



Qty: 100 µg/400 µL

Rabbit anti-mGluR4

Catalog No. 51-3100

Lot No. See product label

Rabbit anti-mGluR4

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. The antibody is epitope-affinity-purified from rabbit antiserum.

POLYCLONAL ANTIBODY DESIGNATION (PAD): ZTS4

ISOTYPE: Rabbit Ig

IMMUNOGEN

Synthetic peptide derived from the C-terminal 200 amino acids of the rat mGluR4 protein.

SPECIFICITY

This antibody is specific for the mGluR4a splice variant ($M_r=93,000-110,00$). Reactivity with other related proteins has not been detected.

REACTIVITY

This antibody reacts with human, mouse, and rat mGluR4. Reactivity was confirmed on Western blots using rat brain cell lysate.

Sample	ELISA	Immuno-histochemistry (paraffin)	Western blotting
Human		nd	+
Mouse		nd	+
Rat		+	+
Immunogen	+		

nd: not determined

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 following concentration ranges are recommended starting points for this product.

ELISA: 0.1-1.0 µg/ml

Western Blotting: 1-2 µg/ml

Immunohistochemistry (paraffin): 5-10 µg/ml

The suitability of this antibody for applications other than those listed here has not been determined.

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

PI513100

(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

The metabotropic glutamate receptors mGluRs can be placed into three groups based on sequence homology, coupling to second messengers and pharmacological profiles. Group III includes mGluR4 and its two splice variants mGluR4a and mGluR4b. mGluR4a is the principal mGluR4 receptor found in the cerebellum, hippocampus and other regions of the brain. mGluR4 is also believed to be a taste receptor for the transduction of mono-sodium glutamate taste.

GENERAL REFERENCES

1. Bradley, S.R. et. al. (1996) J. Neurosci. 16; 2044-2056.
2. Nakanishi, S. (1992) Science 258; 597-603.
3. Tanabe, Y., et. al. Neuron (1992) 8; 169-179.
4. Chaudhari, N., and Roper, S.D. (1998) Annals NY Acad. Sci. 855; 398-406.

RELATED PRODUCTS**Primary Antibodies**

Product	Clone/PAD	Cat. No.
Rb x NMDAR1	ZNR2	51-3600
Ms x NMDAR1	54.1	32-0500
Ms x NMDAR 2A	A3-2D10	32-0600
Ms x NMDAR 2B	B3-13B11	32-0700
Rb x NMDAR 2B	ZK11	71-8600
Ms x GluR1+2+3	2D8	32-0100
Ms x GluR2	6C4	32-0300
Ms x GluR2+4	3A11	32-0200
Ms xGluR3	3B3	32-0400
Ms x Glutamate Transporter EAAC1	35-A9	32-1000

Protein A-Sepharose® 4B 10-1041

rec-Protein G-Sepharose® 4B 10-1241

Secondary Antibodies

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™3 and Cy™5 are trademarks of Amersham Life Sciences, Inc. Sepharose® is a registered trademark of Pharmacia LKB.

For Research Use Only

www.invitrogen.com

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

PI513100

(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.