

Purified Mouse Anti-human/mouse/xenopus RPSA Antibody
RP-01, monoclonal

Catalog number: V1032080

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, xenopus, cow, chicken
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	RPSA
Clone	RP-01

Biological Properties

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

Applications

40S ribosomal protein SA is a 33 kDa protein that can be located in the plasma membrane, extracellular exosome and cytoplasm of cells. It is also known as Laminin-binding protein precursor p40, Laminin receptor 1 and Colon carcinoma laminin-binding protein. In Homo sapiens, 40S ribosomal protein SA is an integral part of organismal processes, in particular, translational initiation, ribosomal small subunit assembly and SRP-dependent cotranslational protein targeting to membrane. It has been closely linked to important functions such as virus receptor and laminin receptor activity. 40S ribosomal protein SA binds with RNA, ribosome and laminin. Mutations and abnormalities in 40S ribosomal protein SA have been thought to be involved with a number of diseases, namely, isolated congenital asplenia. Isolated congenital asplenia, a disorder characterized by abnormality of metabolism/homeostasis, Howell-Jolly bodies and thrombocytosis, has in particular been of interest to researchers.