



cdc25A (Ab-75) Antibody

Catalog Number 21163

(For Research Use Only)

- Catalog Number:** 21163-1, 21163-2, 21163
Amount: 50µg/50µl, 100µg/100µl, 200µg/200µl
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 cdc25A around the phosphorylation site of serine 75.
Purification: The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
Specificity/Sensitivity: cdc25A (Ab-75) antibody detects endogenous levels of total cdc25A protein.
Reactivity: Human, Mouse, Rat
Applications: WB: 1:500~1:1000 ELISA: 1:10000
Swiss-Prot No. : P30304
References: Jin J, et al. (2003) Genes 17(24): 3062-74.
Shimuta K, et al. (2002) EMBO J 21(14): 3694-703.

