

## ABCA12 Polyclonal Antibody

(Catalog # A69003)

### Background

Probable transporter involved in lipid homeostasis.

### Description

ABCA12 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

### Formulation

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

### Specificity

Human

### Isotype

IgG

### Uniprot ID

Q86UK0

### Purification

>95%, Protein G purified

### Immunogen

Recombinant Human ATP-binding cassette sub-family A member 12 protein (45-158AA)

### Storage

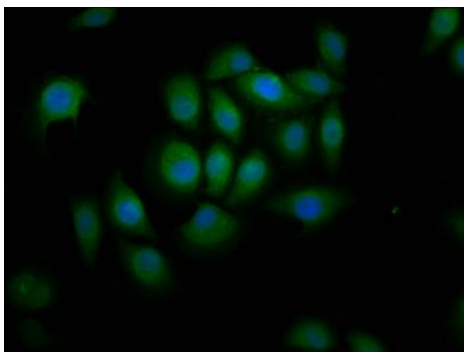
Shipped at 4°C. Upon receipt, store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

### Alternative Names

ATP-binding cassette sub-family A member 12, ATP-binding cassette transporter 12, ATP-binding cassette 12, ABCA12, ABC12

### Application

ELISA, IF; Recommended dilution: IF: 1:50-1:200



Immunofluorescence staining of A549 cells with ABCA12 Polyclonal Antibody at 1:166, counter-stained with DAPI. The cells were fixed in 4% formaldehyde, permeabilized using 0.2% Triton X-100 and blocked in 10% normal Goat Serum. The cells were then incubated with the antibody overnight at 4°C. The secondary antibody was Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L).