

Calnexin (Ab-583) Antibody [#B0463]

Catalog Number:	B0463
Concentration:	1mg/ml
Swiss-Prot No.:	P27824
Other Names:	CALX, CANX, IP90, Major histocompatibility complex class I antigen-binding protein p88, p90
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 Calnexin around the phosphorylation site of serine 583 (N-R-S ^P -P-R).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Calnexin (Ab-583) antibody detects endogenous levels of total Calnexin protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	Shaun P Brothers, J. Mol. Endocrinol., Dec 2006; 37: 479 - 488. Vilasack Thammavongsa, J. Biol. Chem., Sep 2005; 280: 33497 - 33505. Yasushi Okazaki, J. Biol. Chem., Nov 2000; 275: 35751.