

## MICROARRAYS OF CELLS EXPRESSING DEFINED cDNAs

Ziauddin, J. and Sabatini, D.M. (2001) *Nature* **411**, 107.

This article describes a microarray-driven gene expression system. Mammalian cells are cultured on a glass slide onto which a series of cDNAs have been printed. Cells that grow on the printed area are transfected with the DNA, creating localized patches of cells transfected with a particular cDNA. The slides can be screened to identify drug targets or to discover gene products involved in particular cell functions or pathways.

Transfected cell microarrays were used to screen a collection of 192 epitope-tagged cDNAs cloned in expression vectors and enriched for cell signalling molecules. HEK293T cells were transfected and screened with phosphospecific antibodies in order to detect increased phosphotyrosine or activated, phosphorylated forms of the MAPK family. The authors used Promega's Anti-ACTIVE® p38 pAb and Anti-ACTIVE® JNK pAb in the screening. Five of six phosphotyrosine positive cell clusters expressed known tyrosine kinases. Two of the clusters were positive for phosphorylated p38 and JNK (Table 1).

### Ordering Information

Product	Size	Cat.#
Anti-ACTIVE® p38 pAb, Rabbit, (pTEpY)	100µl	V1211
Anti-ACTIVE® JNK pAb, Rabbit, (pTPpY)	40µl	V7931
	120µl	V7932

**Table 1. Phenotypes Induced for cDNAs Expressed in Transfected Cell Microarray.**

Grid: Coordinate	Phenotype	Gene Name
1:A2	P-Tyr	TrkC
1:C7	P-Tyr, p-JNK, p-p38	Syk
1:C9	P-Tyr	unnamed
1:C11	P-Tyr	Syn
1:E3	P-Tyr	Lck
1:F6	P-Tyr, p-JNK, p-p38	Blk
1:C2	cytoskeleton changes	β-PDGF receptor
2:E8	apoptosis	TRAIL receptor 2
2:D10	cell adhesion	NK receptor
2:F7	cell adhesion	CD36

P-Tyr, p-JNK and p-p38 indicates that the cluster had staining of phosphotyrosine, phospho-JNK, and phospho-p38, respectively, above background. Reprinted with permission of the author and *Nature* from Ziauddin, J. and Sabatini, D.M. (2001) *Nature* **411**, 107-110.

### Related Products

Product	Size	Cat.#
Anti-ACTIVE® MAPK pAb, Rabbit, (pTEpY)	40µl	V8031
Anti-Phosphotyrosine, pAb	200µg	V2171

Anti-ACTIVE is a trademark of Promega Corporation and is registered with the U.S. Patent and Trademark Office.

#### Promega Corporation

2800 Woods Hollow Road  
Madison, WI 53711-5399 USA  
Tel: 608-274-4330  
Fax: 608-277-2516  
Toll-Free: 800-356-9526  
Toll-Free Fax: 800-356-1970  
Internet: www.promega.com

#### Promega Biosciences, Inc.

A division of Promega Corporation  
San Luis Obispo, CA

#### Australia, Sydney

Tel: 02-9565-1100  
Fax: 02-9550-4454  
Free Call: 1-800-225-123  
Free Fax: 1-800-626-017  
E-mail: aus\_custserv@au.promega.com

#### China, Beijing

Tel: 10 6849 8287  
Fax: 10 6849 8390  
E-mail: promega@promega.com.cn

#### France, Lyon

Tél: 04-37-22-50-00  
Fax: 04-37-22-50-10  
Numéro Vert: 0-800-48-79-99  
E-mail: fr\_custserv@fr.promega.com

#### Germany/Austria, Mannheim

Tel: (+49) (0) 621-8501-0  
Fax: (+49) (0) 621-8501-222  
Free Tel: 00800-77663422  
Free Fax: 00800-77663423  
E-mail: de\_custserv@de.promega.com

#### Italy, Milan

Tel: 02 290 6651  
Fax: 02 2901 7365  
Numero Verde: 800 69 1818  
E-mail: it\_custserv@it.promega.com

#### Japan, Tokyo

Tel: 03-3669-7981  
Fax: 03-3669-7982  
E-mail: prometec@jp.promega.com

#### The Netherlands, Leiden, NL

Tel: (+31) (0) 71-5324244  
Fax: (+31) (0) 71-5324907  
Free Tel NL: 0800-0221910  
Free Fax NL: 0800-0226545  
E-mail: bnl\_custserv@nl.promega.com

#### Belgium and Luxembourg, Leiden, NL

Tel: (+31) (0) 71-5324244  
Fax: (+31) (0) 71-5324907  
Free Tel BE: 0800-18098  
Free Fax BE: 0800-16971  
E-mail: bnl\_custserv@nl.promega.com

#### Pacific Asia Region

Tel: 65 254 5265  
Fax: 65 254 8645  
E-mail: nng@promega.com

#### Switzerland, Wallisellen

Tel: 01 878 90 00  
Fax: 01 878 90 10  
E-mail: catalys\_custserv@catalys.promega.com

#### United Kingdom, Southampton

Tel: 023 8076 0225  
Fax: 023 8076 7014  
Free Phone: 0800 378994  
Free Fax: 0800 181037  
E-mail: ukcustserve@uk.promega.com



PRINTED IN THE U.S.A.