

**Biotin Mouse Anti-human Cytokeratin 19
Antibody *A53-B/A2, monoclonal***

Catalog number: V103490

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	Cytokeratin 19
Clone	A53-B/A2
Conjugate	Biotin

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with Biotin under optimal conditions
Application	IHC(P), FC (QC TESTED), IP, WB, ICC, ELISA

Applications

Keratin, type I cytoskeletal 19 is a protein with a molecular weight of 44 kDa, located in the the intermediate filament, Z disc and extracellular exosome of cells. It is sometimes referred to as keratin-19 and cytokeratin-19. In humans, cytokeratin-19 binds to protein-containing complex. Cytokeratin-19 is an integral part of organismal processes, in particular, sarcomere organization, viral process and cornification. Sequencing of cytokeratin-19 has demonstrated it contains a primary structural unit, the IF rod domain. Cytokeratin-19 is the subject of extensive application stemming from the fact that it acts in the Notch signaling pathway.