

**Purified Mouse Anti-human Cytokeratin 10
Antibody *VIK-10, monoclonal***Catalog number: V103425
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 10
Clone	VIK-10

Biological Properties

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

Applications

Keratin, type I cytoskeletal 10 (sometimes referred to as Cytokeratin-10, K10, CK-10 or Keratin-10) is a protein with a molecular weight of 69 kDa, located in the cornified envelope, nucleus and cytoplasm of cells. In humans, keratin-10 has been closely linked to vital functions like protein heterodimerization activity, and furthermore, sequencing of Keratin-10 has shown it contains a primary structural unit, the IF rod domain. Moreover, Keratin-10 promotes epidermis development. Keratin-10 is thought to be essential to organismal processes, namely, protein heterotetramerization, cornification and keratinocyte differentiation. Mutations and abnormalities in Keratin-10 have been associated with a number of diseases, for instance, epidermolytic hyperkeratosis (EHK), annular epidermolytic Ichthyosis (AEI) and erythroderma, ichthyosiform, congenital reticular (CRIE). Epidermolytic hyperkeratosis, a disorder characterized by erythroderma, abnormal blistering of the skin and scaling skin, has especially been of interest to researchers.