

Met (Ab-1349) Antibody [#B7152]

Catalog Number: B7152-1, B7152-2, B7152

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

Swiss-Prot No. : P08581

All Names: HGF receptor, HGF-SF receptor, Hepatocyte growth factor receptor precursor, Met proto-oncogene tyrosine kinase, c-met, kinase Met

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 Met around the phosphorylation site of tyrosine 1349 (E-H-Y^P-V-H).

Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.

Specificity/Sensitivity: Met (Ab-1349) antibody detects endogenous levels of total Met protein. This antibody may cross-react with other activated protein tyrosine kinases.

Reactivity: Human, Mouse, Rat

Applications: WB: 1:500~1:1000 ELISA: 1:10000

References: Fan S, et al. (2001) Mol Cell Biol; 21(15): 4968-4984.

Schiering N, et al. (2003) Proc Natl Acad Sci U S A; 100(22): 12654-12659.

Plopper GE, et al. (1995) Mol Biol Cell; 6(10): 1349-1365.

Ponzetto C, et al. (1993) Mol Cell Biol; 13(8): 4600-4608.

Jackson PA, et al. (2001) Plant Physiol; 127(3): 1065-1076.