

Purified Mouse Anti-Benzo[a]pyrene Antibody *BAP-13, monoclonal*

Catalog number: V103165
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

Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Benzo[a]pyrene
Clone	BAP-13

Biological Properties

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

Applications

Benzo[a]pyrene (BaP, BP) is a crystalline, aromatic hydrocarbon consisting of five fused benzene rings and formed during the incomplete combustion of organic matter. It is vital to the research of key cell-based processes like sperm apoptosis, annual basis and mechanistic basis, especially in cell-lines such as Cl41 cells, tumor cells and IgM AFC/spleen cell. And it, furthermore, has been widely used in the examination of key biological proteins like microsomal proteins. It is essential for the study of key cellular enzymes such as flue gases, poisonous gases and signaling by receptor tyrosine kinases.