



Catalog Number	Product Name	Quantity	Clonality, clone (isotype)	Reactive species	Applications	Reg. Status
434000	Mouse anti-human RXR β	100 μ g	mAb clone H7341 (Ms IgG2a)	Hu	WB, ELISA, IP	RUO

Mouse Anti-Human Retinoid X Receptor β

Description Retinoid X receptor β (RXR β ; NR2B2) is a member of the orphan nuclear receptor family. 9-cis retinoic acid can bind to RXR. RXR β is expressed widespread in most tissues of the embryo, including the central nervous system. Along with other members of the RXR family, RXR β plays roles in a variety of processes including embryonic patterning and organogenesis, cell proliferation and differentiation. RXRs commonly function as heterodimers with other members of the nuclear receptor superfamily.

Nomenclature NR2B2

Genbank M84820

Origin Produced in BALB/c mouse ascites after inoculation with hybridoma of mouse myeloma cells (NS-1) and spleen cells derived from a BALB/c mouse immunized with Baculovirus-expressed recombinant human RXR β (9-187 aa).

Specificity This antibody specifically recognizes human RXR β but does not recognize human RXR α and γ . Not yet tested in other species.

Purification Ammonium sulfate fractionation

Formulation Concentration is 1 mg/mL in physiological saline with 0.1% sodium azide as a preservative.

Application	Recommended Concentration*
Western Blot	1 μ g/mL
Non reducing Western Blot	1 μ g/mL
ELISA	1 μ g/mL
Immunoprecipitation	Determine by use
Electrophoretic Mobility Shift Assay	Not tested
Chromatin Immunoprecipitation	Not tested
Immunohistochemistry	Not tested

*In order to obtain the best results, optimal working dilutions should be determined by each individual user.

Storage Store at 2 - 8°C up to one month. For long-term storage, the solution may be frozen at -20°C in working aliquots. Repeated freezing and thawing is not recommended. Storage in a frost-free freezer is not recommended.

Reference Nakamura T, et al., *Nat Genet* 36(5): 528-33, 2004.

www.invitrogen.com

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

PI434000

(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.

Related Products

<i>Product</i>	<i>Conjugate</i>	<i>Cat. No.</i>
Protein A	Sepharose 4B	10-1041
rec-Protein G	Sepharose 4B	10-1241
ZyMAX™ Goat anti-rabbit IgG	Unconjugated	81-6100
ZyMAX™ Goat anti-mouse IgG	Unconjugated	81-6500

Secondary Antibody Conjugates

<i>Conjugate</i>	<i>Goat anti-rabbit IgG (H+L)</i>	<i>Goat anti-mouse IgG (H+L)</i>	<i>Ex/Em*</i>	<i>Fluorescence similar to--</i>
Alexa Fluor® 488	A11008	A11001	495/519	FITC
Alexa Fluor® 555	A21428	A21422	555/565	Cy3
Alexa Fluor® 594	A11012	A11005	590/617	Texas Red®
Alexa Fluor® 647	A21244	A21235	650/668	Cy5
HRP	81-6120	81-6520	NA**	NA
AP	81-6122	81-6522	NA	NA
Biotin	B2770	B2763	NA	NA

*Excitation/emission (nm); **Not applicable

For additional secondary antibody conjugates, visit www.invitrogen.com/antibodies

For Research Use Only

Manufactured by:
Perseus Proteomics, Inc.
4-7-6, Komaba, Meguro-ku
Tokyo 153-0041 Japan
Tel: +81-3-5738-1705
Fax: +81-3-3481-5760



www.invitrogen.com

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

PI434000

(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.