



## Mouse (monoclonal) Anti-Human SLP-76

### PRODUCT ANALYSIS SHEET

---

<b>Catalog Number:</b>	AHO0692
<b>Lot Number:</b>	See product label
<b>Quantity/Volume:</b>	100 µg/200 µL
<b>Clone Number:</b>	AS55
<b>Isotype:</b>	IgG2a
<b>Form of the 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 tissue culture supernatant by Protein G affinity chromatography. Sterile filtered (0.22 µm).
<b>Immunogen:</b>	SLP-76 fusion protein
<b>Myeloma/Fusion Partners:</b>	Hybridoma derived from BALB/c splenocytes.
<b>Specificity:</b>	SH2-domain Leukocyte Protein (SLP-76) is a 76 kDa cytoplasmic protein involved in T-cell receptor mediated signaling. SLP-76 is a substrate for ZAP70 and Syk and is phosphorylated upon T-cell receptor activation. SLP-76 is also known to associate with Grb2. SLP-76 is highly expressed in T-cell and monocytic cell lines and, to a lesser degree in B-cell lines. SLP-76 expression is crucial for receptor-mediated signaling in platelets and in T lymphocytes.
<b>Species Reactivity:</b>	Human. Other species were not tested.
<b>Applications:</b>	This monoclonal antibody is suitable for use in Western blotting and immunoprecipitation.
<b>Suggested Working Dilutions:</b>	For Western blotting, we recommend using the antibody at 1 µg/mL to detect SLP-76 from extract of $1 \times 10^5$ Jurkat cells per lane. For immunoprecipitation, we recommend using 1-5 µg antibody per $1 \times 10^6$ Jurkat cells. The optimal concentration should be determined for each specific application.
<b>Storage:</b>	Store at -20°C. Upon initial thawing, apportion into working aliquots and store at -20°C. Avoid repeated freeze-thaw cycles to prevent denaturing the antibody.
<b>Expiration Date:</b>	Expires one year from date of receipt when stored as instructed.

**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)

PI AHO0692

(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.

**References:**

Jackman, K.J., D.G. Motto, Q. Sun, M. Tanemoto, C.W. Turk, G.A. Peltz, G.A. Koretzky, and P.R. Findell (1995) Molecular cloning of SLP-76, a 76-kDa tyrosine phosphoprotein associated with Grb2 in T cells. *J. Biol. Chem.* 270:7029-7032.

Wardenburg, J.B., C. Fu, J.K. Jackman, H. Flotow, S.E. Wilkinson, D.H. Williams, R. Johnson, G. Kong, A.C. Chan, and P.R. Findel (1996) Phosphorylation of SLP-76 by the ZAP-70 protein-tyrosine kinase is required for T-cell receptor function. *J. Biol. Chem.* 271:19641-19644.

**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)

PI AHO0692

(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.