

# SPECIALIZED ANTIBODIES

## Tag Antibodies

Product Name	Catalog #
Anti-DYKDDDDK	F51-61M
Anti-DYKDDDDK	F51-63R
Anti-GFP	G46-60M
Anti-GST	G52-61M

Product Name	Catalog #
Anti-HA	H98-61M
Anti-HA	H98-63R
Anti-His	H99-61M
Anti-His	H99-63R

Product Name	Catalog #
Anti-Myc	M86-61M
Anti-RFP	R46-61M
Anti-V5	V05-61M

## Phospho Specific Antibodies

Product Name	Catalog #
Anti-phospho-14-3-3 beta (Ser58)	Y71-365R
Anti-phospho-4EBP1 (Thr45)	E35-365R
Anti-phospho-AKT1/2/3 (Thr308)	A16-65R
Anti-phospho-AKT1/2/3 (Ser473)	A16-65BR
Anti-phospho-ASK1 (Ser966)	M13-65R
Anti-phospho-ATF2 (Ser112)	A10-565R
Anti-phospho-ATF2 (Ser490 Ser498)	A10-565BR
Anti-phospho-BRCA1 (Ser1423)	B87-365R
Anti-phospho-CAMK2A/2B (Thr286)	C11-65R
Anti-phospho-Catenin beta (Ser33 Ser37)	C06-365R
Anti-phospho-CDK1 (Tyr15)	C22-65R
Anti-phospho-CHK2 (Thr68)	C48-65R
Anti-phospho-CREB1 (Ser133)	C49-565R
Anti-phospho-EGFR (Thr678)	E10-65BR
Anti-phospho-EGFR (Thr693)	E10-65CR
Anti-phospho-EGFR (Tyr1092)	E10-65DR
Anti-phospho-EGFR (Tyr1110)	E10-65ER
Anti-phospho-EGFR (Tyr1197)	E10-65FR
Anti-phospho-EIF2AK2 (Thr446)	P80-65R
Anti-phospho-ELK1 (Ser383)	E64-65R
Anti-phospho-EPHB2 (Tyr469)	E22-65R
Anti-phospho-EPHB2 (Tyr473)	E22-65BR
Anti-phospho-ERK1/2 (Thr202)	M29-65R
Anti-phospho-ERK1/2 (Thr202/Tyr204)	M29-652R
Anti-phospho-ERK1/2 (Tyr204)	M29-65BR
Anti-phospho-GSK3 beta (Ser9)	G09-65R

Product Name	Catalog #
Anti-phospho-GSK3 beta (Ser9)	G09-652R
Anti-phospho-HER2 (Tyr877)	E27-65R
Anti-phospho-IRS1 (Ser312)	I40-365R
Anti-phospho-IRS1 (Ser639)	I40-365BR
Anti-phospho-JAK1 (Tyr1022)	J01-65R
Anti-phospho-KDR (Tyr951)	K01-65R
Anti-phospho-KDR (Tyr1214)	K01-65BR
Anti-phospho-MARCKS (Ser152 Ser156)	M02-565R
Anti-phospho-MEK1 (Thr292)	M02-65BR
Anti-phospho-MEK1 (Thr386)	M02-65CR
Anti-phospho-MEK1/2 (Ser218 Ser222)	M02-652R
Anti-phospho-MEK1/2 (Ser222)	M02-65R
Anti-phospho-MEK5 (Ser311 Ser315)	M06-65R
Anti-phospho-p38 alpha (Thr180 Tyr182)	M39-652R
Anti-phospho-p38 alpha (Tyr182)	M39-65R
Anti-phospho-p53 (Ser392)	P05-363R
Anti-phospho-PAK1 (Thr212)	P02-65BR
Anti-phospho-PAK1/2/3 (Thr402)	P02-652R
Anti-phospho-Parkin (Ser101)	P215-65R
Anti-phospho-Parkin (Ser378)	P215-65BR
Anti-phospho-PDK1 (Ser241)	P14-65R
Anti-phospho-PLCG1 (Tyr783)	P40-465R
Anti-phospho-PLCG2 (Tyr753)	P41-465R
Anti-phospho-PLK1 (Thr210)	P41-65R
Anti-phospho-RAF1 (Ser301)	R01-65R
Anti-phospho-RAF1 (Ser642)	R01-65BR

# Antibodies

## Phospho Specific Antibodies

Product Name	Catalog #
Anti-phospho-SNCA (Ser129)	S22-365R
Anti-phospho-STAT3 (Tyr705)	S54-365R
Anti-phospho-STAT3 (Ser727)	S54-365BR
Anti-phospho-STAT5A (Ser780)	S56-365R
Anti-phospho-STK10-a (Ser482 Ser486 Ser490)	S29-165R
Anti-phospho-Tau (Ser214)	T08-365GR
Anti-phospho-Tau (Ser235)	T08-365FR
Anti-phospho-Tau (Ser262)	T08-365IR
Anti-phospho-Tau (Ser356)	T08-365MR

Product Name	Catalog #
Anti-phospho-Tau (Ser396)	T08-365ER
Anti-phospho-Tau (Ser404)	T08-365HR
Anti-phospho-Tau (Ser416)	T08-365DR
Anti-phospho-Tau (Ser422)	T08-365KR
Anti-phospho-Tau (Thr181)	T08-365R
Anti-phospho-Tau (Thr205)	T08-365BR
Anti-phospho-Tau (Thr212)	T08-365JR
Anti-phospho-Tau (Thr231)	T08-365CR

## Modification Antibodies

Product Name	Catalog #
Anti-5-Methylated Cytosine (5-methyl cytosine)	C92-67S
Anti-Acetylated Lysine	L95-66R
Anti-Acetylated Lysine, Biotin Conjugated	L95-66BR
Anti-Acetylated Lysine, FITC Conjugated	L95-66CR
Anti-Acetylated Lysine, HRP Conjugated	L95-66DR
Anti-Methylated Arginine (NG,NG-Di-methyl)	A96-67R
Anti-Methylated Arginine (NG-Mono-methyl)	A95-67R
Anti-Methylated Lysine (Tri-epsilon-N-methyl)	L95-67R
Anti-Methylated Lysine (Tri-epsilon-N-methyl), Biotin Conjugated	L95-67BR
Anti-Methylated Lysine (Tri-epsilon-N-methyl), FITC Conjugated	L95-67CR

Product Name	Catalog #
Anti-Methylated Lysine (Tri-epsilon-N-methyl), HRP Conjugated	L95-67DR
Anti-Phosphoserine (Anti-pS)	S95-65R
Anti-Phosphoserine (Anti-pS), Biotin Conjugated	S95-65BR
Anti-Phosphothreonine (Anti-pT)	T95-65R
Anti-Phosphothreonine (Anti-pT), Biotin Conjugated	T95-65BR
Anti-Phosphotyrosine (Anti-pY)	T96-65R
Anti-Phosphotyrosine (Anti-pY), Biotin Conjugated	T96-65BR

## Secondary Antibodies

Product Name	Catalog #
Anti-Mouse HRP	G32-62G

Product Name	Catalog #
Anti-Rabbit HRP	G33-62G

### Signaling Related Antibodies



- » Characterized and Validated
- » Wide Range of Applications
- » Study Post-Translational Modifications

For additional details and antibody applications, please visit [signalchem.com](http://signalchem.com)

20140115