

CYSLTR1 Polyclonal Antibody

(Catalog # A62370)

Background

Receptor for cysteinyl leukotrienes mediating bronchoconstriction of individuals with and without asthma. Stimulation by LTD4 results in the contraction and proliferation of smooth muscle, edema, eosinophil migration and damage to the mucus layer in the lung. This response is mediated via a G-protein that activates a phosphatidylinositol-calcium second messenger system. The rank order of affinities for the leukotrienes is LTD4 >> LTE4 = LTC4 >> LTB4.

Description

CYSLTR1 Polyclonal Antibody. Unconjugated. Raised in: Rabbit.

Formulation

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

Specificity

Human, Mouse

Isotype

IgG

Uniprot ID

Q9Y271

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Cysteinyl leukotriene receptor 1 protein (298-337AA)

Storage

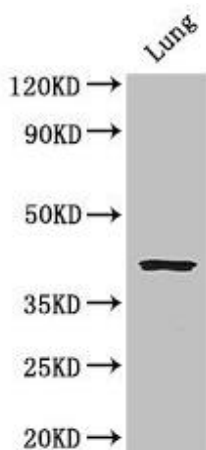
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

Cysteinyl leukotriene receptor 1, CysLTR1, Cysteinyl leukotriene D4 receptor, LTD4 receptor, G-protein coupled receptor HG55, HMTMF81, CYSLTR1, CYSLT1

Application

ELISA, WB, IHC, IF; Recommended dilution: WB:1:1000-1:5000, IHC:1:20-1:200, IF:1:50-1:200



Western Blot

Positive WB detected in: Mouse lung tissue

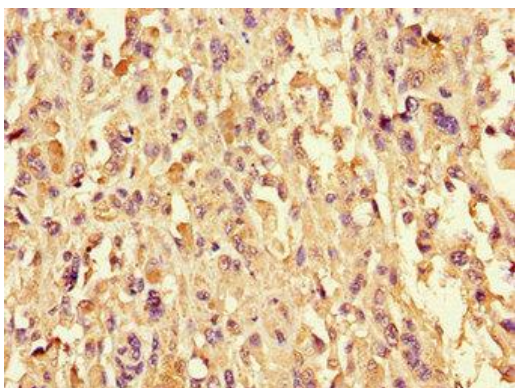
All lanes: CYSLTR1 Polyclonal Antibody at 3ug/ml

Secondary

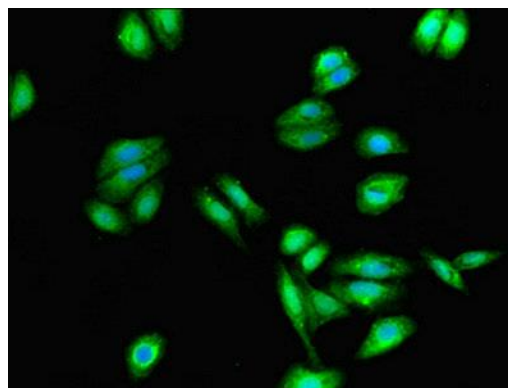
Goat polyclonal to rabbit IgG at 1/50000 dilution

Predicted band size: 39 kDa

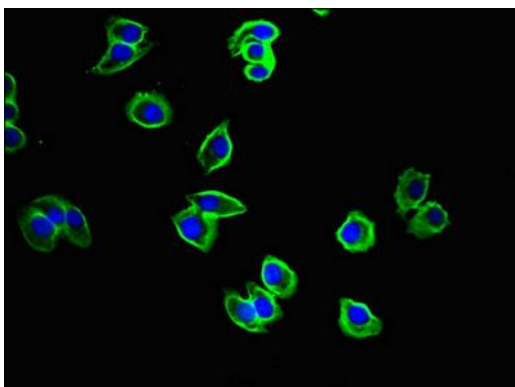
Observed band size: 39 kDa



Immunohistochemistry of paraffin-embedded human melanoma tissue using CYSLTR1 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of HepG2 cells using CYSLTR1 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)



Immunofluorescent analysis of A375 cells using CYSLTR1 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)