

---

## LSR Polyclonal Antibody, Biotin Conjugated

(Catalog #A69018)

---

### Background

Probable role in the clearance of triglyceride-rich lipoprotein from blood. Binds chylomicrons, LDL and VLDL in presence of free fatty acids and allows their subsequent uptake in the cells.

### Description

LSR Polyclonal Antibody, Biotin Conjugated. Biotin. Raised in: Rabbit.

### Formulation

Liquid. 0.03% Proclin 300, 50% Glycerol, 0.01M PBS, pH 7.4.

### Specificity

Human

### Isotype

IgG

### Uniprot ID

Q86X29

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human Lipolysis-stimulated lipoprotein receptor protein (112-233AA)

### Storage

Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

Lipolysis-stimulated lipoprotein receptor, LSR, LISCH

### Application

ELISA; Recommended dilution: ELISA 1:500 - 1:1000

---