS-PAGE (4-12% acrylamide) of 3 µg of RBC BRPF3 (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 BRPF3 (GST)**

Thermal denaturation of BRPF3 (GST) is detected (CFX384™ Touch thermal cycler, 'FRET' channel; Bio-Rad) by increased binding and fluorescence of the dye SYPRO® Orange (Life Technologies). Apo form of BRPF3(GST) displays a T<sub>m</sub> of 43°C and is not stabilized in the presence of various known bromodomain ligands (JQ1, PF1, CBP112, Bromosporine, SGC-CBP30, BET151 and RVX-208; all tested at 25 µM; not shown).

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