

**PE/Texas Red® Anti-human CD19 Antibody**  
**\*HI19a\***Catalog number: 101901S0, 101901S1, 101901S2  
Unit size: 25 tests, 100 tests, 500 tests**Product Details**

---

Storage Conditions	2-8°C with minimized light exposure. Do not freeze.
Expiration Date	12 months upon receiving
Concentration	0.1 mg/mL
Formulation	Phosphate-buffered saline (PBS, pH 7.2), 0.09% sodium azide, 0.2% (w/v) BSA

**Antibody Properties**

---

Species Reactivity	Human
Class	Primary
Clonality	Monoclonal
Host	Mouse
Isotype	Mouse IgG1
Immunogen	CD19 (B4)
Clone	HI19a
Conjugate	PE/Texas Red®

**Biological Properties**

---

Preparation	Antibody purified by affinity chromatography and then conjugated with PE/Texas Red® under optimal conditions
Application	Flow Cytometry (FACS)

**Spectral Properties**

---

Conjugate	PE/Texas Red®
Excitation Wavelength	567 nm
Emission Wavelength	615 nm

**Applications**

---

HI19a is an anti-human monoclonal antibody that targets the CD19 antigen. CD19 (also known as T-cell surface antigen Leu-12 or B-lymphocyte surface antigen B4) is a 95 kD transmembrane glycoprotein that is found on the surface of cells like B cells and stem cells. CD19 plays a role in critical cellular pathways, in particular, the antigen receptor-mediated signaling pathway and B cell receptor signaling pathway. Furthermore, in certain organisms, it plays a role in the upregulation of phosphatidylinositol 3-kinase activity, positively regulates release of sequestered calcium ion into cytosol and plays a role in the upregulation of protein kinase B signaling. From a research standpoint, it is of biological interest due to its

association with vital macromolecules/ligands such as CD81 and Leu-13. CD19 is a very popular antibody target, with over 36000 publications in the last decade. CD19 is frequently used in flow cytometry applications as a phenotypic marker for differentiation of cell types, specifically in the study of immunology and costimulatory molecules. This antibody was purified through affinity chromatography and conjugated to PE/Texas Red® (ex/em = 567/615 nm). It is compatible with the 561 nm laser and 610/20 nm bandpass filter (for example, as in the BD LSRFortessa™ X-20).