



Qty: 100 µg/200 µL

Mouse anti-TRA-1-60

Catalog No. 41-1000

Lot No.

Mouse anti-TRA-1-60

FORM

This monoclonal antibody is supplied as a 200 µL aliquot at a concentration of 0.5 mg/mL in PBS, pH 7.4, containing 0.1% sodium azide. This antibody is highly purified from mouse ascites by protein A chromatography.

CLONE: cl.A

ISOTYPE: Mouse IgM

IMMUNOGEN

2102Ep cl.2A6 human embryonal carcinoma cells

SPECIFICITY

This antibody is specific for the TRA-1-60 antigen. On Western blots, it identifies the target band at ~200-400 kDa.⁽¹⁾

REACTIVITY

Reactivity has been confirmed with human EC1003.7 and NTERA-2⁽⁴⁾ cell lysates by Western blotting, with frozen normal human mammary ducts, stomach, small and large intestine, endometrial and cervical glands, bronchial epithelium, bile ducts, and distal tubules of the kidney by indirect immunofluorescence⁽¹⁾, and with lysates from ¹²⁵I surface-labeled 2102Ep cl.2A6 human EC cells by immunoprecipitation⁽¹⁾.

Sample	Immuno-precipitation (native)	Immuno-fluorescence	Western Blotting
Human	+++ ⁽¹⁾	+++ ⁽¹⁾	+++

(Excellent +++, Good++, Poor +, No reactivity 0, Not applicable N/A, Not Determined ND)

USAGE

Working concentrations for specific applications should be determined by the investigator. Appropriate concentrations will be affected by several factors, including secondary antibody affinity, antigen concentration, sensitivity of detection method, temperature and length of incubations, etc. The suitability of this antibody for applications other than those listed below has not been determined. The following concentration ranges are recommended starting points for this product.

Immunoprecipitation⁽¹⁾: 2-3 µg/mL

Immunofluorescence⁽¹⁾: 5-10 µg/mL

Western Blotting: 1-5 µg/mL

STORAGE

Store at 2-8°C for up to one month. Store at -20°C for long-term storage. Avoid repeated freezing and thawing.

(cont'd)

BACKGROUND

The TRA-1-60 antigen is expressed by human embryonal carcinoma (EC) cells and is associated with a polydisperse glycoprotein with a molecular weight in excess of 200,000.^{1,3} TRA-1-60 recognizes an antigen of similar size as detected by TRA-1-81, GCTM2, K4, and K21, and all these antibodies detect similar but distinct patterns of expression on panels of tumor-derived cell lines.² TRA-1-60 and TRA-1-81 epitopes are associated with distinct molecular species. TRA-1-60, TRA-1-81, GCTM2, and K21 antibodies all recognize distinct epitopes formed by modifications of a common keratan sulfate core molecule that is characteristically expressed by human EC cells.⁴ TRA-1-60, clone cl.A, likely recognizes the carbohydrate part of the molecule as TRA-1-60 activity is destroyed by neuraminidase.⁴

REFERENCES

1. Andrews PW, et al. *Hybridoma* 3(4):347-361, 1984.
2. Andrews PW, et al. *Int J Cancer* 66(6):806-816, 1996.
3. Andrews PW, et al. *Rec Res Cancer Res* 123: 63-83, 1991.
4. Badcock G, et al. *Cancer Research* 59:4715-4719, 1999.

RELATED PRODUCTS

Product	Conjugate	Cat. No.
Protein A	Sepharose [®] 4B	10-1041
rec-Protein G	Sepharose [®] 4B	10-1241

Conjugate	ZyMAX[™] Goat x Rabbit IgG (H+L)	ZyMAX[™] Goat x Mouse IgG (H+L)
Purified	81-6100	81-6500
FITC	81-6111	81-6511
TRITC	81-6114	81-6514
Cy [™] 3	81-6115	81-6515
Cy [™] 5	81-6116	81-6516
HRP	81-6120	81-6520
AP	81-6122	81-6522
Biotin	81-6140	81-6540

Zymed[®] and ZyMAX[™] are trademarks of Zymed Laboratories Inc. Cy[™] and SEPHAROSE[®] are trademarks of GE Healthcare.

For Research Use Only

www.invitrogen.com

Invitrogen Corporation • 542 Flynn Rd • Camarillo • CA 93012 • Tel: 800.955.6288 • E-mail: techsupport@invitrogen.com

PI411000

(Rev 08/09) DCC-08-1089

Important Licensing Information - These products may be covered by one or more Limited Use Label Licenses (see the Invitrogen Catalog or our website, www.invitrogen.com). By use of these products you accept the terms and conditions of all applicable Limited Use Label Licenses. Unless otherwise indicated, these products are for research use only and are not intended for human or animal diagnostic, therapeutic or commercial use.