

## IARS2 Antibody [#C11759]

<b>Catalog Number:</b>	C11759
<b>Concentration:</b>	1mg/ml
<b>Swiss-Prot No.:</b>	Q9NSE4
<b>Other Names:</b>	EC 6.1.1.5, isoleucine-tRNA synthetase 2, mitochondrial
<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 IARS2.
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
<b>Specificity:</b>	IARS2 antibody detects endogenous levels of total IARS2 protein.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	WB: 1:500~1:1000 IHC: 1:50~1:100 IF: 1:500~1:1000 ELISA: 1:5000
<b>References:</b>	Gregory S.G., Nature 441:315-321(2006). Ota T., Nat. Genet. 36:40-45(2004). Olsen J.V., Cell 127:635-648(2006).