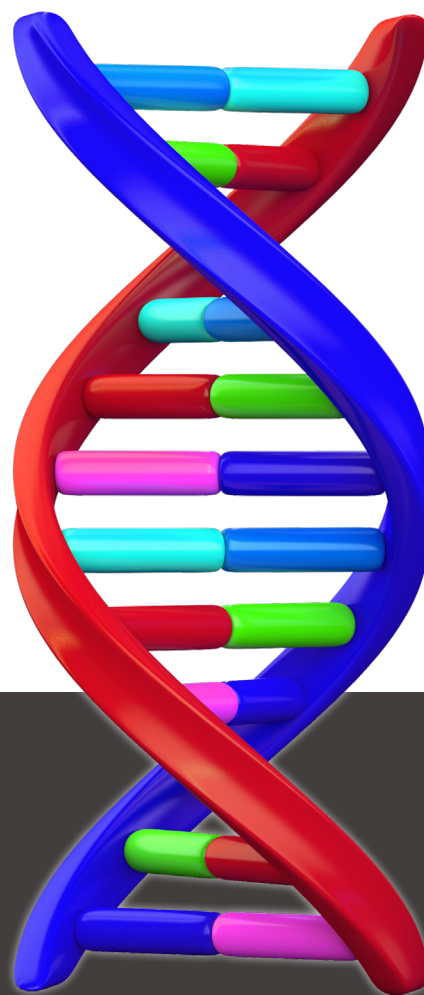


siRNA Catalog



siRNA

In the absence of mutants or as a complement to traditional genetic studies, RNA interference (RNAi) is an excellent tool for studying protein function in cells. RNAi is effective for silencing genes in a highly specific manner and many organisms such as plants and nematodes are amenable to gene silencing via the introduction of long double-stranded RNA (dsRNA) fragments. However, in mammalian systems, the effectiveness of long dsRNAs is limited because secondary effects are prevalent such as the induction of interferons and cytokines. Ultimately, these side-effects lead to unwanted changes to cellular responses. To avoid many of the issues of resulting from the application of long dsRNAs, small interfering RNAs (siRNAs) are typically used in mammalian cells and since siRNAs are downstream mediators of RNAi, the same desired targeted gene silencing is achieved.

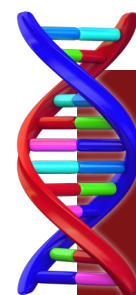
Gene knockdown studies are an excellent complement to most research programs. As a genetic research tool, siRNAs enable scientists to effectively study cell signaling pathways in vivo. For instance, by using siRNA to deplete signaling pathway components, scientists are able to understand the effects of decreased signaling. This information is relevant to understanding the overall effects of pharmaceutical agents that are intended modulate the activity of the aforementioned pathway.

As a provider of complete solutions for researchers, SignalChem also offers the appropriate positive and negative control siRNAs. Our alpha-Actin and GAPDH positive controls are among the prevailing industry standard targets. SignalChem's negative controls consist of scrambled nucleotide sequences, which have no known homology to coding sequences in humans. Among our negative controls, researchers can choose from unlabeled, FAM-labeled and Cy3 labeled products. FAM and Cy3 are fluorescent markers that enable qualitative assessment of cellular uptake, distribution and intracellular localization of siRNAs. Thus, these FAM and Cy3 Labeled Negative Control siRNAs are commonly used to optimize transfection conditions, which are monitored by fluorescence microscopy or fluorescence-activated cell sorting (FACS).

siRNA Sets

Product Name	Cat. No.	Product Name	Cat. No.	Product Name	Cat. No.
14-3-3 alpha/beta siRNA Set I	Y71-911	AMPK A1 siRNA Set I	P47-911	ATF1 siRNA Set I	A09-911B
14-3-3 epsilon siRNA Set I	Y75-911	AMPK A2 siRNA Set I	P48-911	ATF2 siRNA Set I	A10-911
14-3-3 sigma siRNA Set I	Y86-911	AMPK B1 siRNA Set I	P49-911	ATG13 siRNA Set I	A113-911
14-3-3 theta siRNA Set I	Y84-911	AMPK B2 siRNA Set I	P50-911	AURORA A siRNA Set I	A28-911
14-3-3 zeta siRNA Set I	Y92-911	AMPK G1 siRNA Set I	P51-911B	AURORA B siRNA Set I	A31-911
4EBP1 siRNA Set I	E35-911	AMPK G2 siRNA Set I	P52-911C	AURORA C siRNA Set I	A33-911
AATK siRNA Set I	A02-911	AMPK G3 siRNA Set I	P53-911C	AXL siRNA Set I	A34-911
ABL1 siRNA Set I	A03-911	Amyloid beta 42 siRNA Set I	A06-911B	BAD siRNA Set I	B03-911
ABL2 siRNA Set I	A04-911	Annexin 5 siRNA Set I	A61-911	BAG1 siRNA Set I	B05-911B
ACK siRNA Set I	A05-911	APBA1 siRNA Set I	A86-911	BAX siRNA Set I	B14-911
ACVR2A siRNA Set I	A35-911	APC1 siRNA Set I	A87-911	BCKDK siRNA Set I	B15-911
ACVR2B siRNA Set I	A36-911	APC2 siRNA Set I	A88-911	BCL2 siRNA Set I	B25-911
ADD1 siRNA Set I	A38-911	APC3 siRNA Set I	A89-911	BCL2L1 siRNA Set I	B26-911
AFX siRNA Set I	A42-911	APC4 siRNA Set I	A90-911	BCR siRNA Set I	B01-911
AKT1/PKB alpha siRNA Set I	A16-911	APC5 siRNA Set I	A91-911	beta-Actin siRNA Set I	A03-911C
AKT2/PKB beta siRNA Set I	A17-911	APC6 siRNA Set I	A92-911	BID siRNA Set I	B59-911
AKT3 siRNA Set I	A18-911	APC7 siRNA Set I	A93-911	BIRC3 siRNA Set I	B280-911
ALK1 siRNA Set I	A09-911	APC8 siRNA Set I	A94-911	BLK siRNA Set I	B02-911
ALK2 siRNA Set I	A06-911	APC10 siRNA Set I	A95-911	BMP-2 siRNA Set I	B851-911
ALK3 (BMPR1A) siRNA Set I	B04-911	APC11 siRNA Set I	A96-911	BMP2K siRNA Set I	B03-911B
ALK6 (BMPR1B) siRNA Set I	B05-911	APC13 siRNA Set I	A97-911	BMP-4 siRNA Set I	B853-911
ALK7 siRNA Set I	A08-911	ARAF siRNA Set I	A24-911	BMPR2 siRNA Set I	B06-911
alpha 1-Actin siRNA Set I	A03-911B	ASK1 siRNA Set I	M13-911	BMX siRNA Set I	B07-911

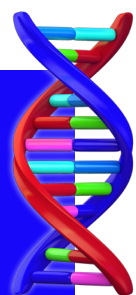
Product Name	Cat. No.	Product Name	Cat. No.	Product Name	Cat. No.
BOK siRNA Set I	B70-911	CK1 gamma 2 siRNA Set I	C68-911B	EPHA3 siRNA Set I	E15-911
BRAF siRNA Set I	B08-911	CK1 gamma 3 siRNA Set I	C68-911C	EPHA4 siRNA Set I	E16-911
BRC A1 siRNA Set I	B87-911	CK2 alpha 1 siRNA Set I	C70-911	EPHA6 siRNA Set I	E18-911
BRK siRNA Set I	P94-911	CK2 alpha 2 siRNA Set I	C71-911	EPHA7 siRNA Set I	E19-911
BRSK1 siRNA Set I	B13-911	c-KIT siRNA Set I	K06-911	EPHA8 siRNA Set I	E20-911
BRSK2 siRNA Set I	B14-911B	CLK1 siRNA Set I	C57-911	EPHB1 siRNA Set I	E21-911
BTK siRNA Set I	B10-911	CLK2 siRNA Set I	C58-911	EPHB2 siRNA Set I	E22-911B
BUB1 siRNA Set I	B11-911	CLK3 siRNA Set I	C60-911	EPHB3 siRNA Set I	E23-911B
BUB1B siRNA Set I	B12-911	c-MER siRNA Set I	M51-911	EPHB4 siRNA Set I	E24-911
Calmodulin1 siRNA Set I	C01-911	Cofilin 1 siRNA Set I	C37-911	ERK1 siRNA Set I	M29-911
CAMK1 siRNA Set I	C07-911	COT siRNA Set I	M16-911	ERK2 siRNA Set I	M28-911
CAMK1 beta siRNA Set I	C08-911	CREB1 siRNA Set I	C49-911	ERK5 siRNA Set I	M32-911
CAMK1 delta siRNA Set I	C09-911	CRKL siRNA Set I	C53-911	EZH1 siRNA Set I	E95-911
CAMK1 gamma siRNA Set I	C10-911	CRYAA siRNA Set I	C256-911	EZH2 siRNA Set I	E96-911
CAMK2 alpha siRNA Set I	C11-911	CSK siRNA Set I	C63-911	FADD siRNA Set I	F02-911
CAMK2 beta siRNA Set I	C12-911	CyclinA2 siRNA Set I	C78-911	FAK siRNA Set I	P91-911
CAMK2 delta siRNA Set I	C13-911	CyclinC siRNA Set I	C80-911	FANCL siRNA Set I	F289-911
CAMK2 gamma siRNA Set I	C14-911	CyclinD1 siRNA Set I	C81-911	FASTK siRNA Set I	F01-911
CAMK4 siRNA Set I	C15-911	CyclinE1 siRNA Set I	C82-911	FER siRNA Set I	F16-911
CAMKK1 siRNA Set I	C17-911	CyclinH siRNA Set I	C83-911	FES siRNA Set I	F03-911
CAMKK2 siRNA Set I	C18-911	CyclinK siRNA Set I	C84-911	FGF4 siRNA Set I	F824-911
CaMKPase siRNA Set I	P06-911	CyclinO siRNA Set I	C85-911	FGFR1 siRNA Set I	F04-911
CASK siRNA Set I	C19-911	CyclinT1 siRNA Set I	C86-911	FGFR2 siRNA Set I	F05-911
CASKIN1 siRNA Set I	C13-911B	CyclinY siRNA Set I	C87-911	FGFR3 siRNA Set I	F06-911
Catenin alpha siRNA Set I	C05-911	DAPK1 siRNA Set I	D01-911	FGFR4 siRNA Set I	F07-911
Catenin beta siRNA Set I	C06-911	DAPK2 siRNA Set I	D02-911	FGR siRNA Set I	F10-911
CBP siRNA Set I	C07-911B	DAPK3 siRNA Set I	D03-911	FKBP1A siRNA Set I	F37-911
CD5L (CT-2) siRNA Set I	C61-911	DCAMKL1 siRNA Set I	D14-911	FLT1 siRNA Set I	F11-911
CDC7 siRNA Set I	C26-911	DCAMKL2 siRNA Set I	D15-911	FLT3 siRNA Set I	F12-911
CDC25A siRNA Set I	C04-911	DCAMKL3 siRNA Set I	D16-911	FLT4 siRNA Set I	F13-911
CDC25B siRNA Set I	C04-911B	DDR1 siRNA Set I	D05-911	FMS siRNA Set I	C74-911
CDC25C siRNA Set I	C04-911C	DDR2 siRNA Set I	D06-911	FOS siRNA Set I	F66-911
CDC42 siRNA Set I	C08-911B	DLK siRNA Set I	M20-911	FOXO1 siRNA Set I	F19-911
CDK1 siRNA Set I	C22-911	DNMT1 siRNA Set I	D351-911	FRK siRNA Set I	F14-911
CDK2 siRNA Set I	C29-911	DNMT3A siRNA Set I	D353-911	FRS3 siRNA Set I	F80-911
CDK3 siRNA Set I	C30-911	DNMT3B siRNA Set I	D353-911B	FYN A siRNA Set I	F15-911
CDK4 siRNA Set I	C31-911	DNMT3L siRNA Set I	D353-911C	FYN C siRNA Set I	F15-911B
CDK5 siRNA Set I	C33-911	DOK1 siRNA Set I	D42-911	GAB1 siRNA Set I	G03-911
CDK5P35 siRNA Set I	C34-911	DRAK1 (STK17A) siRNA Set I	S33-911	GAPDH siRNA Set I	G13-911
CDK6 siRNA Set I	C35-911	DRAK2 (STK17B) siRNA Set I	S34-911	GATA1 siRNA Set I	G19-911
CDK7 siRNA Set I	C36-911	DUSP22 siRNA Set I	D22-911	GCK siRNA Set I	M24-911
CDK8 siRNA Set I	C38-911	DYRK1A siRNA Set I	D09-911	GLK siRNA Set I	M25-911
CDK9 siRNA Set I	C40-911	DYRK2 siRNA Set I	D10-911	GRB2 siRNA Set I	G45-911
CDK11 siRNA Set I	C42-911	DYRK3 siRNA Set I	D11-911	GRB7 siRNA Set I	G45-911B
CDK11B (CDC2L1) siRNA Set I	C23-911	EEF2K siRNA Set I	E01-911	GRK1 siRNA Set I	R03-911
CDKL1 siRNA Set I	C43-911	EGFR siRNA Set I	E10-911	GRK2 siRNA Set I	A14-911
CDKL2 siRNA Set I	C44-911	EHMT2 siRNA Set I	E318-911	GRK3 siRNA Set I	A15-911
CDKL3 siRNA Set I	C45-911	EIF2AK1 (HRI) siRNA Set I	H07-911	GRK5 siRNA Set I	G03-911B
CDKL4 siRNA Set I	C69-911	EIF2AK2 siRNA Set I	P80-911		
CDKN3 siRNA Set I	C01-911B	EIF2AK3 siRNA Set I	E11-911		
CHK1 siRNA Set I	C47-911	EIF2AK4 siRNA Set I	E12-911		
CHK2 siRNA Set I	C48-911	EIF2S1 siRNA Set I	E23-911		
c-Jun siRNA Set I	J05-911	EIF4E siRNA Set I	E34-911		
CK1 alpha 1 siRNA Set I	C64-911	ELK1 siRNA Set I	E64-911		
CK1 delta siRNA Set I	C65-911	ENPP2 siRNA Set I	E22-911		
CK1 epsilon siRNA Set I	C66-911	EPHA1 siRNA Set I	E13-911		
CK1 gamma 1 siRNA Set I	C68-911	EPHA2 siRNA Set I	E14-911		



Product Name	Cat. No.	Product Name	Cat. No.	Product Name	Cat. No.
GRK6 siRNA Set I	G04-911	HSP27 siRNA Set I	H30-911	KDM3A siRNA Set I	K423-911
GRK7 siRNA Set I	G05-911	HSP40 siRNA Set I	H32-911	KDM3B siRNA Set I	K423-911B
GSK3 alpha siRNA Set I	G08-911	HSP60 siRNA Set I	H33-911	KDM4A siRNA Set I	K424-911
GSK3 beta siRNA Set I	G09-911	HSP70 siRNA Set I	H34-911	KDM4B siRNA Set I	K424-911B
Haspin (GSG2) siRNA Set I	G10-911	HSP90 alpha siRNA Set I	H36-911	KDM4C siRNA Set I	K424-911C
HCK siRNA Set I	H02-911	ICK siRNA Set I	I01-911	KDM5B siRNA Set I	K425-911
HDAC1 siRNA Set I	H83-911	IFN-gamma siRNA Set I	I837-911	KDM5C siRNA Set I	K425-911B
HDAC2 siRNA Set I	H84-911	IGF1R siRNA Set I	I02-911	KDM6B siRNA Set I	K426-911
HDAC3 siRNA Set I	H85-911	IKK beta siRNA Set I	I03-911	KDM8 siRNA Set I	K428-911
HDAC4 siRNA Set I	H86-911	IKK epsilon siRNA Set I	I04-911	KDR siRNA Set I	K01-911
HDAC5 siRNA Set I	H87-911	IKK alpha siRNA Set I	C51-911	KHS1 siRNA Set I	M27-911
HDAC6 siRNA Set I	H88-911	IKK gamma siRNA Set I	I25-911	KSR1 siRNA Set I	K07-911
HDAC7 siRNA Set I	H89-911	IL-2 siRNA Set I	I862-911	LATS1 siRNA Set I	L01-911
HDAC8 siRNA Set I	H90-911	IL-8 (72AA) siRNA Set I	I868-911	LATS2 siRNA Set I	L02-911
HDAC9 siRNA Set I	H91-911	IL-18 siRNA Set I	I878-911	LC20 siRNA Set I	M89-911
HDAC10 siRNA Set I	H92-911	INCENP siRNA Set I	I30-911	LCK siRNA Set I	L03-911
HDAC11 siRNA Set I	H93-911	InsR siRNA Set I	I08-911	Leptin siRNA Set I	L833-911
HER2 siRNA Set I	E27-911	IP-10 siRNA Set I	I838-911	LIMK1 siRNA Set I	L04-911
HER4 siRNA Set I	E29-911	IRAK1 siRNA Set I	I09-911	LKB1 (STK11) siRNA Set I	L15-911
HERC4 siRNA Set I	H265-911	IRAK2 siRNA Set I	I10-911	LMTK3 siRNA Set I	L08-911
HGK siRNA Set I	M26-911	IRAK3 siRNA Set I	I11-911	LOK siRNA Set I	S29-911
HIPK1 siRNA Set I	H03-911	IRAK4 siRNA Set I	I12-911	LOXL1 siRNA Set I	L261-911



Product Name	Cat. No.	Product Name	Cat. No.	Product Name	Cat. No.
HIPK2 siRNA Set I	H04-911	IRR siRNA Set I	I07-911	LOXL2 siRNA Set I	L262-911
HIPK3 siRNA Set I	H05-911	IRS1 siRNA Set I	I40-911	LOXL3 siRNA Set I	L263-911
HIPK4 siRNA Set I	H06-911	ITK siRNA Set I	I13-911	LRRK2 siRNA Set I	L10-911
Histone H1 siRNA Set I	H10-911	IκB alpha siRNA Set I	I20-911	LTK siRNA Set I	L11-911
Histone H3 siRNA Set I	H12-911	IκB beta siRNA Set I	I21-911	LYN A siRNA Set I	L13-911
HO1 siRNA Set I	H20-911	JAK2 siRNA Set I	J02-911	LYN B siRNA Set I	L13-911B
HO2 siRNA Set I	H21-911	JAK3 siRNA Set I	J03-911	MAPKAPK2 siRNA Set I	M40-911
HPK1 siRNA Set I	M23-911	JMJD1C siRNA Set I	J48-911	MAPKAPK3 siRNA Set I	M41-911
HRAS1 siRNA Set I	R05-911	JNK1 siRNA Set I	M33-911	MAPKAPK5 siRNA Set I	M42-911
HSD17B10 siRNA Set I	H23-911	JNK2 siRNA Set I	M34-911	MARCKS siRNA Set I	M02-911
HSF1 siRNA Set I	H25-911	JNK3 siRNA Set I	M35-911	MARK1 siRNA Set I	M43-911
HSP10 siRNA Set I	H29-911	KAT2A (GCN5) siRNA Set I	K311-911	MARK2 siRNA Set I	M44-911
		KAT2B (PCAF) siRNA Set I	K311-911B	MARK3 siRNA Set I	M45-911
		KAT3B (EP300) siRNA Set I	K312-911B	MARK4 siRNA Set I	M46-911
		KAT4 siRNA Set I	K313-911	MDM2 siRNA Set I	M45-911B
		KAT5 (TIP60) siRNA Set I	K314-911	MDM4 siRNA Set I	M47-911
		KAT6B (MYST4) siRNA Set I	K315-911	MEK1 siRNA Set I	M02-911B
		KAT7 (MYST2) siRNA Set I	K316-911	MEK2 siRNA Set I	M03-911
		KAT8 (MYST1) siRNA Set I	K317-911	MEKK1 siRNA Set I	M09-911
		KAT9 (ELP3) siRNA Set I	K318-911	MEKK2 siRNA Set I	M10-911
		KDM1A siRNA Set I	K421-911	MEKK3 siRNA Set I	M11-911
		KDM2A siRNA Set I	K422-911	MEKK6 siRNA Set I	M14-911



Product Name	Cat. No.	Product Name	Cat. No.	Product Name	Cat. No.
MELK siRNA Set I	M50-911	NIK siRNA Set I	M22-911	PDE1A siRNA Set I	P89-911
MET siRNA Set I	M52-911	NIM1 siRNA Set I	N14-911	PDE1B siRNA Set I	P89-911B
MGMT siRNA Set I	M354-911	NLK siRNA Set I	N15-911	PDE1C siRNA Set I	P89-911C
MGRN1 siRNA Set I	M287-911	NUAK1 siRNA Set I	N19-911	PDE2A siRNA Set I	P90-911
MINK1 siRNA Set I	M53-911	NUAK2 siRNA Set I	N20-911	PDE3A siRNA Set I	P91-911B
MLCK siRNA Set I	M62-911	OPG siRNA Set I	O878-911	PDE3B siRNA Set I	P91-911C
MLK1 siRNA Set I	M17-911	OSR1 siRNA Set I	O09-911	PDE4A siRNA Set I	P92-911
MLK2 siRNA Set I	M18-911	p16INK4A siRNA Set I	C30-911B	PDE4B siRNA Set I	P92-911B
MLK3 siRNA Set I	M19-911	p18INK4C siRNA Set I	C32-911	PDE4C siRNA Set I	P92-911C
MLK4 siRNA Set I	M48-911	p19INK4D siRNA Set I	C33-911B	PDE4D siRNA Set I	P92-911D
MNK1 siRNA Set I	M54-911	p21CIP1 siRNA Set I	C27-911B	PDE5A siRNA Set I	P93-911
MNK2 siRNA Set I	M55-911	p27KIP1 siRNA Set I	C28-911B	PDE6A siRNA Set I	P94-911B
MO25 alpha siRNA Set I	M51-911B	p33ING1 siRNA Set I	I32-911	PDE6B siRNA Set I	P94-911C
Moesin siRNA Set I	M52-911B	p33ING2 siRNA Set I	I32-911B	PDE6C siRNA Set I	P94-911D
Mortalin siRNA Set I	M54-911B	p38 alpha siRNA Set I	M39-911	PDE7A siRNA Set I	P95-911
MOS siRNA Set I	M57-911	p38 beta siRNA Set I	M36-911	PDE7B siRNA Set I	P95-911B
MRCK alpha siRNA Set I	C27-911	p38 gamma siRNA Set I	M37-911	PDE8A siRNA Set I	P96-911
MRCK beta siRNA Set I	C28-911	p38 delta siRNA Set I	M38-911	PDE8B siRNA Set I	P96-911B
MSK1 siRNA Set I	R19-911	p53 siRNA Set I	P05-911	PDE9A siRNA Set I	P97-911
MSK2 siRNA Set I	R18-911	p57KIP2 siRNA Set I	C29-911B	PDE10A siRNA Set I	P98-911
MSSK1 siRNA Set I	S41-911	p63 siRNA Set I	P06-911B	PDE12 siRNA Set I	P100-911
MST1 siRNA Set I	S25-911	p70S6K siRNA Set I	R21-911	PDGF siRNA Set I	P841-911



Product Name	Cat. No.	Product Name	Cat. No.	Product Name	Cat. No.
MST3 siRNA Set I	S42-911	p70S6Kb siRNA Set I	R22-911	PDGFR alpha siRNA Set I	P12-911
MST4 siRNA Set I	M59-911	p70S6KL1 siRNA Set I	R25-911	PDGFR beta siRNA Set I	P13-911
mTOR siRNA Set I	F17-911	p73 alpha siRNA Set I	P08-911	PDHA1 siRNA Set I	P57-911
MUSK siRNA Set I	M60-911	p73 beta siRNA Set I	P08-911B	PDHK1 siRNA Set I	P57-911B
MYLK2 siRNA Set I	M63-911	p73 gamma siRNA Set I	P08-911C	PDHK2 siRNA Set I	P58-911
MYLK3 siRNA Set I	M64-911	PAD1 siRNA Set I	P312-911	PDHK3 siRNA Set I	P59-911
MYO3 beta siRNA Set I	M66-911	PAD2 siRNA Set I	P312-911B	PDHK4 siRNA Set I	P60-911
MYT1 siRNA Set I	M67-911	PAD3 siRNA Set I	P312-911C	PDK1 siRNA Set I	P14-911
N6AMT2 siRNA Set I	N355-911	PAD4 siRNA Set I	P312-911D	PEAK1 siRNA Set I	P18-911
NCK siRNA Set I	N06-911	PAD6 siRNA Set I	P312-911E	PFKFB1 siRNA Set I	P321-911
NDR siRNA Set I	S49-911	PAK1 siRNA Set I	P02-911	PFKFB3 siRNA Set I	P323-911
NDR2 siRNA Set I	S50-911	PAK2 siRNA Set I	P03-911	PFKFB4 siRNA Set I	P324-911
NEK1 siRNA Set I	N02-911	PAK3 siRNA Set I	P04-911	PFTK1 (CDK14) siRNA Set I	P15-911
NEK2 siRNA Set I	N03-911	PAK4 siRNA Set I	P05-911B	PHKG1 siRNA Set I	P19-911
NEK3 siRNA Set I	N05-911	PAK6 siRNA Set I	P06-911C	PHKG2 siRNA Set I	P20-911
NEK5 siRNA Set I	N07-911	PAK7 siRNA Set I	P07-911B	PHLPP2 siRNA Set I	P72-911
NEK6 siRNA Set I	N08-911	PARK7 (DJ-1) siRNA Set I	P219-911	PI3K (p55 gamma) siRNA Set I	P31-911B
NEK7 siRNA Set I	N09-911	Parkin siRNA Set I	P215-911	PI3K (p85 alpha) siRNA Set I	P31-911C
NEK9 siRNA Set I	N11-911	PASK siRNA Set I	P08-911D	PI3K (p85 beta) siRNA Set I	P31-911D
NEK11 siRNA Set I	N13-911	PCNA siRNA Set I	P11-911	PI3K (p110 alpha) siRNA Set I	P27-911
NFATC1 siRNA Set I	N12-911	PCTK1 siRNA Set I	P09-911B	PI3K (p110 beta) siRNA Set I	P28-911
NFKB2 siRNA Set I	N13-911B	PCTK3 siRNA Set I	P11-911B	PI3K (p110 delta) siRNA Set I	P30-911

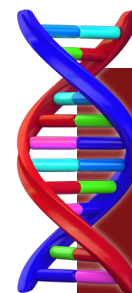
Product Name	Cat. No.	Product Name	Cat. No.	Product Name	Cat. No.
PI3K (p120 gamma) siRNA Set I	P29-911B	PP2C gamma siRNA Set I	P07-911C	RSK2 siRNA Set I	R17-911
PI4K2A siRNA Set I	P21-911	PRKAR1A siRNA Set I	P57-911C	RSK3 siRNA Set I	R16-911
PI4K2B siRNA Set I	P22-911B	PRKAR1B siRNA Set I	P59-911B	RSK4 siRNA Set I	R20-911
PI4KA siRNA Set I	P31-911	PRKG1 siRNA Set I	P78-911	RYK siRNA Set I	R28-911
PI4KB siRNA Set I	P32-911	PRKG2 siRNA Set I	P79-911	SCYL1 siRNA Set I	S04-911
PIAS1 siRNA Set I	P280-911	PRKX siRNA Set I	P81-911	SCYL3 siRNA Set I	S09-911
PIAS2 siRNA Set I	P281-911	PRKY siRNA Set I	P82-911	SDF-1 siRNA Set I	S821-911
PIAS3 siRNA Set I	P282-911	PRMT1 siRNA Set I	P365-911	SETD7 siRNA Set I	S344-911
PIAS4 siRNA Set I	P283-911	PRMT3 siRNA Set I	P365-911B	SETD8 siRNA Set I	S345-911
PIK3C2A siRNA Set I	P23-911	PRMT5 siRNA Set I	P365-911C	SETDB2 siRNA Set I	S346-911
PIK3C2B siRNA Set I	P24-911	PRPF19 siRNA Set I	P270-911	SETMAR siRNA Set I	S347-911
PIK3C2G siRNA Set I	P25-911	PSKH1 siRNA Set I	P89-911D	SGK1 siRNA Set I	S06-911
PIK3C3 (Vps34) siRNA Set I	P26-911	PTEN siRNA Set I	P23-911B	SGK2 siRNA Set I	S07-911
PIM1 siRNA Set I	P35-911	PTK7 siRNA Set I	P95-911C	SGK3 siRNA Set I	S08-911
PIM2 siRNA Set I	P36-911	PTK9 siRNA Set I	P97-911B	SH2D1A (SAP) siRNA Set I	S03-911
PIM3 siRNA Set I	P37-911	PTPN1 siRNA Set I	P28-911C	SIK siRNA Set I	S14-911
PIP4K2A siRNA Set I	P76-911	PTPN2 siRNA Set I	P29-911	SIK3 siRNA Set I	S12-911
PIP4K2B siRNA Set I	P76-911B	PTPN6 siRNA Set I	P33-911	SIRT1 siRNA Set I	S35-911
PIP4K2C siRNA Set I	P76-911C	PTPN7 siRNA Set I	P34-911	SIRT2 siRNA Set I	S36-911
PIP5K1A siRNA Set I	P16-911	PTPN11 siRNA Set I	P38-911E	SIRT3 siRNA Set I	S37-911
PIP5K1B siRNA Set I	P16-911B	PTPN12 siRNA Set I	P39-911	SIRT4 siRNA Set I	S38-911
PIP5K1C siRNA Set I	P16-911C	PTPN13 siRNA Set I	P40-911B	SIRT5 siRNA Set I	S39-911
PIP5K3 siRNA Set I	P17-911	PTPRA siRNA Set I	P48-911C	SIRT6 siRNA Set I	S40-911
PKAc alpha siRNA Set I	P51-911	PTPRC (CD45) siRNA Set I	P50-911B	SIRT7 siRNA Set I	S41-911B
PKAc beta siRNA Set I	P52-911	PTPRE siRNA Set I	P52-911B	SLK siRNA Set I	S11-911
PKAc gamma siRNA Set I	P53-911	PTPRF siRNA Set I	P53-911B	SMAD1 siRNA Set I	S10-911
PKC alpha siRNA Set I	P61-911	PTPRM siRNA Set I	P58-911B	SMAD2 siRNA Set I	S11-911B
PKC beta I siRNA Set I	P62-911	PTPRS siRNA Set I	P64-911B	SMAD3 siRNA Set I	S12-911B
PKC beta II siRNA Set I	P63-911	PYK2 siRNA Set I	P92-911E	SMAD4 siRNA Set I	S13-911
PKC delta siRNA Set I	P64-911	QIK siRNA Set I	S15-911	SMAD5 siRNA Set I	S14-911B
PKC epsilon siRNA Set I	P65-911	RAC1 siRNA Set I	R03-911B	SMAD9 siRNA Set I	S17-911
PKC eta siRNA Set I	P67-911	RAC2 siRNA Set I	R04-911	SMC1 siRNA Set I	S19-911
PKC gamma siRNA Set I	P66-911	RAC3 siRNA Set I	R02-911	SMYD2 siRNA Set I	S348-911
PKC iota siRNA Set I	P68-911	RAF1 (EE) siRNA Set I	R01-911	SMYD3 siRNA Set I	S348-911B
PKC mu siRNA Set I	P72-911B	RANTES (CCL5) siRNA Set I	R801-911	SNCA siRNA Set I	S22-911
PKC nu siRNA Set I	P73-911	Raptor siRNA Set I	R47-911	SNCB siRNA Set I	S22-911B
PKC theta siRNA Set I	P74-911	RAS siRNA Set I	R05-911B	SNCG siRNA Set I	S22-911C
PKC zeta siRNA Set I	P75-911	Rb siRNA Set I	R05-911C	SOD1 siRNA Set I	S26-911
PKD2 siRNA Set I	P76-911D	REL A siRNA Set I	R29-911	SOD2 siRNA Set I	S27-911
PKN2/PRK2 siRNA Set I	P71-911B	REL B siRNA Set I	R30-911	SOK1 siRNA Set I	S43-911
PKN3/PRK3 siRNA Set I	P69-911	RET siRNA Set I	R02-911B	SPHK1 siRNA Set I	S17-911B
PLAC1 siRNA Set I	P38-911	RGS1 siRNA Set I	R39-911	SPHK2 siRNA Set I	S18-911
PLAC2 siRNA Set I	P38-911B	RHEB siRNA Set I	R42-911	SRC siRNA Set I	S19-911B
PLAC3 siRNA Set I	P38-911C	RhoA siRNA Set I	R44-911	SRMS siRNA Set I	S20-911
PLAC4 siRNA Set I	P38-911D	RhoC siRNA Set I	R44-911B	SRPK1 siRNA Set I	S21-911
PLC gamma 1 siRNA Set I	P40-911	RhoD siRNA Set I	R44-911C	SRPK2 siRNA Set I	S22-911D
PLK1 siRNA Set I	P41-911	RIPK1 siRNA Set I	R07-911	STAT1 alpha siRNA Set I	S52-911
PLK2 siRNA Set I	P42-911	RIPK2 siRNA Set I	R08-911	STAT2 siRNA Set I	S53-911
PLK3 siRNA Set I	P43-911	RIPK3 siRNA Set I	R09-911	STAT3 siRNA Set I	S54-911
PLK4 siRNA Set I	P44-911	RIPK5 siRNA Set I	R27-911	STAT4 siRNA Set I	S55-911
POR siRNA Set I	P242-911	ROCK1 siRNA Set I	R10-911	STAT5 siRNA Set I	S56-911
PP1A siRNA Set I	P13-911B	ROCK2 siRNA Set I	R11-911	STAT5B siRNA Set I	S56-911B
PP1B siRNA Set I	P14-911B	RON siRNA Set I	M58-911	STAT6 siRNA Set I	S57-911
PP2A alpha siRNA Set I	P16-911D	ROR1 siRNA Set I	R12-911	Stathmin siRNA Set I	S71-911
PP2A beta siRNA Set I	P17-911B	ROR2 siRNA Set I	R13-911	STK3 siRNA Set I	S24-911
PP2B siRNA Set I	P18-911B	ROS1 siRNA Set I	R14-911	STK9 siRNA Set I	C46-911
PP2C alpha siRNA Set I	P02-911B	RSK1 siRNA Set I	R15-911	STK16 siRNA Set I	S31-911

Product Name	Cat. No.	Product Name	Cat. No.	Product Name	Cat. No.
STK19 siRNA Set I	S35-911B	TLK2 siRNA Set I	T11-911	UBE2F siRNA Set I	U220-911
STK21 (CIT) siRNA Set I	C52-911	TNF-alpha siRNA Set I	T860-911	UBE2G1 (UBC7) siRNA Set I	U221-911
STK31 siRNA Set I	S45-911	TNFB siRNA Set I	T861-911	UBE2H siRNA Set I	U223-911
STK32 (YANK2) siRNA Set I	Y04-911	TNFK siRNA Set I	T27-911	UBE2I siRNA Set I	U224-911
STK32C (YANK3) siRNA Set I	Y05-911	TOPK siRNA Set I	T14-911	UBE2J1 siRNA Set I	U225-911
STK33 siRNA Set I	S46-911	TP53RK siRNA Set I	T29-911	UBE2K (UBC1) siRNA Set I	U227-911
STK35 (CLIK1) siRNA Set I	S47-911	TPX2 siRNA Set I	T40-911	UBE2L6 (UBCH8) siRNA Set I	U229-911
STK36 siRNA Set I	S48-911	TRAF2 siRNA Set I	T57-911	UBE2N (UBC13) siRNA Set I	U231-911
STK39 (STLK3) siRNA Set I	S51-911B	TRAF3 siRNA Set I	T58-911	UBE2O siRNA Set I	U232-911
STK40 (SHIK) siRNA Set I	S54-911B	TRAIL siRNA Set I	T880-911	UBE2R1 siRNA Set I	U234-911
STUB1 (CHIP) siRNA Set I	C268-911	TRDMT1 (DNMT2) siRNA Set I	T352-911	UBE2R2 (UBC3B) siRNA Set I	U235-911
STYK1 (NOK) siRNA Set I	S53-911B	TRIM32 siRNA Set I	T291-911	UBE2U siRNA Set I	U237-911
SUMO1 siRNA Set I	S293-911	TRIM37 siRNA Set I	T292-911	UBE2V1 (UEV1A) siRNA Set I	U238-911
SUMO2 siRNA Set I	S294-911	TRKA siRNA Set I	N16-911	UBE2V2 siRNA Set I	U239-911
SUMO3 siRNA Set I	S295-911	TRKB siRNA Set I	N17-911	UBE2W (UBC16) siRNA Set I	U240-911
SUV39H1 siRNA Set I	S350-911	TRKC siRNA Set I	N18-911	UBE2Z (USE1) siRNA Set I	U241-911
SYK siRNA Set I	S52-911B	Troponin1 siRNA Set I	T68-911	UBLE1A (SAE1) siRNA Set I	U208-911
TAB1 siRNA Set I	T01-911	TSC1 siRNA Set I	T71-911	UCHL1 siRNA Set I	U22-911
TAK1 siRNA Set I	M15-911	TSC2 siRNA Set I	T72-911	UCHL3 siRNA Set I	U24-911
TAOK1 siRNA Set I	T24-911	TSSK1B siRNA Set I	S40-911B	UCHL5 siRNA Set I	U26-911
TAOK2 siRNA Set I	T25-911	TSSK2 siRNA Set I	S38-911B	UHMK1 (KIS) siRNA Set I	K05-911
TAOK3 siRNA Set I	T26-911	TSSK3 siRNA Set I	S39-911B	ULK1 siRNA Set I	U01-911
Tau-316 siRNA Set I	T02-911	TTBK1 siRNA Set I	T17-911	ULK2 siRNA Set I	U02-911
Tau-352 siRNA Set I	T03-911	TTBK2 siRNA Set I	T18-911	ULK3 siRNA Set I	U03-911
Tau-381 siRNA Set I	T04-911	TTK siRNA Set I	T20-911	ULK4 siRNA Set I	U04-911
Tau-383 siRNA Set I	T05-911	TUBB1 siRNA Set I	T92-911	VASP siRNA Set I	V12-911
Tau-410 siRNA Set I	T06-911	Tubulin alpha siRNA Set I	T91-911	VAV1 siRNA Set I	V14-911
Tau-412 siRNA Set I	T07-911	TXK siRNA Set I	T19-911	VEGF-C siRNA Set I	V812-911
Tau-441 siRNA Set I	T08-911	TYK2 (JTK1) siRNA Set I	T21-911	Vimentin siRNA Set I	V49-911
TBK1 siRNA Set I	T02-911B	TYRO3 siRNA Set I	T22-911	VLK siRNA Set I	V05-911
TEC siRNA Set I	T03-911B	Ubiquitin siRNA Set I	U06-911	VRK1 siRNA Set I	V01-911
TESK1 siRNA Set I	T05-911B	UBA1 (UBE1) siRNA Set I	U201-911	VRK2 siRNA Set I	V02-911
TESK2 siRNA Set I	T06-911B	UBA2 (SAE2) siRNA Set I	U202-911	VRK3 siRNA Set I	V03-911
TET1 siRNA Set I	T636-911	UBA3 (UBE1C) siRNA Set I	U203-911	WEE1 siRNA Set I	W01-911
TET2 siRNA Set I	T637-911	UBA7 (UBE1L) siRNA Set I	U207-911	WNK1 siRNA Set I	W02-911
TET3 siRNA Set I	T638-911	UBE2A siRNA Set I	U210-911	WWP1 siRNA Set I	W296-911
TGFB1 siRNA Set I	T831-911	UBE2B siRNA Set I	U211-911	WWP2 siRNA Set I	W297-911
TGFB3 siRNA Set I	T833-911	UBE2C siRNA Set I	U212-911	XIAP siRNA Set I	X08-911
TGFBR1 siRNA Set I	T07-911B	UBE2D1 (UBCH5A) siRNA Set I	U213-911	YES1 siRNA Set I	Y01-911
TGFBR2 siRNA Set I	T08-911B	UBE2D2 (UBC4) siRNA Set I	U214-911	ZAK siRNA Set I	Z01-911
TIE 2 siRNA Set I	T04-911B	UBE2D3 (UBCH5C) siRNA Set I	U215-911	ZAP70 siRNA Set I	Z02-911
TLK1 siRNA Set I	T10-911	UBE2E3 (UBCH9) siRNA Set I	U219-911	ZIP11 siRNA Set I	Z45-911

siRNA Controls

Product Name	Cat. No.	Product Name	Cat. No.
beta-Actin Positive Control siRNA	A03-911D	GAPDH Positive Control siRNA	G13-911B
Cy3 Labeled Negative Control siRNA	S02-910N	Unlabeled Negative Control siRNA	S01-910N
FAM Labeled Negative Control siRNA	S03-910N		

Visit our website for more product offerings



Tel: 1.604.232.4600 or
1.866.9KINASE (54.6273)
Fax: 1.604.232.4601
orders@signalchem.com
or order directly at:
<https://www.signalchem.com/ordering>



www.signalchem.com