



Qty: 3 x 20 µg

Phosphoprotein Antibody

Sampler Pack

Catalog No. 90-0200

Lot No.

Phosphoprotein Antibody Sampler Pack

FORM

All antibodies are supplied separately in phosphate buffered saline, pH 7.4, containing 0.1% sodium azide. All antibodies in this sampler pack are specifically purified by antigen-affinity-purification.

ANTIBODY DESCRIPTIONS

Catalog Number	61-8100	71-8200	13-6600
Antibody	Rabbit anti-Phosphoserine	Rabbit anti-Phosphothreonine	Mouse anti-Phosphotyrosine
Host	Rabbit (polyclonal)	Rabbit (polyclonal)	Mouse (monoclonal cocktail with clones PY-7E1, PY20)
Quantity	20 µg/80 µl	20 µg/80 µl	20 µg/40 µl
Cell lysates tested	NIH 3T3 (+/- TPA) cells, K562 cells, and EGF-stimulated A431 cells.	NIH 3T3 (+/- TPA) cells, K562 cells, and EGF-stimulated A431 cells.	NIH 3T3 cells (+/- TPA), K562 cells, EGF-stimulated A431 cells, MCF-7 cells, mouse brain.
Application	Suggested Starting Amounts*		
ELISA	0.1-1.0 µg/ml	0.1-1.0 µg/ml	0.1-0.2 µg/ml
Immunoprecipitation	ND	5-10 µg	2-5 µg
Tissue Staining	ND	ND	ND
Western blotting ⁽⁸⁻¹⁹⁾	1.0-2.0 µg/ml	1.0-2.0 µg/ml	0.5-1.0 µg/ml

* Optimal concentrations should be determined by the researcher for each specific application. ND: Not determined.

NOTE

Milk-derived blocking solutions reportedly contain phosphoproteins that may inhibit phosphoamino acid antibody binding and therefore should be avoided. Invitrogen's Membrane Blocking Solution (Cat. No. 00-0105) is optimized for use with anti-phosphoamino acid antibodies and provides enhanced blocking of non-specific signal. A 3% BSA (bovine serum albumin) solution may also be used.

STORAGE

Store at 2-8°C for up to one month. Store at -20°C for long term storage. Avoid repeated freezing and thawing.

(cont'd)

www.invitrogen.com

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

PI900200

(Rev 12/08) 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.

CITATIONS

Polyclonal Rabbit anti-pSer (Cat. No. 61-8100)Application

Glogauer, M., et al., <i>J. Biol. Chem.</i> 273: 1689-1698 (1998)	WB
Yagi, K., et al., <i>J. Biol. Chem.</i> 274: 703-709 (1997)	WB
Jin, D-Y, et al., <i>Cell</i> 93: 81-91 (1998)	WB
Fleming, I., et al., <i>J. Biol. Chem.</i> 272: 33105-33110 (1997)	WB
Nishimura, R., et al., <i>J. Biol. Chem.</i> 273: 1872-1879 (1998)	WB
Kamiguti, A., et al., <i>J. Biol. Chem.</i> 272: 32599-32605 (1997)	WB
Kawabata, M., et al., <i>EMBO J.</i> 17: 4056-4065 (1998)	WB
Mahoney, C.W., et al., <i>Anal. Biochem.</i> 268:371-6 (1999)	ELISA
Reilein, A.R., et al., <i>J. Cell Biol.</i> , 142: 803-813 (1998)	WB
Tuvia, S., et al., <i>J. Cell Biol.</i> , 141:1551-1561 (1998)	WB
Musarò, S., et al., <i>Nature</i> , 400:581-585 (1999)	WB

Polyclonal Rabbit anti-pThr (Cat. No. 71-8200)

Wilson, J.R., and T. Biden, <i>J. Biol. Chem.</i> 273: 22729-22737 (1998)	WB
Kamiguti, A., et al., <i>J. Biol. Chem.</i> 272: 32599-32605 (1997)	WB
Fleming, I., et al., <i>J. Biol. Chem.</i> 272: 33105-33110 (1997)	WB, IP
Fleming, I., et al., <i>J. Biol. Chem.</i> 274: 12753-12758 (1997)	WB
Rintamäki, E., et al., <i>J. Biol. Chem.</i> 272: 30476-30482 (1997)	WB
Mahoney, C.W., et al., <i>Anal. Biochem.</i> 268:371-6 (1999)	ELISA

PY-Plus™ Mouse Monoclonal Cocktail (clones PY-7E1, PY-1B2, and PY20, cat. no. 13-6600)

Bourguignon, L.Y., et al., <i>J. Biol. Chem.</i> 272: 27913-27918 (1997)	IP
Conus, N.M., et al., <i>J. Biol. Chem.</i> 273: 4776-4782 (1998)	WB
Hua, C.T., et al., <i>J. Biol. Chem.</i> 273: 28332-28340 (1998)	WB

RELATED PRODUCTS

Product	PAD*/clone	Cat. No.	
Rb x Phosphoserine	Z-PS1	61-8100	
Phosphoserine Ab Inhibitor	----	79-0001	
Rb x PS/PT/PY (pan)	polyclonal	61-8300	
Phosphotyrosine Sampler Pack	6 antibodies	90-0100	
Rb x Phosphothreonine	Z-PT1	71-8200	
Ms x Phosphothreonine	PT-5H5	13-9200	
Phosphothreonine Ab Inhibitor		79-0002	
Rb x Phosphotyrosine	Z-PY1	61-5800	
Rb x Phosphotyrosine-HRP	Z-PY1	61-5820	
Rb x Phosphotyrosine-Sepharose®	Z-PY1	61-5841	
Ms x Phosphotyrosine	PY-7E1	13-5900	
Ms x Phosphotyrosine-HRP	PY-7E1	13-5920	
Ms x Phosphotyrosine	PY-1B2	13-6300	
Ms x Phosphotyrosine	PY20	03-7700	
Ms x Phosphotyrosine (1 mg size)	PY20	03-7799	Excellent Value!
Ms x Phosphotyrosine-HRP	PY20	03-7720	
Ms x Phosphotyrosine-AP	PY20	03-7722	
Ms x Phosphotyrosine-Biotin	PY20	03-7740	
Ms x Phosphotyrosine-Sepharose®	PY20	03-7742	
Ms x Phosphotyrosine	Z027	03-5800	
Phosphotyrosine Ab inhibitor	---	79-0003	
PY-Plus™ Cocktail	3 Clones	13-6600	
PY-Plus™ Cocktail-HRP	3 Clones	13-6620	

PAD: Polyclonal Antibody Designation

Protein A	Sepharose® 4B	10-1041
rec-Protein G	Sepharose® 4B	10-1241

Zymed® and ZyMAX™ are trademarks of Zymed Laboratories Inc. Cy™3 and Cy™5 are trademarks of Amersham Life Sciences, Inc. Sepharose® is a registered trademark of Pharmacia LKB.

For Research Use Only

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

PI900200

(Rev 12/08) 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.