



Catalog Number	Product Name	Quantity	Clonality, clone (isotype)	Reactive species	Applications	Reg. Status
434300	Mouse anti-human SHP	100 µg	mAb clone N7519 (Ms IgG1)	Hu	WB, ELISA	RUO

## Mouse Anti-Human Small Heterodimer Partner

**Description** Small heterodimer partner SHP (NR0B2) is an orphan receptor that is expressed in liver and at lower level in heart, adrenal gland, spleen and pancreas. Since SHP binds to several receptors, including retinoid receptors RAR, RXR and thyroid hormone receptor, the function is suggested as a negative regulator of receptor-dependent signaling pathways.

**Nomenclature** NR0B2

**Genbank** L76571

**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 *E. coli*-expressed recombinant human SHP (1-257 aa).

**Specificity** This antibody specifically recognizes human SHP. 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 µg/mL
Non reducing Western Blot	3 µg/mL
ELISA	0.5 µg/mL
Immunoprecipitation	Not tested
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.

[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)

PI434300

(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](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.

## 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](http://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](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)

PI434300

(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](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.