

## HIRA(Phospho-Thr555) Antibody [#8A1037]

<b>Catalog Number:</b>	8A1037
<b>Concentration:</b>	1mg/ml
<b>Origin of Antibody:</b>	Rabbit
<b>Swiss-Prot No.:</b>	P54198
<b>NCBI Gene Symbol:</b>	HIRA
<b>Other Names:</b>	DGCR1, TUP1 like enhancer of split protein 1, TUPLE1
<b>All Sites:</b>	Human: Thr555; Mouse: Thr554
<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 phosphopeptide derived from human HIRA around the phosphorylation site of threonine 555 (L-T-T <sub>P</sub> -P-S).
<b>Purification:</b>	The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific phosphopeptide. The antibody against non-phosphopeptide was removed by chromatography using non-phosphopeptide corresponding to the phosphorylation site.
<b>Specificity:</b>	HIRA (Phospho-Thr555) antibody detects endogenous levels of HIRA only when phosphorylated at threonine 555.
<b>Reactivity:</b>	Human, Mouse
<b>Applications:</b>	IF: 1:500~1:1000 , ELISA: 1:5000
<b>References:</b>	Lamour V., Hum. Mol. Genet. 4:791-799(1995). Lorain S., Genome Res. 6:43-50(1996). Halford S., Hum. Mol. Genet. 2:2099-2107(1993).