



Qty: 100 µg/400 µL

Rabbit anti-Kinesin

Catalog No. 40-8100

Lot No.

Rabbit anti-Kinesin

FORM

This polyclonal antibody is supplied as a 400 µL aliquot at a concentration of 0.25 mg/mL in phosphate buffered saline (pH 7.4) containing 0.1% sodium azide. This antibody is epitope-affinity purified from rabbit antiserum.

PAD: ZMD.502

IMMUNOGEN

Bacterially expressed cDNA encoding full-length human neuronal kinesin heavy chain, which shares 96% homology with mouse

SPECIFICITY

This antibody is specific for the FKHC3 coiled-coil stalk domain of kinesin protein. On Western blots, it identifies the target band at ~120-130 kDa.

REACTIVITY

Reactivity has been confirmed with pig and mouse brain homogenates and HEK 293T lysates by Western blotting, and with mouse 3T3 cells by immunocytochemistry.

| Sample | Western Blotting | Immunocytochemistry |
|-----------|------------------|---------------------|
| Human | +++ | ND |
| Mouse | +++ | +++ |
| Pig | +++ | ND |
| Immunogen | N/A | N/A |

(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.

Western Blotting: 1–3 µg/mL

Immunocytochemistry: 0.5 – 1.0 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)

www.invitrogen.com

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

PI408100

(Rev 10/08) 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.

BACKGROUND

Active transport in cells is driven by three distinct classes of molecular motors: the kinesin family that moves toward the plus-end of microtubules, the dynein family that moves toward the minus-end of microtubules, and myosin motors that move along actin filaments¹.

Kinesins are microtubule-based motor proteins involved in the transport of organelles in eukaryotic cells. They typically consist of 2 identical, ~110-120 kDa heavy chains and 2 identical, ~60-70 kDa light chains. The heavy chain contains 3 domains: a globular N-terminal motor domain, which converts the chemical energy of ATP into a motile force along microtubules in one fixed direction, a central alpha-helical rod domain (stalk domain) which enables the 2 heavy chains to dimerize, and a globular C-domain which interacts with light chains and organelle receptors²⁻³. Kinesin proteins operate as a dimer, traveling in a coordinated manner along microtubule protofilaments⁴⁻⁵. Human kinesins can be classified into three groups based on the position of their motor domains: N-terminal, C-terminal, and internal. There are three known kinesin family members denoted as A, B and C. Kinesins 5A and 5C are exclusively neuronal, whereas 5B appears to be ubiquitous in expression.

REFERENCES

1. Mallik R, et al. *Curr Biol* 14(22):R971-982, 2004.
2. Pfister K, et al. *J Cell Biol* 108(4):1453-1463, 1989.
3. Skoufias DA, et al. *J Biol Chem* 269(2):1477-1485, 1994.
4. Kull FJ, et al. *Essays Biochem* 35:61-73, 2000.
5. Diefenbach RJ, et al. *Biochem Biophys Res Commun* 319(3):987-992, 2004.

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 Amersham Biosciences Ltd.

For Research Use Only

AC311005

www.invitrogen.com

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

PI408100

(Rev 10/08) 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.