

PE Mouse Anti-human/mouse/rabbit Acidic Cytokeratins Antibody *AE1, monoclonal*Catalog number: V103010
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, mouse, rabbit, cat, rat, dog, chicken, non-human primates
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Acidic Cytokeratins
Clone	AE1
Conjugate	PE

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with PE under optimal conditions
Application	FC (QC TESTED)

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. Sequencing of keratin-10 has shown it contains a primary structural unit, the IF rod domain. Keratin-10 promotes epidermis development. It 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.