

# APOPTOSIS/AUTOPHAGY

www.signalchem.com

| Product Name                | Cat #    | Product Name                      | Cat #      | Product Name                | Cat #      |
|-----------------------------|----------|-----------------------------------|------------|-----------------------------|------------|
| AATK Protein                | A02-35G  | DNMT1 Protein                     | D351-31G   | JNK1, Unactive              | M33-14G    |
| AKT1, Active                | A16-10G  | DNMT3A, Active                    | D353-380G  | JNK2, Active                | M34-10BG   |
| AKT1 Mutant (E17K), Active  | A16-12G  | DNMT3A Protein                    | D353-30G   | JNK2, Unactive              | M34-14G    |
| AKT1, Unactive              | A16-14G  | DNMT3B, Active                    | D353-380BG | JNK3, Active                | M35-10BG   |
| AKT2, Active                | A17-10G  | DNMT3B Protein                    | D353-30BG  | JNK3, Unactive              | M35-14G    |
| AKT2, Active                | A17-10H  | DNMT3L, Active                    | D353-380CG | KAT2A (GCN5), Active        | K311-381G  |
| AKT2 Mutant (E17K), Active  | A17-12G  | DNMT3L Protein                    | D353-30CG  | KAT2B (PCAF), Active        | K311-381BG |
| AKT2 Mutant (R274H), Active | A17-12BG | DRAK1 (STK17A), Active            | S33-10G    | KAT3A (CREBBP), Active      | K312-381G  |
| AKT2, Unactive              | A17-14G  | DRAK2 (STK17B), Active            | S34-10G    | KAT3A (CREBBP) Protein      | K312-31G   |
| AKT2, Unactive              | A17-15H  | DUSP22 (MKPX), Active             | D22-20G    | KAT3B (EP300) Protein       | K312-31BG  |
| AKT3, Active                | A18-10G  | DYRK2, Active                     | D10-10G    | KAT4 Protein                | K313-31G   |
| AKT3 Mutant (G171R), Active | A18-12BG | DYRK3, Active                     | D11-11G    | KAT5 (TIP60), Active        | K314-380G  |
| AKT3, Unactive              | A18-14G  | EEF2K, Active                     | E01-10G    | KAT6B (MYST4), Active       | K315-381BG |
| AMPK (A1/B1/G1), Active     | P47-10H  | EGFR, Active                      | E10-11G    | KAT7 (MYST1), Active        | K316-380G  |
| AMPK (A1/B1/G2), Active     | P55-10H  | EGFR Mutant (T790M), Active       | E10-12EG   | KAT8 (MYST1) Protein        | K317-31G   |
| AMPK (A1/B1/G3), Active     | P56-10H  | EGFR Mutant (T790M L858R), Active | E10-12DG   | KAT9 (ELP3) Protein         | K318-30G   |
| AMPK (A1/B2/G1), Active     | P50-10H  | EGFR Mutant (L858R), Active       | E10-12BG   | KDR, Active                 | K01-11G    |
| AMPK (A1/B2/G2), Active     | P45-10G  | EGFR Mutant (L861Q), Active       | E10-12CG   | KSR1, Active                | K07-11G    |
| AMPK (A1/B2/G3), Active     | P83-10G  | EHMT2, Active                     | E318-381G  | KSR2, Active                | K08-11G    |
| AMPK (A2/B1/G1), Active     | P48-10H  | ELK1 Protein                      | E64-30G    | KSR2 Mutant (R676S), Active | K08-12BG   |
| AMPK (A2/B1/G2), Unactive   | P48-14H  | ENPP2 Protein                     | E22-35G    | LATS1, Active               | L01-11G    |
| AMPK (A2/B1/G2), Active     | P84-10GH | ERK1, Active                      | M29-10G    | LATS2, Active               | L02-11G    |
| AMPK (A2/B1/G3), Active     | P85-10GH | ERK1, Active                      | M29-10U    | LATS2, Unactive             | L02-35G    |
| AMPK (A2/B2/G1), Active     | P49-10H  | ERK1, Unactive                    | M29-14G    | LKB1 (STK11) Protein        | L15-30G    |
| AMPK (A2/B2/G2), Active     | P54-10H  | ERK1, Unactive                    | M29-14U    | MAPKAPK2, Active            | M40-11G    |
| AMPK (A2/B2/G3), Active     | P46-10GH | ERK1 (K71A) Protein               | M29-16H    | MAPKAPK2, Active            | M40-11H    |
| AMPK (A2/B2/G3), Unactive   | P46-14GH | ERK2, Active                      | M28-10G    | MAPKAPK2, Unactive          | M40-14G    |
| Annexin 5 Protein           | A61-30H  | ERK2, Unactive                    | M28-14G    | MAPKAPK3, Active            | M41-10G    |
| ASK1, Active                | M13-11G  | ERK2, Unactive                    | M28-14U    | MAPKAPK3, Unactive          | M41-14G    |
| ATF1 Protein                | A09-54G  | ERK5, Active                      | M32-10G    | MAPKAPK5, Active            | M42-10G    |
| ATF2 Protein                | A10-55G  | FADD Protein                      | F02-30H    | MDM2 Protein                | M45-31G    |
| ATG13 Protein               | A113-30G | FANCL Protein                     | F289-30G   | MDM2 (1-118) Protein        | M45-31BH   |
| BAD Protein                 | B03-30G  | FASTK, Active                     | F01-10G    | MDM4 (1-137) Protein        | M47-31BH   |
| BAG1 Protein                | B05-31G  | FOS Protein                       | F66-30G    | MEK1, Active                | M02-10G    |
| BAX Protein                 | B14-30G  | GAB1 Protein                      | G03-30H    | MEK1 (EE), Active           | M02-12G    |
| BCL2 Protein                | B25-30G  | GFP Protein                       | E620-30H   | MEK1, Unactive              | M02-14BG   |
| BID Protein                 | B59-30G  | GRB2 Protein                      | G45-30G    | MEK1, Unactive              | M02-14H    |
| BIRC3 Protein               | B280-30G | GRB7 Protein                      | G45-30GG   | MEK1 (K97A), Unactive       | M02-16H    |
| BLK, Active                 | B02-10G  | GSK3 alpha, Active                | G08-10G    | MEK2, Active                | M03-10G    |
| BRAF, Active                | B08-11G  | GSK3 beta, Active                 | G09-10G    | MEK2, Unactive              | M03-14G    |
| BRAF Mutant (V599E), Active | B08-13G  | GSK3 beta, Active                 | G09-10H    | MEKK1, Active               | M09-11G    |
| BRSK1, Active               | B13-10G  | GSK3 beta, Unactive               | G09-14G    | MEKK2, Active               | M10-10G    |
| BRSK2, Active               | B14-10G  | HGK, Active                       | M26-11G    | MEKK3, Active               | M11-10G    |
| c-JUN Protein               | J05-55G  | HRAS1 (Mature form) Protein       | R05-31H    | MEKK6, Active               | M14-11G    |
| CBP Protein                 | C07-31G  | HRAS1 (G12V) Protein              | R05-32H    | MGMT Protein                | M354-30H   |
| CREB1 Protein               | C49-55G  | HSP70 Protein                     | H34-54G    | MGRN1 Protein               | M287-30G   |
| CRKL Protein                | C53-30H  | HSP70 Protein                     | H34-54H    | MLK1, Active                | M17-11G    |
| DAPK1, Active               | D01-11G  | HSP90 alpha Protein               | H36-50H    | MLK2, Active                | M18-11G    |
| DAPK2, Active               | D02-10G  | HSP90 beta Protein                | H36-54H    | MLK3, Active                | M19-11G    |
| DAPK3, Active               | D03-10G  | IGF1R, Active                     | I02-11H    | MLK4, Active                | M48-11G    |
| DDR1, Active                | D05-11G  | IkBA Protein                      | I20-30G    | Mortalin Protein            | M54-31G    |
| DDR2, Active                | D06-11BG | IkBA Protein                      | I20-31G    | MST1, Active                | S25-10G    |
| DDR2 Mutant (T654M), Active | D06-12G  | IkBB Protein                      | I21-30G    | MST3, Active                | S42-11G    |
| DLK Protein                 | M20-31G  | JNK1, Active                      | M33-10G    | MST4, Active                | M59-10G    |

# APOPTOSIS/AUTOPHAGY

| Product Name                                      | Cat #     |
|---|-----------|
| NGAMT2 Protein                                    | N355-30BH |
| NDR, Active                                       | S49-10G   |
| NDR, Unactive                                     | S49-14G   |
| NDR2, Active                                      | S50-10G   |
| NFKB2 Protein                                     | N13-31G   |
| NUAK1, Active                                     | N19-10G   |
| NUAK2, Active                                     | N20-10G   |
| P300 Protein                                      | P07-31G   |
| p38 alpha, Active                                 | M39-10BG  |
| p38 alpha (K53A), Unactive                        | M39-16H   |
| p38 alpha, Unactive                               | M39-14G   |
| p38 alpha, Unactive                               | M39-14U   |
| p38 beta, Active                                  | M36-10BG  |
| p38 beta, Unactive                                | M36-14G   |
| p38 delta, Active                                 | M38-10BG  |
| p38 delta, Unactive                               | M38-14G   |
| p38 gamma, Active                                 | M37-10BG  |
| p38 gamma, Unactive                               | M37-14G   |
| p53 Protein                                       | P05-30BG  |
| p53 Protein                                       | P05-30G   |
| p63 Protein                                       | P06-30G   |
| p70S6K, Active                                    | R21-10H   |
| p70S6K, Unactive                                  | R21-14H   |
| p70S6Kb, Active                                   | R22-10G   |
| p70S6Kb, Unactive                                 | R22-14G   |
| p70S6K11 Protein                                  | R25-34G   |
| p73 alpha Protein                                 | P08-30G   |
| p73 beta Protein                                  | P08-30BG  |
| p73 gamma Protein                                 | P08-30CG  |
| PARK7 (DJ-1) Protein                              | P219-31H  |
| Parkin Protein                                    | P215-30G  |
| PDHA1 Protein                                     | P57-55H   |
| PDK1, Active                                      | P14-10H   |
| PI3K (p55 gamma) Protein                          | P31-30CH  |
| PI3K (p65 alpha) Protein                          | P31-31H   |
| PI3K (p85 alpha) Protein                          | P31-30H   |
| PI3K (p85 beta) Protein                           | P31-30BH  |
| PI3K (p110 alpha/p55 gamma), Active               | P27-10CH  |
| PI3K (p110 alpha/p85 alpha), Active               | P27-10H   |
| PI3K (p110 alpha/p85 alpha), Active               | P27-18H   |
| PI3K (p110 alpha(E545K)/p85 alpha) Mutant, Active | P27-15H   |
| PI3K (p110 beta/p85 alpha), Active                | P28-10H   |
| PI3K (p110 beta/p85 beta), Active                 | P28-10BH  |
| PI3K (p110 delta/p85 alpha), Active               | P30-10H   |
| PI3K (p120 gamma), Active                         | P29-10H   |
| PI4K2A, Active                                    | P21-10G   |
| PI4K2B, Active                                    | P22-10G   |
| PI4KA Protein                                     | P31-30G   |
| PI4KB, Active                                     | P32-10G   |
| PIK3C2A, Active                                   | P23-11G   |
| PIK3C2B Protein                                   | P24-31BG  |
| PIK3C2G, Active                                   | P25-11G   |
| PIK3C3 (Vps34), Active                            | P26-10G   |
| PIM1, Active                                      | P35-10G   |
| PIM2, Active                                      | P36-10G   |
| PIM3, Active                                      | P37-10G   |
| PIP4K2A, Active                                   | P76-10AG  |
| PIP4K2B Protein                                   | P76-30BG  |
| PIP4K2C, Active                                   | P76-10CG  |

| Product Name                | Cat #      |
|-----------------------------|------------|
| PIP5K1A, Active             | P16-10AG   |
| PIP5K1B, Active             | P16-10BG   |
| PIP5K1C, Active             | P16-10CG   |
| PIP5K3 Protein              | P17-31G    |
| PKAc alpha, Active          | P51-10G    |
| PKAc beta, Active           | P52-10G    |
| PKAc gamma, Active          | P53-10G    |
| PKC alpha, Active           | P61-10G    |
| PKC alpha, Active           | P61-18G    |
| PKC beta I, Active          | P62-10G    |
| PKC beta II, Active         | P63-10G    |
| PKC delta, Active           | P64-10G    |
| PKC epsilon, Active         | P65-10G    |
| PKC eta, Active             | P67-10G    |
| PKC gamma, Active           | P66-10G    |
| PKC iota, Active            | P68-10G    |
| PKC mu, Active              | P72-10G    |
| PKC nu, Active              | P73-10G    |
| PKC theta, Active           | P74-10G    |
| PKC zeta, Active            | P75-10G    |
| PKD2, Active                | P76-10G    |
| PKD2 Mutant (G848E), Active | P76-12BG   |
| PKD2 Mutant (G870E), Active | P76-12CG   |
| PKN2/PRK2, Active           | P71-10G    |
| PKN3/PRK3, Active           | P69-10G    |
| POR Protein                 | P242-30G   |
| PRKAR1A Protein             | P57-30H    |
| PRKAR1B Protein             | P59-30H    |
| PRKX, Active                | P81-10G    |
| PRMT1, Active               | P365-380G  |
| PRMT3, Active               | P365-380CG |
| PRMT5 Protein               | P365-30EG  |
| PRPF19 Protein              | P270-30G   |
| PTEN, Active                | P23-20G    |
| PTEN Protein                | P23-34G    |
| RAF1 (EE), Active           | R01-13G    |
| RAS Protein                 | R05-30G    |
| Rel B Protein               | R30-30H    |
| RHEB Protein                | R42-30H    |
| RhoA Protein                | R44-30H    |
| RhoC Protein                | R44-30CH   |
| RhoD Protein                | R44-30DH   |
| RIPK1 Protein               | R07-34G    |
| RIPK2, Active               | R08-11G    |
| RIPK3, Active               | R09-10G    |
| RIPK5, Active               | R27-10G    |
| ROCK1, Active               | R10-11G    |
| ROCK2, Active               | R11-11G    |
| ROCK2, Active               | R11-11H    |
| RSK1, Active                | R15-10G    |
| RSK2, Active                | R17-10G    |
| RSK2 Mutant (I416V), Active | R17-12BG   |
| RSK2 Mutant (Y483C), Active | R17-12DG   |
| RSK2 Mutant (L608F), Active | R17-12CG   |
| RSK3, Active                | R16-10G    |
| RSK4, Active                | R20-10G    |
| SETD7, Active               | S344-380G  |

| Product Name           | Cat #     |
|------------------------|-----------|
| SETD8 Protein          | S345-30G  |
| SETDB2 Protein         | S346-30BG |
| SETMAR Protein         | S347-30G  |
| SGK1, Active           | S06-11G   |
| SGK1, Unactive         | S06-15G   |
| SGK2, Active           | S07-10G   |
| SGK2, Unactive         | S07-14G   |
| SGK3, Active           | S08-10G   |
| SGK3, Unactive         | S08-14G   |
| SMYD2 Protein          | S348-31BG |
| SOD1 Protein           | S26-30G   |
| SOD2 Protein           | S27-30G   |
| SOK1, Active           | S43-10G   |
| SOK1, Unactive         | S43-14G   |
| SPHK1, Active          | S17-10G   |
| SPHK2, Active          | S18-10G   |
| SRC, Active            | S19-10G   |
| SRC, Active            | S19-18G   |
| SRMS, Active           | S20-11G   |
| STAT1 alpha Protein    | S52-50G   |
| STAT1 beta Protein     | S52-54G   |
| STAT2 Protein          | S53-54G   |
| STAT3 Protein          | S54-54G   |
| STAT3 Protein          | S54-54H   |
| STAT4 Protein          | S55-54H   |
| STAT5 Protein          | S56-54H   |
| STAT5B Protein         | S56-54BG  |
| STAT6 Protein          | S57-30G   |
| STK3, Active           | S24-10G   |
| SUMO1 (UBL1) Protein   | S293-31H  |
| SUMO2 Protein          | S294-31H  |
| SUMO3 Protein          | S295-31H  |
| SUV39H1 Protein        | S350-30G  |
| TESK2, Active          | T06-10G   |
| TP53RK Protein         | T29-34G   |
| TPX2 Protein           | T40-30H   |
| TRAF2 Protein          | T57-30H   |
| TRDMT1 (DNMT2) Protein | T352-30G  |
| TRIM32 Protein         | T291-30G  |
| TRIM37 Protein         | T292-30G  |
| Ubiquitin Protein      | U06-54H   |
| UCLH1 Protein          | U22-30H   |
| UCLH3 Protein          | U24-30H   |
| UCLH5 Protein          | U26-30H   |
| ULK1, Active           | U01-11G   |
| ULK2, Active           | U02-11G   |
| ULK3, Active           | U03-10H   |
| ULK4 Protein           | U04-35G   |
| VAV1 Protein           | V14-31G   |
| Vimentin Protein       | V49-30H   |
| VRK1, Active           | V01-10G   |
| VRK2, Active           | V02-11G   |
| VRK3 Protein           | V03-30G   |
| XIAP Protein           | X08-30G   |
| YES1, Active           | Y01-10G   |
| ZAK, Active            | Z01-10G   |