



PRIMARY ANTIBODIES

Isoform Specific Antibodies

Product Name	Catalog #
Anti-AKT3	A18-63R
Anti-AMPK alpha 1	P47-63R
Anti-CAMK1	C07-60M
Anti-CHK2	C48-60M
Anti-ERK1	M29-60BM
Anti-ERK1	M29-60M
Anti-GSK3 alpha	G08-63R
Anti-GSK3 beta	G09-60M
Anti-HSP90 beta	H36-63R
Anti-JAK2	J02-63R

Product Name	Catalog #
Anti-JNK1	M33-63R
Anti-JNK2	M34-60M
Anti-MEK1	M02-60M
Anti-MEK2	M03-63R
Anti-p38 alpha	M39-63R
Anti-p38 beta	M36-63R
Anti-p38 gamma	M37-63R
Anti-PAK1	P02-63R
Anti-PKC delta	P64-63CR
Anti-PKC delta	P64-63R

Product Name	Catalog #
Anti-PKC mu	P72-63R
Anti-PKC theta	P74-63R
Anti-SMAD2	S11-63R
Anti-STAT1 beta	S52-63R
Anti-STAT3	S54-63BR
Anti-STAT3	S54-63R
Anti-STAT4	S55-63R
Anti-STAT5	S56-63BR
Anti-STAT5	S56-63R

Primary Antibodies

Product Name	Catalog #
Anti-4EBP1	E35-563R
Anti-ADD1	A38-63R
Anti-AFX	A42-63R
Anti-AKT (pan)	A16-63BR
Anti-AKT (pan)	A16-63R
Anti-AKT1	A16-60M
Anti-AKT1	A16-63CR
Anti-AKT2/3	A17-63BR
Anti-ALK2	A06-63R
Anti-ALK4	A07-63R
Anti-ALK7	A08-63R
Anti-alpha 1-Actin	A03-360M
Anti-AMPK alpha 1/2	P47-63BR
Anti-AOS1	A80-63R
Anti-APBA1	A86-363R
Anti-APC1	A87-363R
Anti-APC2	A88-363R
Anti-APC3	A89-363R
Anti-APC4	A90-363R
Anti-APC5	A91-363R
Anti-APC6	A92-363R
Anti-APC7	A93-363R
Anti-APC8	A94-363R
Anti-APC10	A95-363R
Anti-APC11	A96-363R
Anti-APC13	A97-363R
Anti-ASK1	M13-60BM
Anti-ASK1	M13-60M
Anti-ASK1	M13-63BR
Anti-ASK1	M13-63CR
Anti-ASK1	M13-63R

Product Name	Catalog #
Anti-ATF2	A10-63R
Anti-BAD	B03-363R
Anti-BAG1	B05-363BR
Anti-BAG1	B05-363R
Anti-BAK1	B09-363R
Anti-BAX	B14-363R
Anti-BCL2	B25-63BR
Anti-BCL2	B25-63CR
Anti-BCL2	B25-63R
Anti-BCL2L1	B26-363R
Anti-BCL2L10	B26-363JR
Anti-BCL2L14	B26-363NR
Anti-beta-Actin	A03-60BM
Anti-beta-Actin	A03-63BR
Anti-BID	B59-363BR
Anti-BID	B59-363R
Anti-BIRC2	B279-63R
Anti-BOK	B70-363R
Anti-BRAF	B08-63R
Anti-BRSK1	B13-63BR
Anti-BRSK1	B13-63R
Anti-BRSK2	B14-63R
Anti-BTK	B10-63R
Anti-BUB1	B11-63BR
Anti-BUB1	B11-63R
Anti-c-JUN	C05-63R
Anti-CASK	C19-63R
Anti-CASKIN1	C13-363R
Anti-CD5L (CT-2)	C61-363BR
Anti-CD5L (CT-2)	C61-363R
Anti-CDC25A	C04-63R

Product Name	Catalog #
Anti-CDC42	C08-363R
Anti-CDK5	C33-60M
Anti-CHK2	C48-63R
Anti-CREB	C49-63R
Anti-CyclinO	C85-363BR
Anti-CyclinO	C85-363R
Anti-CyclinT1	C86-363BR
Anti-DAPK2	D02-63R
Anti-DAPK3	D03-63R
Anti-DCAMKL1	D14-63R
Anti-DCAMKL2	D15-63R
Anti-DCAMKL3	D16-63R
Anti-DLK1	M20-63R
Anti-DOK1	D42-63R
Anti-DRAK2 (STK17B)	S34-63R
Anti-DYRK1A	D09-63R
Anti-DYRK2	D10-63R
Anti-EGFR	E10-63BR
Anti-EGFR	E10-63R
Anti-EIF2AK2 (PKR)	P80-63R
Anti-EIF2AK2 (PKR)	P80-63BR
Anti-EIF2AK2 (PKR)	P80-63CR
Anti-eIF4E	E34-63R
Anti-ELK1	E64-63R
Anti-ERK1	M29-60CM
Anti-ERK1/2	M29-60EM
Anti-ERK2	M28-63R
Anti-EZH1	E95-363R
Anti-EZH2	E96-363R
Anti-FGF4	F824-63R
Anti-FLT3	F12-63R

Antibodies

Primary Antibodies

Product Name	Catalog #
Anti-FOXO1	F19-63R
Anti-GAB1	G03-63R
Anti-GAPDH	G13-61M
Anti-GAPDH	G13-63R
Anti-GATA1	G19-63R
Anti-GSK3 alpha/ beta	G09-63R
Anti-HDAC1	H83-63R
Anti-HDAC2	H84-63R
Anti-HDAC4	H86-63R
Anti-HDAC5	H87-63R
Anti-HDAC6	H88-63R
Anti-HDAC7	H89-63R
Anti-HDAC8	H90-63R
Anti-HDAC9	H91-63R
Anti-HDAC11	H93-63R
Anti-HER2	E27-63R
Anti-HSF1	H25-63R
Anti-HSP27	H30-63R
Anti-IKK alpha	C51-63BR
Anti-IKK alpha	C51-63CR
Anti-IKK alpha	C51-63R
Anti-IKK beta	I03-63R
Anti-IKK epsilon	I04-63R
Anti-IKK gamma	I25-63R
Anti-IRAK1	I09-63R
Anti-IRAK2	I10-63BR
Anti-IRAK2	I10-63CR
Anti-IRAK2	I10-63R
Anti-IRAK3	I11-63R
Anti-IRAK4	I12-63R
Anti-IRS1	I40-363R
Anti-JMJD1C	J48-363R
Anti-JNK2/3	M34-63R
Anti-KDM1A	K421-63BR
Anti-KDM1A	K421-63R
Anti-KDM3A	K423-63R
Anti-KDM3B	K423-63BR
Anti-KDM4A	K424-63R
Anti-KDM4B	K424-63BR
Anti-KDM6B	K426-63BR
Anti-KDM8	K428-63R
Anti-LKB1 (STK11)	L15-63R
Anti-MAPKAPK2/3	M40-63R
Anti-MARCKS	M02-562R
Anti-MBP	M42-563C
Anti-MEK1/2	M02-63R
Anti-mTOR	F17-63BR
Anti-mTOR	F17-63R
Anti-MYT1L	M92-363R
Anti-NF-kB p105	N12-63R
Anti-NIK	N22-63R
Anti-p16INK4A	C30-363R
Anti-p21CIP1	C27-63R

Product Name	Catalog #
Anti-p38 alpha	M39-60M
Anti-p53	P05-63R
Anti-p57KIP2	C29-363R
Anti-p70S6K	R21-63BR
Anti-p70S6K	R21-63R
Anti-PAK2	P03-63BR
Anti-PAK2	P03-63R
Anti-PAK4	P05-163R
Anti-PAK6	P06-63R
Anti-PAK7	P07-63R
Anti-PDGFB	P842-63R
Anti-PHLP1	P71-263R
Anti-PHLP2	P72-263R
Anti-PIAS1	P280-63R
Anti-PIAS2	P281-63R
Anti-PIAS3	P282-63R
Anti-PIAS4	P283-63R
Anti-PKAc	P51-63R
Anti-PKC delta/ theta	P64-63DR
Anti-PKC zeta	P75-63R
Anti-PLAC1	P38-363R
Anti-PLAC2	P38-363BR
Anti-PLAC3	P38-363CR
Anti-PLAC4	P38-363DR
Anti-PLC gamma 1	P40-63R
Anti-PTEN	P23-263BR
Anti-PTEN	P23-263R
Anti-PTK7	P95-63BR
Anti-PTK7	P95-63R
Anti-PTPN11 (SHP2)	P38-63BR
Anti-PTPN11 (SHP2)	P38-63CR
Anti-PYK2	P92-63R
Anti-QIK	S15-63R
Anti-RAF1	R01-63R
Anti-Raptor	R47-363BR
Anti-Raptor	R47-363R
Anti-Rel A	R29-63BR
Anti-Rel A	R29-63R
Anti-RHEB	R42-363BR
Anti-RHEB	R42-363R
Anti-RIPK1	R07-63R
Anti-RIPK2	R08-63BR
Anti-RIPK2	R08-63R
Anti-RIPK3	R09-63R
Anti-ROCK1	R10-63R
Anti-ROCK2	R11-63R
Anti-RSK (pan)	R15-63BR
Anti-RSK1	R15-63CR
Anti-RSK1	R15-63R
Anti-RSK3	R16-63R
Anti-SGK1	S06-63R
Anti-SH2D1A (SAP)	S03-63R
Anti-SHP2	P38-63R

Product Name	Catalog #
Anti-SIK1	S14-63R
Anti-SIK3	S12-63R
Anti-SIRT1	S35-463R
Anti-SIRT2	S36-463BR
Anti-SIRT2	S36-463R
Anti-SIRT3	S37-463R
Anti-SIRT4	S38-463R
Anti-SIRT5	S39-463R
Anti-SIRT6	S40-463R
Anti-SIRT7	S41-463R
Anti-SMC1	S19-63R
Anti-SNCA	S22-360M
Anti-STAT1 alpha	S52-63BR
Anti-STAT1 beta	S52-63CR
Anti-STAT3	S54-63CR
Anti-Stathmin	S71-363BR
Anti-Stathmin	S71-363R
Anti-STK39 (STLK3)	S51-63BR
Anti-STK39 (STLK3)	S51-63R
Anti-SUMO1	S293-63R
Anti-SUMO2/3	S294-63BR
Anti-SUMO3	S295-63R
Anti-TAB1	T01-363R
Anti-TAK1	M15-63R
Anti-TBK1	T02-63R
Anti-TET3	T638-63R
Anti-TRAF2	T57-363BR
Anti-TRAF2	T57-363R
Anti-TRAF3	T58-363BR
Anti-TRAF3	T58-363R
Anti-Trail	T880-63R
Anti-TSC1	T71-363BR
Anti-TSC1	T71-363R
Anti-TSC2	T72-363BR
Anti-TSC2	T72-363R
Anti-TTBK1	T18-63R
Anti-TTBK2	T18-63BR
Anti-Tubulin alpha	T91-63R
Anti-Tubulin beta	T92-63R
Anti-TYK2 (JTK1)	T21-63R
Anti-UBA2 (SAE2)	U202-63R
Anti-UBC9	U19-63R
Anti-UBE2N (UBC13)	U231-63R
Anti-UBE2V1 (UEV1A)	U238-63R
Anti-UBLE1A (SAE1)	U208-63R
Anti-VASP	V12-63R
Anti-VAV	V14-63R
Anti-Vimentin	V49-362C
Anti-VLK	V05-163R
Anti-XIAP	X08-63R
Anti-ZIP11	Z45-363R

For additional details and antibody applications, please visit signalchem.com

20140115