

GSK3 α / β (Ab-279/216) Antibody [#B0012]

- Catalog Number:** B0012-1, B0012-2, B0012
- Amount:** 50 μ g/50 μ l, 100 μ g/100 μ l, 200 μ g/200 μ l
- Swiss-Prot No. :** P49840
- All Names:** EC 2.7.11.26, Factor A, GSK-3 alpha, GSK3-alpha, Glycogen synthase kinase-3 alpha, kinase GSK3-alpha, GSK-3 beta, GSK3B, Glycogen synthase kinase-3 beta, Protein kinase GSK-3-beta, kinase GSK3-beta
- 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 GSK3 α / β around the phosphorylation site of tyrosine 279/216 (V-S-Y-I-C).
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** GSK3 α / β (Ab-279/216) antibody detects endogenous levels of total GSK3 α / β protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100
ELISA: 1:20000
- References:** Minsub Shim, J. Biol. Chem., May 2003; 278: 19674 -19681.
Junichi Gotoh Carcinogenesis, Mar 2003; 24: 435 - 442.
Katrina MacAulay, J. Biol. Chem., Mar 2005; 280: 9509 - 9518.
Victor H. Bustos, PNAS, Dec 2006; 103: 19725 - 19730.