

**Purified Mouse Anti-human BG8, Lewis^y
Antibody *AHC0018, Monoclonal, IgG1,
kappa***

Catalog number: V104015
Unit size: 100 µg

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), 0.02% sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	IgG1, kappa
Immunogen	BG8, Lewis ^y
Clone	AHC0018

Biological Properties

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	IHC

Applications

Identifies BG8, Lewis y blood group antigens.