

Thyroid Hormone Receptor α Antibody [#C0345]

- Catalog Number:** C0345-1, C0345-2, C0345
- Amount:** 50 μ g/50 μ l, 100 μ g/100 μ l, 200 μ g/200 μ l
- Swiss-Prot No. :** P10827
- All Names:** C-erbA-alpha, EAR-7, EAR7, ERBA1, NR1A1, THA, THRA, THRA1, TR-alpha-2, Thyroid hormone receptor alpha, c-erbA-1
- Form of Antibody:** Rabbit IgG in phosphate buffered saline (without Mg²⁺ and Ca²⁺), pH 7.4, 150mM NaCl, 0.02% sodium azide and 50% glycerol.
- Storage/Stability:** Store at -20°C/1 year
- Immunogen:** The antiserum was produced against synthesized peptide derived from human thyroid hormone receptor α .
- Purification:** The antibody was affinity-purified from rabbit antiserum by affinity-chromatography using epitope-specific immunogen.
- Specificity/Sensitivity** Thyroid Hormone Receptor α antibody detects endogenous levels of total thyroid hormone receptor α protein.
- Reactivity:** Human, Mouse, Rat
- Applications:** WB: 1:500~1:1000 IHC: 1:50~1:100 ELISA: 1:10000
- References:** D E Banker, Mol. Cell. Biol., Oct 1991; 11: 5079 - 5089.
V. Krishna K. Chatterjee, PNAS, Dec 1989; 86: 9114 - 9118.
GE Muscat, Cell Growth Differ, Apr 1993; 4: 269.
Yan-Yun Liu, J. Biol. Chem., Oct 2003; 278: 38913 - 38920.