

Purified Mouse Anti-human Cytokeratin 18 Antibody *DC-10, monoclonal, Cross Adsorbed*

Catalog number: V103440

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 18
Clone	DC-10

Biological Properties

Preparation	Antibody purified by affinity chromatography, cross-adsorbed against mouse, dog, rat, pig, hamster, sheep, cow serum and then conjugated with under optimal conditions
Application	IP, WB, IHC(P), ICC, ELISA, FC

Applications

Keratin, type I cytoskeletal 18 is a 56 kDa protein that can be found in the nucleolus, cytoplasm and microtubule organizing center of cells. It is sometimes referred to as cell proliferation-inducing gene 46 protein, Keratin-18 or Cytokeratin-18. In Homo sapiens, keratin, type I cytoskeletal 18 recognizes scaffold protein, cadherin involved in cell-cell adhesion and RNA. It is an integral part of organismal processes, for example, intermediate filament cytoskeleton organization, anatomical structure morphogenesis and Golgi to plasma membrane CFTR protein transport. It is the subject of intensive study stemming from the fact that it is a member of the extrinsic apoptotic signaling pathway and tumor necrosis factor-mediated signaling pathway. Sequencing of Keratin, type I cytoskeletal 18 has shown evidence that it contains a primary structural unit, the IF rod domain. It has been closely linked to essential functions like structural molecule activity. Keratin, type I cytoskeletal 18 is a negative regulator of apoptotic process. Mutations and abnormalities in Keratin, type I cytoskeletal 18 have been associated with a number of diseases, for example, Cirrhosis (CIRRH).