

## c-PLA2 (Ab-505) Antibody [#B0065]

<b>Catalog Number:</b>	B0065
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	P47712
<b>Other Names:</b>	CPLA2, Cytosolic phospholipase A2, EC 3.1.1.4, EC 3.1.1.5, PA24, PA24A, PLA2G4, PLA2G4A, Phosphatidylcholine 2-acylhydrolase
<b>Storage/Stability:</b>	Store at -20°C/1 year
<b>Form of Antibody:</b>	Rabbit IgG in phosphate buffered saline (without Mg <sup>2+</sup> and Ca <sup>2+</sup> ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
<b>Immunogen:</b>	The antiserum was produced against synthesized non-phosphopeptide derived from human c-PLA2 around the phosphorylation site of serine 505 (P-L-S <sup>P</sup> -P-L).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	c-PLA2 (Ab-505) antibody detects endogenous levels of total c-PLA2 protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:10000
<b>References:</b>	D. D. Denson, J Neurophysiol, Jun 2005; 93: 3146 - 3156. Diane M. Duffy, J. Clin. Endocrinol. Metab., Oct 2005; 90: 5858 - 5865. Andrei I. Ivanov, Am J Physiol Regulatory Integrative Comp Physiol, Nov 2002; 283: 1104. Miriam Kolko, Invest. Ophthalmol. Vis. Sci., Mar 2007; 48: 1401 - 1409.