

**Biotin Mouse Anti-mammalian Cytokeratins
Antibody *C-11, monoclonal***

Catalog number: V103535

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	Mammalian
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Cytokeratins
Clone	C-11
Conjugate	Biotin

Biological Properties

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

Applications

Keratin, type II cytoskeletal 7 (sometimes referred to as CK-7, Type-II keratin Kb7, Keratin-7, Sarcolectin, K7 or Cytokeratin-7) is a protein with a molecular weight of 60 kDa, found in the keratin filament, cytosol and cytoplasm of cells. Sequencing of cytokeratin-7 has exemplified it contains a primary structural unit, the IF rod domain. It is thought to be essential to organismal processes, in particular, viral process, cornification and keratinization.