

LSMEM1 Polyclonal Antibody

(Catalog # A66202)

Background

cytoplasm

Description

LSMEM1 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

Q8N8F7

Purification

>95%, Protein G purified

Immunogen

Recombinant Human Leucine-rich single-pass membrane protein 1 protein (1-131AA)

Storage

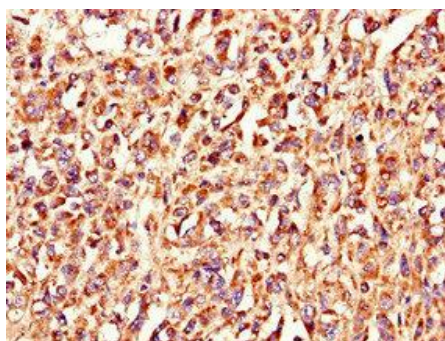
Shipped at 4°C. Upon delivery aliquot and store at -20°C (short-term) or -80°C (long-term). Avoid repeated freeze.

Alternative Names

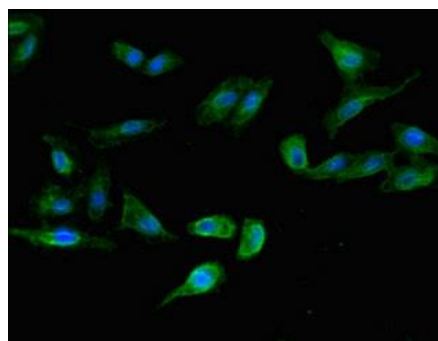
Leucine-rich single-pass membrane protein 1, LSMEM1, C7orf53

Application

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



Immunohistochemistry of paraffin-embedded human melanoma cancer using LSMEM1 Polyclonal Antibody at dilution of 1:100



Immunofluorescent analysis of HeLa cells using LSMEM1 Polyclonal Antibody at a dilution of 1:100 and Alexa Fluor 488-conjugated AffiniPure Goat Anti-Rabbit IgG (H+L)