

FAK (Ab-407) Antibody [#B0479]

Catalog Number:	B0479
Concentration:	1mg/ml
Swiss-Prot No.:	Q05397
Other Names:	EC 2.7.10.2, FADK 1, FAK1, Focal adhesion kinase 1, PTK2, Protein- tyrosine kinase 2, pp125FAK
Storage/Stability:	Store at -20°C/1 year
Form of Antibody:	Rabbit IgG in phosphate buffered saline (without Mg ²⁺ and Ca ²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
Immunogen:	The antiserum was produced against synthesized non-phosphopeptide derived from human FAK around the phosphorylation site of tyrosine 407 (D-T-Y ^P -T-M).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	FAK (Ab-407) antibody detects endogenous levels of total FAK protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
References:	Graham Jones. Cancer Res., Aug 2001; 61: 5688 - 5691. Chuanhai Zheng. J. Biol. Chem., Jan 1998; 273: 2384. Pamela J. Bruce-Staskal. J. Cell Sci., Jul 2002; 115: 2689.