

**Culture supernatant Mouse Anti-human
Blood Group B Antibody *HEB-29,
monoclonal***

Catalog number: V103280

Unit size: 1.0 ml

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	Culture supernatant, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Blood Group B
Clone	HEB-29
Conjugate	Culture supernatant

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with Culture supernatant under optimal conditions
Application	AGG, IHC(P)

Applications

Identifies erythrocytes and glycoproteins of blood group B.