

**Purified Chicken Anti-Digoxigenin Antibody**  
**\*Polyclonal\***Catalog number: V102065  
Unit size: 1.0 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	, 0.2% (w/v) BSA

**Antibody Properties**

Class	Primary
Clonality	Polyclonal
Host	Chicken
Immunogen	Digoxigenin

**Biological Properties**

Preparation	Antibody purified by affinity chromatography and then conjugated with under optimal conditions
Application	Flow Cytometry (FACS)

**Applications**

Digoxigenin is a metabolite of digoxin(1,2). It has a variety of applications in the study of essential biological pathways such as TNF receptor signaling pathway, FAS (CD95) signaling pathway and versatile pathway, particularly in cell-lines such as multinucleated odontoclasts, structural cell and investigations using hepatocytes. Digoxigenin, also, is vital to the application of critical biological receptors like adenosine receptor, androgen receptor and CX3CL1 receptor, and it has a variety of applications in the research of critical cellular diseases such as demyelinating disease, preventing disease and genetic disease.