

**Purified Mouse Anti-human Thyroglobulin
Antibody *2H11, monoclonal***Catalog number: V1032330
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	Thyroglobulin
Clone	2H11

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	IHC(P)

Applications

Thyroglobulin (Tg) is a 305 kDa protein that can be found in the extracellular region of cells. In humans, Tg has been associated with vital functions such as hormone activity. It recognizes identical protein. Tg aids in myelination, and furthermore, plays an important role in organismal processes, for example, signal transduction, thyroid gland development and hormone biosynthetic process. Sequencing of Tg has shown it contains 11 types of conserved structural units: Thyroglobulin type-1 1, Thyroglobulin type-1 2, Thyroglobulin type-1 3, Thyroglobulin type-1 4, Thyroglobulin type-1 5, Thyroglobulin type-1 6, Thyroglobulin type-1 7, Thyroglobulin type-1 8, Thyroglobulin type-1 9, Thyroglobulin type-1 10 and Thyroglobulin type-1 11 domain. Mutations and abnormalities in Tg have been thought to be involved with a number of diseases, namely, thyroid dysmorphogenesis 3 (TDH3) and autoimmune thyroid disease type 3 (AITD3). Thyroid dysmorphogenesis 3, an autosomal recessive inheritance disorder characterized by Intellectual disability, goiter and Increased T3/T4 ratio, has especially been of interest to scientists.