

SREBP-1 (Ab-439) Antibody [#B8425]

Catalog Number:	B8425
Concentration:	1mg/ml
Swiss-Prot No.:	P36956
Other Names:	SRBP1; SREBF1; SREBP1; sterol regulatory element-binding protein 1; sterol regulatory element-binding transcription factor 1
Storage/Stability:	Store at -20°C/1 year
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.
Immunogen:	The antiserum was produced against synthesized non-phosphopeptide derived from human SREBP-1 around the phosphorylation site of serine 439 (Q-S ^P -P-L).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	SREBP-1 (Ab-439) antibody detects endogenous levels of total SREBP-1 protein.
Reactivity:	Human (Identities = 100%, Positives = 100%); Rat (Identities = 90%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:20000
References:	Yokoyama C., Cell 75:187-197(1993). Hua X., Genomics 25:667-673(1995). Olsen J.V., Cell 127:635-648(2006).