



Bcr (Ab-177) Antibody

Catalog Number 21197

(For Research Use Only)

- Catalog Number:** 21197-1, 21197-2, 21197
Amount: 50µg/50µl, 100µg/100µl, 200µg/200µl
Swiss-Prot No. : P11274
All Names: BCR, BCR protein, BCR1, Breakpoint cluster region protein
Form of Antibody: Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Storage/Stability: Store at -20 °C /1 year
Immunogen: The antiserum was produced against synthesized non-phosphopeptide derived from human Bcr around the phosphorylation site of tyrosine 177 (P-F-Y^P-V-N).
Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity/Sensitivity Bcr (Ab-177) antibody detects endogenous levels of total Bcr protein.
Reactivity: Human, Mouse, Rat
Applications: WB: 1:500~1:1000 ELISA: 1:20000
References: ML Campbell, Cell Growth Differ., Jul 1993; 4: 581.
Asok Mukhopadhyay, J. Biol. Chem., Aug 2002; 277: 30622 - 30628.
Jeanne Wilson-Rawls, Cancer Res., Aug 1996; 56: 3426 - 3430.
Lori Rink, Blood, Apr 2007; 10.1182.

