

**Purified Mouse Anti-human/pig/cow
Cytokeratin 7/17 Antibody *C-46,
monoclonal, Cross Adsorbed***

Catalog number: V103505
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, pig, cow
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Cytokeratin 7/17
Clone	C-46

Biological Properties

Preparation	Antibody purified by affinity chromatography, cross-adsorbed against mouse, rabbit, rat serum and then conjugated with under optimal conditions
Application	IP, WB, 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.