

**Purified Mouse Anti-human/mouse/rabbit
Basic Cytokeratins Antibody *AE3,
monoclonal***

Catalog number: V103140

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, cow, non-human primates, chicken, rat
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Basic Cytokeratins
Clone	AE3

Biological Properties

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

Applications

Keratin, type II cytoskeletal 2 epidermal is a 77 kDa protein that can be found in the nucleus, cytosol and extracellular exosome of cells. It is also known as keratin-2 epidermis, type-II keratin Kb2 and keratin-2e. In Homo sapiens, type-II keratin Kb2 recognizes cytoskeletal protein. Sequencing of type-II keratin Kb2 has shown it contains a primary structural unit, the IF rod domain. Type-II keratin Kb2 plays a role in the upregulation of epidermis development. It aids in organismal processes, in particular, peptide cross-linking, intermediate filament organization and keratinocyte migration. Type-II keratin Kb2 is clinically significant because abnormalities in its function have been associated with diseases such as Ichthyosis bullosa of Siemens (IBS).