

## AASDHPPT Antibody [#C16447]

<b>Catalog Number:</b>	C16447
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q9NRN7
<b>Other Names:</b>	4'-phosphopantetheinyl transferase; AASD-PPT; ADPPT; alpha-aminoadipic semialdehyde dehydrogenase-phosphopantetheinyl transferase; EC 2.7.8.-; L-aminoadipate-semialdehyde dehydrogenase-phosphopantetheinyl transferase; LYS2; LYS5
<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 peptide derived from internal of human AASDHPPT.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	AASDHPPT antibody detects endogenous levels of total AASDHPPT protein.
<b>Reactivity:</b>	Human, Mouse, Rat
<b>Applications:</b>	WB: 1:500~1:1000 ELISA: 1:40000
<b>References:</b>	Praphanphoj V., Mol. Genet. Metab. 72:336-342(2001). Lai C.-H., Genome Res. 10:703-713(2000). Zhang Q.-H., Genome Res. 10:1546-1560(2000).