

Rhodopsin (Ab-334) Antibody [#B0569]

Catalog Number:	B0569
Concentration:	1mg/ml
Swiss-Prot No.:	P08100
Other Names:	OPN2, OPSD, RHO, opsin 2
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 Rhodopsin around the phosphorylation site of serine 334 (D-A-S ^P -A-T).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	Rhodopsin (Ab-334) antibody detects endogenous levels of total Rhodopsin protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:5000
References:	Edwin S. Lee, Invest. Ophthalmol. Vis. Sci., Jun 2007; 48: 2868 - 2876. DJ Roof, Invest. Ophthalmol. Vis. Sci., Nov 1994; 35: 4049. Angelika Schmitt, J. Exp. Biol., Apr 2005; 208: 1247 - 1256. Paul A. Sieving, J. Neurosci., Aug 2001; 21: 5449.