

c-Kit (Ab-721) Antibody [#B7138]

- Catalog Number:** B7138-1, B7138-2, B7138
Amount: 50µg/50µl, 100µg/100µl, 200µg/200µl
Swiss-Prot No. : P10721
All Names: CD117, Mast/stem cell growth factor receptor precursor, SCFR, SL, c-kit, kinase Kit
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 c-Kit around the phosphorylation site of tyrosine 721 (N-E-Y^P-M-D).
Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity/Sensitivity: c-Kit (Ab-721) antibody detects endogenous levels of total c-Kit protein.
Reactivity: Human, Mouse
Applications: WB: 1:500~1:1000 ELISA: 1:10000
References: G. Sakuntala Warshamana-Greene Mol, et al. (2004) Cancer Ther. 3: 527-536
Y Kubota, et al. (1998) Cell Growth Differ. 9: 247-256
Pramod Pandey, et al. (1995) Cancer Res. 55: 4000-4003.
Nancy K. Pryer, Clin. et al. (2003) Cancer Res. 9: 5729-5734