

Fibrillin-1 Antibody [#C0194]

- Catalog Number:** C0194-1, C0194-2, C0194
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** P35555
- All Names:** MASS, MFS1, OCTD, SGS, WMS, 350 kda glycoprotein component extracellular microfibril, Weill-Marchesani syndrome, fibrillin 1 (Marfan syndrome), fibrillin 15
- 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.
- Storage/Stability:** Store at -20°C/1 year
- Immunogen:** The antiserum was produced against synthesized peptide derived from human Fibrillin-1.
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** Fibrillin-1 antibody detects endogenous levels of total Fibrillin-1 protein.
- Reactivity:** Human (Identities = 100%, Positives = 100%);
Mouse (Identities = 92%, Positives = 100%);
Rat (Identities = 92%, Positives = 100%)
- Applications:** IHC: 1:50~1:100 ELISA: 1:40000
- References:** Ehab El-Hallous, J. Biol. Chem., Mar 2007; 282: 8935 - 8946.
Shazia S. Chaudhry, J. Cell Biol., Jan 2007; 176: 355 - 367.
Kerstin Tiedemann, J. Biol. Chem., Sep 2001; 276: 36035 - 36042.
Liliana Schaefer, Am. J. Pathol., Aug 2004; 165: 383 - 396.