

## LYPLA1 Antibody [#C14279]

|                           |                                                                                                                                                               |
|---------------------------|---------------------------------------------------------------------------------------------------------------------------------------------------------------|
| <b>Catalog Number:</b>    | C14279                                                                                                                                                        |
| <b>Concentration:</b>     | 1mg/ml                                                                                                                                                        |
| <b>Swiss-Prot No.:</b>    | O75608                                                                                                                                                        |
| <b>Other Names:</b>       | Acyl-protein thioesterase 1; EC 3.1.2.-; Lysophospholipase 1; Lysophospholipase I; APT1, LPL1                                                                 |
| <b>Storage/Stability:</b> | Store at -20°C/1 year                                                                                                                                         |
| <b>Form of Antibody:</b>  | Rabbit IgG in phosphate buffered saline (without $Mg^{2+}$ and $Ca^{2+}$ ), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.                          |
| <b>Immunogen:</b>         | The antiserum was produced against synthesized peptide derived from internal of human LYPLA1.                                                                 |
| <b>Purification:</b>      | The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.                                         |
| <b>Specificity:</b>       | LYPLA1 antibody detects endogenous levels of total LYPLA1 protein.                                                                                            |
| <b>Reactivity:</b>        | Human, Mouse, Rat                                                                                                                                             |
| <b>Applications:</b>      | WB: 1:500~1:1000   ELISA: 1:40000                                                                                                                             |
| <b>References:</b>        | Hu G.; Submitted (AUG-1998) to the EMBL/GenBank/DDBJ databases.<br>Devedjiev Y., Structure 8:1137-1146(2000).<br>Zhang Q.-H., Genome Res. 10:1546-1560(2000). |