

PKR (Ab-258) Antibody [#B8174]

Catalog Number:	B8174
Concentration:	1mg/ml
Swiss-Prot No.:	P19525
Other Names:	ADRB2; E2AK2; EC 2.7.11.1; EIF2aK; EIF2AK2; Interferon-induced, double-stranded RNA-activated protein kinase; Interferon-inducible RNA-dependent protein kinase; kinase PKR; P1/eIF-2A protein kinase; p68 kinase; PRKR; TIK
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 PKR around the phosphorylation site of threonine 258 (K-Y-T ^P -V-D).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PKR (Ab-258) antibody detects endogenous levels of total PKR protein.
Reactivity:	Human (Identities = 100%, Positives = 100%)
Applications:	WB: 1:500~1:1000 ELISA: 1:40000
References:	Meurs E., Cell 62:379-390(1990). Thomis D.C., Virology 188:33-46(1992). Kuhlen K.L., Gene 178:191-193(1996).