

**Purified Mouse Anti-human Sialyl Lewis a
Antibody *121SLE, monoclonal***Catalog number: V1032110
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	Tris buffered saline (TBS) solution, 15 mM sodium azide, 0.2% (w/v) BSA

Antibody Properties

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Immunogen	Sialyl Lewis a
Clone	121SLE

Biological Properties

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

Applications

Identifies Sialyl Lewis X, also known as CD15s.