



Profilin polyclonal antibody

Cat. # APUF01

Lot: 035

Source: Profilin antibody is provided as an affinity purified rabbit polyclonal antibody. The antibody has been raised against purified human profilin and it is predicted to have broad species cross reactivity. The positive control protein consists of an extract of bovine thymus that identifies a characteristic profilin band at 14 kDa. The control protein (bovine thymus extract, Cat. # EXT03) is supplied as a lyophilized powder; this should be reconstituted in 500 μ l of cold Milli-Q water to 2 mg/ml. After addition of water the protein should be left at room temperature for 5 minutes with occasional pipetting up and down to make sure that the protein has completely dissolved into solution, it can then be stored at -20°C . We recommend running 10 μ g of positive control protein in Western blots.

Materials: The profilin antibody (50 μ g) is supplied as a lyophilized powder. It is recommended to resuspend the powder in 200 μ l of 30% glycerol/water. This gives a concentration of 250 μ g/ml in PBS plus 30% glycerol with 0.01% chlorhexidine for preservative.

Recommended Working Dilution The profilin antibody should be used for Western blots at a 1:1000 in TBST.

Specificity: The profilin antibody will cross-react with profilin of mouse, rat, and human origin, and there are no non-specific bands in thymus and platelet tissue extracts (see figure 1).



Figure 1: **Western blot probed with profilin antibody.** Lane 1; 100 ng profilin protein, Lane 2; 10 μ g thymus extract protein, Lane 3; 10 μ g of platelet extract protein. Primary antibody used at 1:1000 dilution, secondary antibody used at 1/50 000 dilution anti-rabbit HRP, Supersignal West Dura™ (Pierce Co.) detection method with film exposure for 30 seconds.

Storage: Store antibody at 4°C . Store the control protein at -20°C or -70°C .