

PEA-15 (Ab-116) Antibody [#B0545]

Catalog Number:	B0545
Concentration:	1mg/ml
Swiss-Prot No.:	Q15121
Other Names:	Astrocytic phosphoprotein PEA-15, PE15, PEA15, PED, Phosphoprotein enriched in diabetes
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 PEA-15 around the phosphorylation site of serine 116 (Q-P-S ^P -E-E).
Purification:	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity:	PEA-15 (Ab-116) antibody detects endogenous levels of total PEA-15 protein.
Reactivity:	Human, Mouse, Rat
Applications:	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:40000
References:	Joseph Krueger, Mol. Biol. Cell, Aug 2005; 16: 3552 - 3561. Alessandra Trencia, Mol. Cell. Biol., Jul 2003; 23: 4511 - 4521. Daniel Kitsberg, J. Neurosci., Oct 1999; 19: 8244. Angela Glading, Cancer Res., Feb 2007; 67: 1536 - 1544.