

PKR (Ab-446) Antibody [#B7198]

- Catalog Number:** B7198
- Amount:** 50µg/50µl, 100µg/100µl, 200µg/200µl
- Swiss-Prot No. :** P19525
- All Names:** EIF2AK2 , Interferon-induced double-stranded RNA-activated protein kinase
- Form of Antibody:** Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
- Storage/Stability:** Store at -20°C/1 year
- Immunogen:** The antiserum was produced against synthesized non-phosphopeptide derived from human PKR around the phosphorylation site of threonine 446 (K-R-T^P-R-S).
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity:** PKR (Ab-446) antibody detects endogenous levels of total PKR protein.
- Reactivity:** Human
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:20000
- References:** Abujiang Pataer, et.al. (2002) Cancer Res; 62: 2239.
K. D. Ryman, et.al. (2005) J. Virol; 79: 1487-1499.
Susana Guerra, et.al. (2006) J. Biol. Chem; 281: 18734-18745.