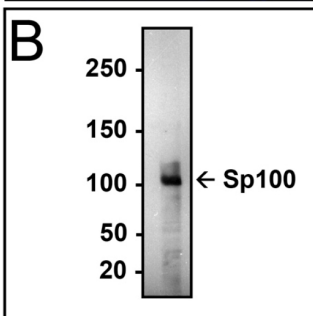
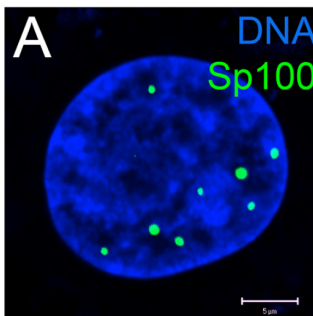


anti-Sp100 Peptide anti-human Sp100 rabbit, polyclonal

Cat. No.	Amount
ABD-031	50 μ l



For *in vitro* use only
Quality guaranteed for 12 months
Store in aliquots at -20°C

Avoid freeze / thaw cycles

Form

Liquid. Supplied as serum.

Description

Polyclonal Antibody pABahSp100(209-227)#2 to human Sp100.

Immunogen

The antibody binds to a peptide corresponding to amino acids 209-227 (CETEQINAKRKDTTSDKDD) of the human Sp100 protein.

Specificity

Antibody specificity was evaluated by indirect immunofluorescence and western blotting (see Figure 1A and B)

Application

Indirect Immunofluorescence, concentration: 1:1000 to 1:2000.

Western Blot, concentration: 1:500.

Selected References:

Szostecki *et al.* (1990) Isolation and characterization of cDNA encoding a human nuclear antigen predominantly recognized by autoantibodies from patients with primary biliary cirrhosis. *J Immunol.* **145**:4338.