



# Mouse (monoclonal) Anti-IRS-2 Unconjugated

## PRODUCT ANALYSIS SHEET

<b>Catalog Number:</b>	AHO1532
<b>Lot Number:</b>	See product label
<b>Quantity/Volume:</b>	100 µg/0.2 mL
<b>Clone Number:</b>	8S15
<b>Isotype:</b>	IgG1 κ (mouse)
<b>Form of Antibody:</b>	Purified immunoglobulin in phosphate buffered saline, pH 7.4.
<b>Preservation:</b>	0.1% sodium azide (Caution: sodium azide is a poisonous and hazardous substance. Handle with care and dispose of properly.)
<b>Purification:</b>	Purified from ascites by affinity chromatography.
<b>Immunogen:</b>	Recombinant human IRS-2 N-terminal fragment.
<b>Specificity:</b>	This antibody recognizes an 180 kDa protein identified as insulin receptor substrate 2 (IRS-2). Insulin exerts its effects by binding to the insulin receptor, disulfide-linked heterotetrameric protein comprised of two α subunits and two β subunits. The α subunits each contain insulin binding sites and are entirely extracellular in localization. The β subunits each possess an extracellular domain, a single transmembrane domain, and a cytoplasmic tyrosine kinase domain. Binding of insulin to the α subunits induces a conformation change in the receptor which activates the kinase domain, stimulating tyrosine autophosphorylation of the receptor and tyrosine phosphorylation of at least five different insulin receptor substrates designated IRS-1-4, and Shc. This tyrosine phosphorylation produces docking sites for proteins bearing SH2 domains, such as PI3-K, Grb-2, Nck, and Crk. While tyrosine phosphorylation of insulin receptor substrates propagates signaling from insulin, serine and threonine phosphorylation by MAPK's, PKA, and PKC may reduce insulin signaling.
<b>Species Reactivity:</b>	Human and mouse. Other species were not tested.
<b>Applications:</b>	This antibody is suitable for use in Western blotting.
<b>Suggested Working Dilutions:</b>	For Western blotting, the recommended concentration is 1 µg/mL. The optimal antibody concentration should be determined for each specific application.
<b>Recommended Positive Control:</b>	Human Jurkat and HeLa cells, and mouse NIH3T3 cells.
<b>Storage:</b>	Store at 2-8°C. For long term storage, aliquot into small volumes and store at -20°C. Avoid repeated freeze-thaw cycles to prevent denaturing the antibody.

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

[www.invitrogen.com](http://www.invitrogen.com)

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

PIAHO1532 (Rev 10/08)

(Page 1 of 2) 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](http://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.

**References:**

Dearth, R.K. et al, (2006) Mammary tumorigenesis and metastasis caused by overexpression of insulin receptor substrate 1 (IRS-1) or IRS-2. *Mol Cell Biol.* 26:9302-9314.

Greene, M.W., et al. (2003) Modulation of insulin stimulated degradation of human insulin receptor substrate-1 by serine 312 phosphorylation. *J. Biol. Chem.* 278:8199-8211.

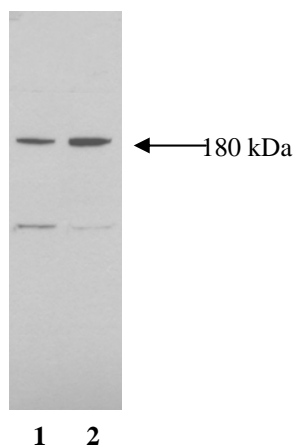
Sun, X.J., et al. (1995) Role of IRS-2 in insulin and cytokine signalling. *Nature* 377:173-177.

Mothe I., and E. van Obberghen (1996) Phosphorylation of insulin receptor substrate-1 on multiple serine residues, 612, 632, 662, and 731, modulates insulin action. *J. Biol. Chem.* 271:11222-11227.

White, M.F. (1998) The IRS-signalling system: a network of docking proteins that mediate insulin action. *Mol. Cell. Biochem.* 182(1-2):3-11.

**Related Products:**

Anti IRS-1 monoclonal antibody	Cat. # AHO1322
IRS-1 (total) ELISA	Cat. # KHO0511
IRS-1 pS312 ELISA	Cat. # KHO0521
Anti-IRS-2 [pS721] antibody	Cat. # 44-828

**Western Blot Analysis**

Proteins from cell extracts of human Jurkat and HeLa cells were resolved by SDS-PAGE and transferred to PVDF. The membranes were incubated with this IRS-2 monoclonal antibody (clone 8S15) at a concentration of 1 µg/mL for two hours at room temperature (Lane 1 – Jurkat, Lane 2 – HeLa). After washing, the membranes were incubated with a goat F(ab')<sub>2</sub> anti-mouse IgG HRP conjugated antibody (Cat. #65-6420) at a 1:5000 dilution. Bands were detected with ECL Chemiluminescent Substrate and Pierce CL-X Posure™ film.

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

[www.invitrogen.com](http://www.invitrogen.com)

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