



## Bcr (Phospho-Tyr177) Antibody

Catalog Number 11199

(For Research Use Only)

**Catalog Number:** 11199-1, 11199-2, 11199

**Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl

**Swiss-Prot No. :** P11274

**Form of Antibody:** The antiserum was produced against synthesized phosphopeptide derived from human Bcr around the phosphorylation site of tyrosine 177 (P-F-Y<sup>P</sup>-V-N).

**Storage/Stability:** Store at -20°C/1 year

**Immunogen:** The antiserum was produced against synthesized phosphopeptide derived from human SMC1 around the phosphorylation site of serine 957 (G-S-S<sup>P</sup>-Q-G).

**Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific phosphopeptide. The antibody against non-phosphopeptide was removed by chromatography using non-phosphopeptide corresponding to the phosphorylation site.

**Specificity/Sensitivity:** Bcr (phospho-Tyr177) antibody detects endogenous levels of Bcr only when phosphorylated at tyrosine 177.

**Reactivity:** Human, Mouse

**Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100

**References:** Sattler M, et al. Mol Cell Biol. 1999 Nov; 19(11): 7473-7480.

Li S et al. EMBO J. 2001 Dec 3; 20(23): 6793-6804.

Million RP, et,al. Mol Cell Biol. 2004 Jun; 24(11): 4685-4695.

Malcolm A. Meyn, et al. . Biol. Chem., Oct 2006; 281: 30907 – 30916..

