

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 »](#)

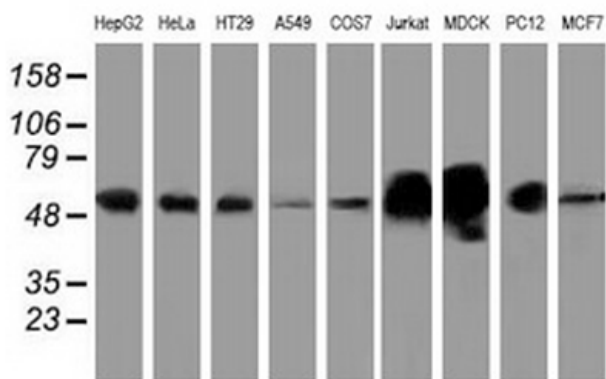
Ordering Information

[Order Online »](#)

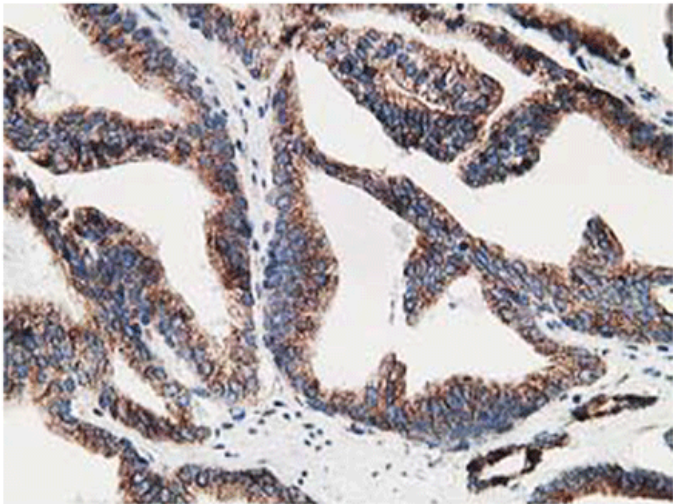
ENZ-ABS300-0100	100µl
-----------------	-------

Manuals, SDS & CofA

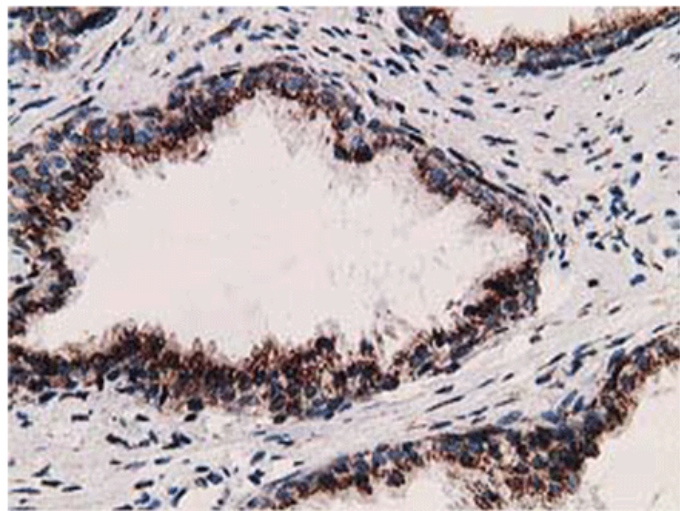
[View Online »](#)



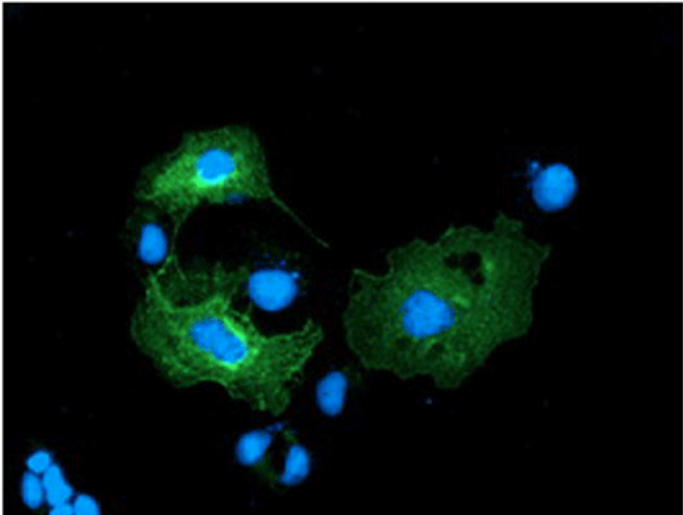
Western blot analysis of extracts (35µg) from 9 different cell lines by using ERGIC-53 monoclonal antibody (OTI1A8).



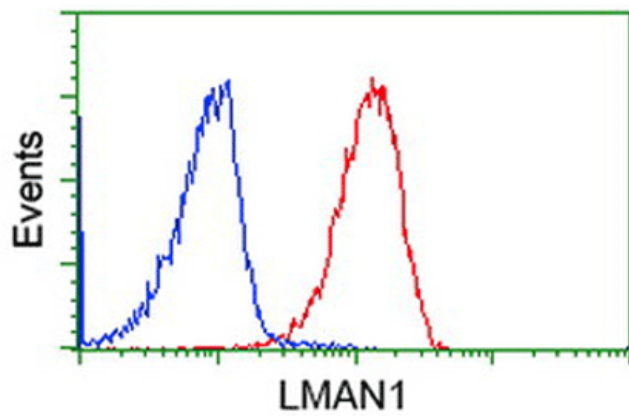
Immunohistochemistry staining of paraffin-embedded carcinoma of human prostate tissue using ERGIC-53 monoclonal antibody (OTI1A8).



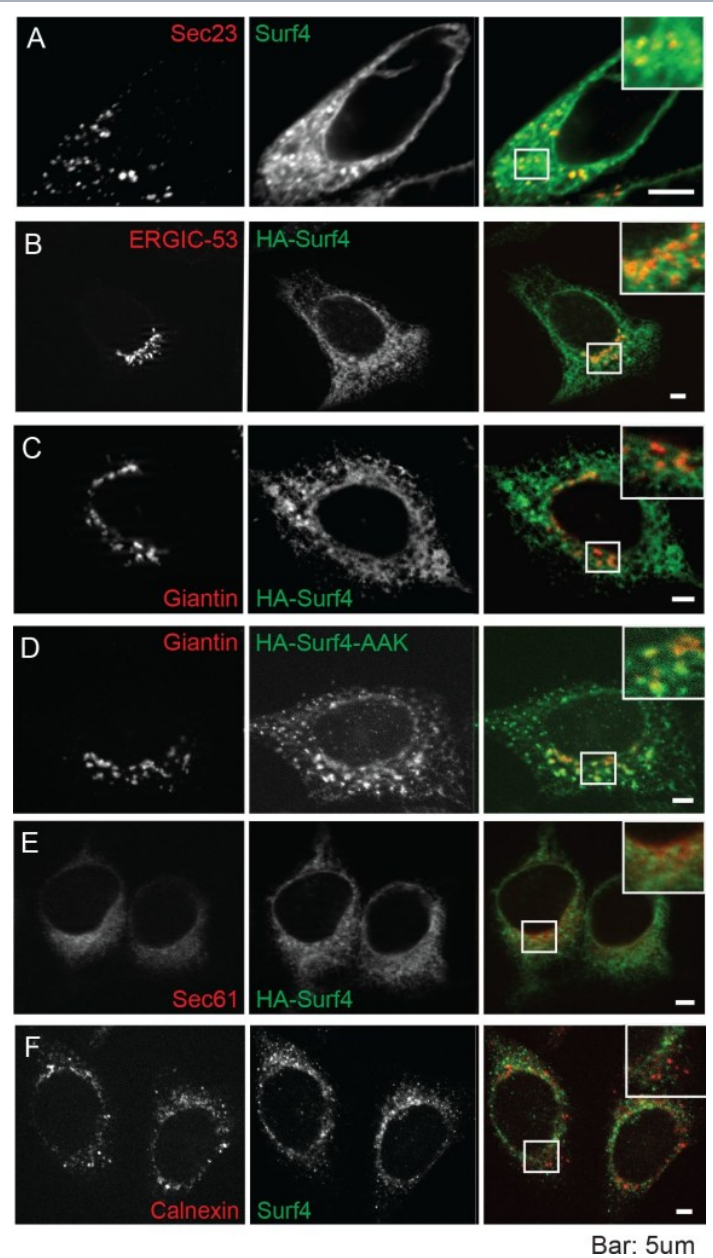
Immunohistochemistry staining of paraffin-embedded human prostate tissue within the normal limits using ERGIC-53 monoclonal antibody (OTI1A8).



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).



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 carboxyl-terminal 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	Avoid freeze/thaw cycles.
Long Term Storage	-20°C
Shipping	Blue Ice

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.
Isotype	IgG1
Purity Detail	Affinity purified.

Species Reactivity Dog, Human, Monkey, Rat

UniProt ID P49257

Worry-free Guarantee This antibody is covered by our [Worry-Free Guarantee](#)



ENZO LIFE SCIENCES,
INC.
Phone: 800.942.0430
[info-
usa@enzolifesciences.com](mailto:info-usa@enzolifesciences.com)

European Sales Office
ENZO LIFE SCIENCES
(ELS) AG
Phone: +41 61 926 8989
[info-
eu@enzolifesciences.com](mailto:info-eu@enzolifesciences.com)

Belgium, The Netherlands
& Luxembourg
Phone: +32 3 466 0420
[info-
be@enzolifesciences.com](mailto:info-be@enzolifesciences.com)

France
Phone: +33 472 440 655
[info-
fr@enzolifesciences.com](mailto:info-fr@enzolifesciences.com)

Germany
Phone: +49 7621 5500 526
[info-
de@enzolifesciences.com](mailto:info-de@enzolifesciences.com)

UK & Ireland
Phone (UK customers):
0845 601 1488
Phone: +44 1392 825900
[info-
uk@enzolifesciences.com](mailto:info-uk@enzolifesciences.com)