

Multiplexable Assay Matrix

Total Target Assay	Tested Antibody Species*	Akt	ATF-2	Bcr-Abl	c-Jun	CREB	ERK2	ERK1/2	HSP27	IκB-α	JNK	MEK1	p38 MAPK	p53	p90RSK
Akt	H, M, R														
ATF-2	H														
Bcr-Abl	H, M, R														
c-Jun	H, M, R														
CREB	H, M, R														
ERK2	H, M, R														
ERK1/2	H, M														
HSP27	H, K														
IκB-α	H														
JNK	H														
MEK1	H, M, R														
p38 MAPK	H, M														
p53	H														
p90RSK	H, M, R														

* H, human; M, mouse; R, rat; K, monkey.

Multiplexing Key**



Multiplexable Not multiplexable

** Representative data are based on lysates provided with the assay (in-house cell lines).