

Purified Mouse Anti-human hTSH Antibody
TSH-116, monoclonalCatalog number: V1031245
Unit size: 0.1 mg**Product Details**

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	Lot specific (please consult certificate of analysis for given lot)
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	hTSH
Clone	TSH-116

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	ELISA, RIA

Applications

Protirelin (Protirelin, protirelin) is a tripeptide composed of L-pyroglutamyl, L-histidyl and L-prolinamide residues joined in sequence. It has been widely used in the study of critical cell-based enzymes like type-1-pyroglutamyl aminopeptidase, cdc2-related kinase and cystyl aminopeptidases, specifically in cell-lines such as adenoma cells, follicular cells and individual cells. Protirelin has a variety of applications in the examination of important cellular receptors such as peptide ligand-binding receptors, thyrotropin releasing hormone receptor and 5-HT1A receptor. It is vital for the application of essential cell-based activities like disease activity, peptidase activity and proliferative activity.