ERGIC-53 monoclonal antibody (OTI1A8)

The ER-Golgi intermediate compartment (ERGIC) marker ERGIC-53 is a mannose-specific membrane lectin operating as a cargo receptor for the transport of glycoproteins from the ER to the ERGIC. ERGIC-53 can also be used for studying aspects of protein trafficking in the secretory pathway.

This antibody is covered by our Worry-Free Guarantee.

Citations: 19

View Online »

Ordering Information

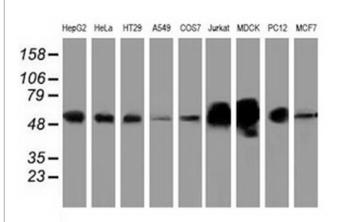
Order Online »

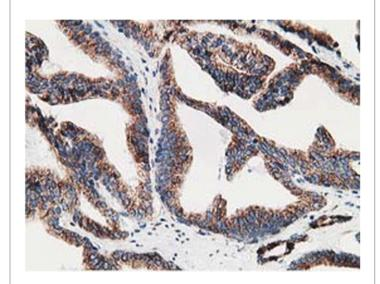
ENZ-ABS300-0100

100µl

Manuals, SDS & CofA

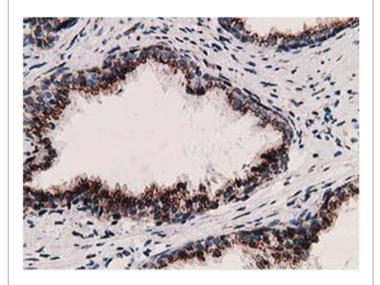
View Online »

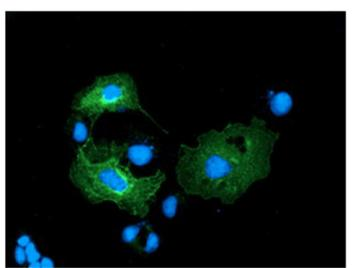




Western blot analysis of extracts (35µg) from 9 different cell lines by using ERGIC-53 monoclonal antibody (OTI1A8).

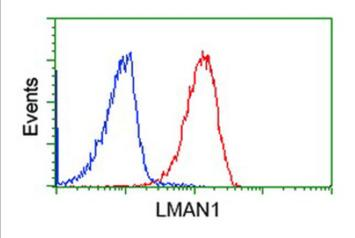
Immunohistochemistry staining of paraffin-embedded carcinoma of human prostate tissue using ERGIC-53 monoclonal antibody (OTI1A8).



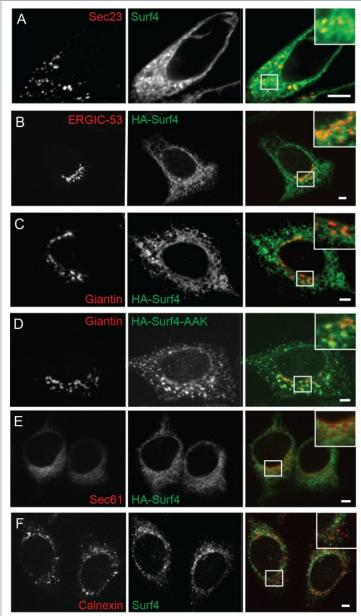


Immunohistochemistry staining of paraffin-embedded human prostate tissue within the normal limits using ERGIC-53 monoclonal antibody (OTI1A8).

ERGIC-53 monoclonal antibody (OTI1A8) immunofluorescent staining of COS7 cells transiently transfected by pCMV6-ENTRY ERGIC-53.



Flow cytometric analysis of Jurkat cells, using ERGIC-53 monoclonal antibody (OTI1A8) (red), compared to a nonspecific negative control antibody (blue).



Bar: 5um

Immunofluorescence microscopy of HEK293A cells shows Surf4 accumulates in and around ERESs.(A) Fluorescent signal for endogenous Surf4 (green) was strongest at punctate structures positive for ERES marker Sec23 (red). Note additional Surf4 fluorescence in weblike structures surrounding ERES. (B) HA-Surf4 signal (green) was observed within the ERGIC (ERGIC-53, red). (C) HA-Surf4 (green) showed only low levels of colocalization with cis-Golgi marker, giantin (red). (D) Mutation of proposed COPI recycling motif by replacement of two of three near-carboxy-terminal lysines to alanines (HA-Surf4-AAK, green) increased colocalization with cis-Golgi marker, giantin (red). (E) Newly synthesized HA-Surf4 (green) was found at low levels in the rER (Sec61 marker, red). (F) Surf4 (green) did not colocalize with chaperone, calnexin (red), in the quality control domain. HEK293A cells were transfected with wild-type HA-tagged Surf4 plasmid (B, C, E), or Surf4KO HEK293A cells were transfected with carboxylterminal di-lysine mutation, HA-Surf4-AAK (D), 18 hr prior to fixation. Bars = 5µm. The cells in each panel are shown 3 times, first with organelle marker, then Surf4

Handling & Storage

Handling Avoid freeze/thaw cycles.

Long Term Storage -20°C

Shipping Blue Ice

Regulatory Status RUO - Research Use Only

Product Details

Alternative Name ER-Golgi intermediate compartment 53kDa protein,

LMAN1

Application Flow Cytometry, IF, IHC, WB

Clone OTI1A8

Formulation Liquid. In PBS, pH 7.3, containing 1% BSA, 50% glycerol

and 0.02% sodium azide.

Host Mouse

Immunogen Full length recombinant human ERGIC-53.

lsotype lgG1

Purity Detail Affinity purified.

Recommendation Dilutions/Conditions Flow Cytometry (1:100)Immunofluorescence

(1:100)Immunohistochemistry (1:150)Western Blot (1:200-1:500)Suggested dilutions/conditions may not be available for all applications.Optimal conditions must be determined

individually for each application.

Species Reactivity Dog, Human, Monkey, Rat

UniProt ID P49257

Worry-free Guarantee This antibody is covered by our Worry-Free Guarantee

•

Last modified: May 29, 2024



info-

eu@enzolifesciences.com