

**Purified Mouse Anti-epstein-barr virus
EBNA-1 (EBV) Antibody *E1-2.5,
monoclonal***

Catalog number: V103610

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	Epstein-barr virus
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	EBNA-1 (EBV)
Clone	E1-2.5

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	ICC, ELISA

Applications

Epstein-Barr nuclear antigen 1 is a protein with a molecular weight of 56 kDa, expressed in the nucleus of the host cells. In human herpesvirus 4, EBNA-1 binds to DNA. It has been thought to be involved with key functions such as DNA-binding transcription factor activity. EBNA-1 aids in DNA replication, and in addition, has been found to be involved in organismal processes, for example, suppression by virus of host antigen processing and presentation, suppression by virus of host adaptive immune response and viral latency. EBNA-1 positively regulates DNA-templated transcription.