

**Mouse (monoclonal)
Anti-Human IGF-1R
(β -Subunit)**

PRODUCT ANALYSIS SHEET

Catalog Number:	AHR0361
Lot Number:	See product label
Quantity/Volume:	100 μ g/0.5 mL
Clone Number:	1-2
Isotype:	Mouse IgG2b kappa
Form of Antibody:	Purified immunoglobulin in phosphate buffered saline, pH 7.4, with 0.2% bovine serum albumin.
Preservation:	0.09% sodium azide (Caution: sodium azide is a poisonous and hazardous substance. Handle with care and dispose of properly.)
Purification:	Purified from ascites by Protein A affinity chromatography.
Immunogen:	A synthetic peptide corresponding to amino acid residues 1323-1337 (YRKNERALPLQSSTC) from the C-terminal region of the β -subunit of human IGF-1R.
Myeloma/Fusion Partners:	Produced by fusion between BALB/c mouse splenocytes and mouse myeloma NSO/1 cells.
Specificity:	<p>This monoclonal antibody recognizes a protein of $M_r=95$ kDa, identified as the β-subunit of type 1 insulin-like growth factor receptor (IGF-1R). As with the receptor for insulin, IGF-1R is synthesized as a single polypeptide which is glycosylated and proteolytically cleaved to yield a receptor composed of two α- and two β-subunits arranged in the following configuration: β-α-α-β. The α-subunits are entirely extracellular. The β-subunits each possess an extracellular domain, a single transmembrane domain, and a cytoplasmic tyrosine kinase domain which bears homology with other tyrosine kinases.</p> <p>The epitope recognized by this antibody is localized between amino acid residues 1323-1337 of the C-terminal region of the β-subunit. This antibody shows no cross-reactivity with insulin receptor (IR).</p>
Species Reactivity:	Human and rat. Other species were not tested.
Applications:	This antibody is suitable for use in Western blot analysis. Preparations from which the BSA has been omitted have been used successfully in the affinity purification of IGF-1R and IGF-1R-containing fusion proteins.
Suggested Working Dilutions:	For Western blotting, the recommended concentration is 1.0 μ g/mL with an incubation of 2 hours at room temperature. The optimal antibody concentration should be determined for each specific application.
Recommended Positive Control:	Placenta or breast carcinoma.

This product is for research use only. Not for use in diagnostic procedures.

www.invitrogen.com

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

PI AHR0361

(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.

Storage:	Store at 2-8°C. For long term storage, apportion into working aliquots and store at -20°C. Avoid repeated freeze-thaw cycles to prevent denaturing the antibody.
Expiration Date:	Expires one year from date of receipt when stored as instructed.
References:	<p>Soos, M.A., C.E. Field, R. Lammers, A. Ullrich, B. Zhang, R.A. Roth, A.S. Andersen, T. Kjeldsen, and K. Siddle (1992) A panel of monoclonal antibodies for the type 1 insulin-like growth factor receptor. Epitope mapping, effects on ligand binding, and biological activity. <i>J. Biol. Chem.</i> 267:12955-12963.</p> <p>Soos, M.A., C.E. Field, and K. Siddle (1993) Purified hybrid insulin/insulin-like growth factor-1 receptors bind insulin-like growth factor-1, but not insulin, with high affinity. <i>Biochem. J.</i> 290:419-426.</p> <p>Soos, M.A., B.T. Nave, and K. Siddle (1993) Immunological studies of type 1 IGF receptors and insulin receptors: characterization of hybrid and atypical receptor subtypes. <i>Adv. Exp. Med. Biol.</i> 343:145-157.</p>

This product is for research use only. Not for use in diagnostic procedures.

www.invitrogen.com

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

PI AHR0361

(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.