

MMP-23 Antibody [#C0271]

Catalog Number: C0271-1, C0271-2, C0271

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

Swiss-Prot No. : Q13561

All Names: Matrix metalloproteinase-23 [Precursor], EC 3.4.24.-, Matrix metalloproteinase 21, MMP-21
Matrix metalloproteinase 22, MMP-22, Femalysin, MIFR-1

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 peptide derived from human MMP-23 antibody.

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

Specificity/Sensitivity MMP-23 antibody detects endogenous levels of total MMP-23 protein.

Reactivity: Human, Mouse, Rat

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

References: Scott S. Grieshaber, J. Cell Sci., Sep 2003; 116: 3793.

Janis K. Burkhardt, J. Cell Biol., Oct 1997; 139: 469.

Karin A. Melkonian, J. Biol. Chem., Apr 2007; 10.1074.

Lili Yue, J. Biol. Chem., Jul 2000; 275: 23948.