

Arrestin 1 (Ab-412) Antibody [#B0455]

Catalog Number:	B0455
Concentration:	1mg/ml
Swiss-Prot No.:	P49407
Other Names:	ARR1, arrestin 2, arrestin beta 1, beta-arrestin-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 Arrestin 1 around the phosphorylation site of serine 412 (T-G-S ^P -P-Q).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Arrestin 1 (Ab-412) antibody detects endogenous levels of total Arrestin 1 protein.
Reactivity:	Human
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
References:	D. Alex Groarke. Mol. Pharmacol., Feb 2001; 59: 375. Karen McConalogue. Mol. Biol. Cell, Aug 1998; 9: 2305. Maria Stella Lombardi. Stroke, Apr 2004; 35: 981 - 986.