

RPPAセルシグナル解析サービス 対象解析ターゲットリスト

2012年 12月 1日 更新

Angiogenesis	p-eNOS (Ser1177)
Apoptosis	p-Bad (Ser112)
Apoptosis	p-Bad (Ser136)
Apoptosis	p-Bad (Ser155)
Apoptosis	p-Bcl-2 (Ser70)
Apoptosis	p-Bcl-2 (Thr56)
Apoptosis	p-Bim (Ser69)
Apoptosis	p-NDRG1 (Thr346)
Cell cycle	p-Aurora A (Thr288)
Cell cycle	p-Aurora A (Thr288)/Aurora B (Thr232)/Aurora C (Thr198)
Cell cycle	p-Cyclin D1 (Thr286)
Cell cycle	p-Rb (Ser780)
Cell cycle	p-Rb (Ser795)
Cell cycle	p-Rb (Ser807/811)
Cell Cycle	p-Wee1 (Ser642)
Cytoskelton	p-AP2M1 (Thr156)
Cytoskelton	p-Cofilin (Ser3)
Cytoskelton	p-Ezrin (Thr567)/Radixin (Thr564)/Moesin (Thr558)
Cytoskelton	p-MARCKS (Ser167/170)
Cytoskelton	p-MYPT (Ser507)
Cytoskelton	p-MYPT (Ser668)
Cytoskelton	p-MYPT (Thr853)
Cytoskelton	p-VASP (Ser157)
Cytoskelton	p-VASP (Ser239)
DNA damage	p-ATM (Ser1981)
DNA damage	p-ATR (Ser428)
DNA damage	p-BAP1 (Ser592)
DNA damage	p-BRCA1 (Ser1524)
DNA damage	p-Chk1 (Ser296)
DNA damage	p-Chk1 (Ser317)
DNA damage	p-Chk1 (Ser345)
DNA damage	p-Chk2 (Ser19)
DNA damage	p-Chk2 (Ser33/35)
DNA damage	p-Chk2 (Thr68)
DNA damage	p-p53 (Ser15)
DNA damage	p-p53 (Ser20)
DNA damage	p-p53 (Ser37)
DNA damage	p-p53 (Ser392)
DNA damage	p-p53 (Ser46)
DNA damage	p-p53 (Ser6)
DNA damage	p-p53 (Ser9)
Energy homeostasis	p-Acetyl-CoA carboxylase (Ser79)
Energy homeostasis	p-AMPKα (Thr172)
Energy homeostasis	p-AMPKβ1 (Ser108)
Energy homeostasis	p-LKB1 (Ser428)
Forkhead	p-FoxO1 (Ser256)
Forkhead	p-FoxO1 (Thr24)/FoxO3a (Thr32)
Forkhead	p-FoxO1 (Thr24)/FoxO3a (Thr32)/FoxO4 (Thr28)
Forkhead	p-FoxO3a (Ser253)
Forkhead	p-FoxO3a (Ser318/321)
Insulin receptor	p-IRS-1(Ser307)
Insulin receptor	p-IRS-1(Ser612)
Insulin receptor	p-IRS-1(Ser636/639)
Mitosis marker	p-Histone H2A.X (Ser139)
Mitosis marker	p-Histone H3 (Ser10)
Mitosis marker	p-Histone H3 (Ser28)
Mitosis marker	p-Histone H3 (Thr11)
Mitosis marker	p-Histone H3 (Thr3)
Mitosis marker	p-NPM (Ser4)
Mitosis marker	p-TACC3 (Ser558)
NFκB	p-IκB-α (Ser32/36)
NFκB	p-NFκB p65 (Ser468)
NFκB	p-NFκB p65 (Ser536)
NFκB	p-RelB (Ser552)
NFκB	p-TBK1/NAK (Ser172)
p38 MAPK	p-HSP27 (Ser82)
p38 MAPK	p-MAPKAPK-2 (Thr222)
p38 MAPK	p-MAPKAPK-2 (Thr334)
p38 MAPK	p-MKK3/MKK6 (Ser189/207)
p38 MAPK	p-MSK1 (Thr581)
p38 MAPK	p-p38 MAPK (Thr180/Tyr182)
PI3K-AKT-mTOR	Non-p-4E-BP1 (Thr46)
PI3K-AKT-mTOR	p-4E-BP1 (Ser65)
PI3K-AKT-mTOR	p-4E-BP1 (Thr37/46)
PI3K-AKT-mTOR	p-4E-BP1 (Thr70)
PI3K-AKT-mTOR	p70-S6 kinase (Thr389)
PI3K-AKT-mTOR	p-Akt (Ser473)
PI3K-AKT-mTOR	p-Akt (Thr308)
PI3K-AKT-mTOR	p-eEF2k (Ser366)
PI3K-AKT-mTOR	p-eIF4B (Ser422)
PI3K-AKT-mTOR	p-eIF4E (Ser209)
PI3K-AKT-mTOR	p-eIF4G (Ser1108)
PI3K-AKT-mTOR	p-GSK-3α (Ser21)
PI3K-AKT-mTOR	p-GSK-3α/β (Ser21/9) GSK-3α preferred
PI3K-AKT-mTOR	p-GSK-3β (Ser9)
PI3K-AKT-mTOR	p-mTOR (Ser2448)
PI3K-AKT-mTOR	p-PDK1 (Ser241)
PI3K-AKT-mTOR	p-PKD/PKC γ (Ser744/748)
PI3K-AKT-mTOR	p-PKD/PKC γ (Ser916)
PI3K-AKT-mTOR	p-PRAS40 (Thr246)

PI3K-AKT-mTOR	p-PEN (Ser380)
PI3K-AKT-mTOR	p-Raptor (Ser792)
PI3K-AKT-mTOR	p-S6Rb (Ser235/236)
PI3K-AKT-mTOR	p-Tuberin/TSC2 (Ser939)
PI3K-AKT-mTOR	p-Tuberin/TSC2 (Thr1462)
PI3K-AKT-mTOR	p-Tuberin/TSC2 (Thr1571)
Ras-Raf	p-A-Raf (Ser299)
Ras-Raf	p-B-Raf (Ser445)
Ras-Raf	p-c-Raf (Ser259)
Ras-Raf	p-c-Raf (Ser289/296/301)
Ras-Raf	p-c-Raf (Ser338)
Ras-Raf	p-CREB (Ser133)
Ras-Raf	p-MEK1/2 (Ser217/221)
Ras-Raf	p-Mnk1 (Thr197/202)
Ras-Raf	p-p44/42 MAP kinase (Thr202/Tyr204)
Ras-Raf	p-p90RSK (Ser380)
Ras-Raf	p-p90RSK (Thr359/Ser363)
Ras-Raf	p-p90RSK (Thr573)
Ras-Raf	p-PAK1 (Ser144)/PAK2 (Ser141)
Ras-Raf	p-PAK1 (Ser199/204)/PAK2 (Ser182/197)
Ras-Raf	p-PAK1 (Thr423)/PAK2 (Thr402)
Ras-Raf	p-PAK2 (Ser20)
Ras-Raf	p-RSK2 (Ser227)
Ras-Raf	p-RSK3(Thr356/Ser360)
RTK	Non-p-Src (Tyr416)
RTK	Non-p-Src (Tyr527)
RTK	p-ALK (Tyr1282/1283)
RTK	p-c-Kit(Tyr703)
RTK	p-EGF Receptor (Tyr1045)
RTK	p-EGF Receptor (Tyr1068)
RTK	p-EGF Receptor (Tyr992)
RTK	p-FLT3 (Tyr591)
RTK	p-Gab1 (Tyr627)
RTK	p-Gab2 (Tyr452)
RTK	p-HER2/ErbB2 (Tyr1221/1222)
RTK	p-HER2/ErbB2 (Tyr1248)
RTK	p-HER4/ErbB4 (Tyr984)
RTK	p-IGF-I Receptor (Tyr1131)/Insulin Receptor (Tyr1146)
RTK	p-IGF-I Receptor (Tyr1135/1136)
RTK	p-IGF-I Receptor (Tyr980)
RTK	p-Met (Tyr1234/1235)
RTK	p-PDGF Receptor-β (Tyr1009)
RTK	p-PDGF Receptor-β (Tyr1021)
RTK	p-PDGF Receptor-β (Tyr740)
RTK	p-PDGF Receptor-β (Tyr751)
RTK	p-PDGF Receptor-β (Tyr771)
RTK	p-PKC(pan) βII (Ser660)
RTK	p-PKCα/β (Thr638/641)
RTK	p-PKCδ (Thr505)
RTK	p-PKCδ/θ (Ser643/676)
RTK	p-PKCζ/λ (Thr410/403)
RTK	p-PKCθ (Thr538)
RTK	p-PKM2 (Tyr105)
RTK	p-PLCγ (Tyr1217)
RTK	p-PLCγ (Tyr759)
RTK	p-PLCγ (Tyr783)
RTK	p-Ret (Tyr905)
RTK	p-Shc (Tyr239)
RTK	p-Shc (Tyr317)
RTK	p-SHIP2 (Tyr1135)
RTK	p-SHP-2 (Tyr580)
RTK	p-Src Family (Tyr416)
RTK	p-Src (Tyr527)
RTK	p-Syk (Tyr323)
RTK	p-Syk (Tyr525/526)
RTK	p-VEGF Receptor 2 (Tyr1059)
RTK	p-VEGF Receptor 2 (Tyr1175)
RTK	p-VEGF Receptor 2 (Tyr951)
RTK	p-Zap-70 (Tyr319)/Syk (Tyr352)
SAPK/JNK	p-ATF-2 (Thr71)
SAPK/JNK	p-c-Fos (Ser32)
SAPK/JNK	p-c-Jun (Ser63)
SAPK/JNK	p-c-Jun (Ser73)
SAPK/JNK	p-SEK1/MKK4 (Thr261)
SAPK/JNK	p-SAPK/JNK(Thr183/Tyr185)
STAT	p-Jak2 (Tyr1007/1008)
STAT	p-Stat 1(Tyr701)
STAT	p-Stat 2 (Tyr690)
STAT	p-Stat 3 (Ser727)
STAT	p-Stat 3 (Tyr705)
STAT	p-Stat 5 (Tyr694)
STAT	p-Stat 6 (Tyr641)
STAT	p-Tyk2 (Tyr1054/1055)
TGF-β	p-Smad1/5 (Ser463/465)
TGF-β	p-Smad2 (Ser465/467)
TGF-β	p-Smad2 (Ser465/467)/Smad3 (Ser423/425)
TGF-β	p-TAK1 (Thr184/187)
Wnt	p-LRP6 (Ser1490)
Wnt	p-β-catenin (Ser33/37/Thr41)
Wnt	p-β-catenin (Ser45)
Housekeeping	γ-tubulin