

Protein Kinase Catalytic Domain Conservation

Most Evolutionary Conserved Protein Kinases

| Kinase Name | Mouse | Rat | Dog | Chicken | Fly | Worm | Mustard Plant | Rice | Budding Yeast | Neurospora | SUM |
|------------------------|-------|-----|-----|---------|-----|------|---------------|------|---------------|------------|-----|
| Average Conserv. | 89 | 89 | 91 | 77 | 50 | 47 | 43 | 44 | 41 | 44 | 615 |
| CK2 α 1/CSNK2A1 | 98 | 98 | 100 | 97 | 88 | 78 | 75 | 76 | 59 | 70 | 839 |
| AMPK α 2/PRKAA2 | 97 | 98 | 98 | 90 | 60 | 58 | 46 | 46 | 42 | 80 | 715 |
| GSK3 β | 96 | 96 | 99 | 95 | 70 | 74 | 63 | | 54 | 62 | 709 |
| p38 α /MAPK14 | 95 | 99 | 95 | 96 | 68 | 62 | 45 | 46 | 52 | 49 | 707 |
| CDK7 | 93 | 95 | 92 | 94 | 66 | 60 | 57 | 58 | 48 | 42 | 705 |
| PKAC β /PRKACB | 97 | 97 | 88 | 90 | 83 | 80 | 35 | | 54 | 55 | 679 |
| CDK5 | 99 | 99 | 99 | 98 | 79 | 74 | | | 55 | 60 | 663 |
| CK1 δ /CSNK1D | 99 | 99 | 99 | 97 | 61 | | 74 | 61 | | 71 | 661 |
| ERK2/MAPK1 | 99 | 99 | 96 | 98 | 81 | 81 | 50 | | 50 | | 654 |
| CK1 α /CSNK1A1 | 92 | 91 | 100 | 99 | 70 | 79 | 59 | 60 | | | 650 |
| MAP2K1/MEK1 | 98 | 99 | 91 | 95 | 63 | 52 | 36 | 33 | 42 | 39 | 648 |
| AurA/STK6 | 82 | 83 | 89 | 71 | 50 | 57 | 57 | 53 | 43 | 55 | 640 |
| PAK1 | 97 | 98 | 96 | 94 | 51 | 53 | | 49 | 45 | 40 | 623 |
| FRAP/FRAP1 | 98 | 98 | 99 | 93 | 53 | 51 | 40 | 46 | 41 | | 619 |
| NDR2/STK38L | 97 | | 99 | 97 | 71 | 67 | 50 | 49 | 47 | 42 | 619 |
| CK1 ϵ /CSNK1E | 98 | 99 | 99 | 95 | 84 | | 63 | 74 | | | 612 |
| PDHK3/PDK3 | 97 | 97 | 99 | 92 | 58 | 53 | 38 | 39 | | 37 | 610 |
| PKAC α /PRKACA | 98 | 98 | 98 | 95 | 82 | 83 | | | 48 | | 602 |
| CLK2 | 96 | 96 | 98 | 88 | 58 | 62 | 49 | | | 46 | 593 |
| CDK11/CDC2L6 | 87 | 96 | 96 | 79 | 85 | 60 | 44 | 41 | | | 588 |
| CDK9 | 98 | 98 | 98 | 90 | 73 | 51 | | | 41 | 34 | 583 |
| NDR1/STK38 | 98 | 97 | 95 | 93 | 80 | 66 | 49 | | | | 578 |
| RIOK2 | 83 | 81 | 84 | 67 | 45 | 41 | 45 | 45 | 40 | 47 | 578 |
| MST2/STK3 | 94 | 96 | 98 | 98 | 46 | 54 | 43 | 43 | | | 572 |
| p70S6K/RPS6KB1 | 99 | 99 | 99 | 98 | 72 | 60 | 43 | | | | 570 |
| PKC ϵ /PRKCE | 97 | 98 | 98 | 82 | 57 | 57 | | | 46 | 33 | 568 |
| PIK3R4 | 95 | 95 | 97 | 91 | 41 | 29 | 30 | 26 | 26 | 30 | 560 |
| PRP4/PRPF4B | 95 | 95 | 98 | 90 | 43 | 41 | 33 | 48 | | | 543 |
| RIOK1 | 80 | 82 | 89 | 75 | 48 | 42 | 46 | | 38 | 41 | 541 |
| TRRAP | 98 | 98 | 98 | 93 | 49 | 24 | 27 | | 27 | 27 | 541 |
| CaMKK2 | 89 | 92 | 91 | 74 | 54 | 55 | 40 | | | 45 | 540 |
| CDK2 | 96 | 72 | 96 | | 70 | | | 71 | 64 | 64 | 533 |
| OSR1/OXSR1 | 95 | 95 | 97 | 86 | 56 | 52 | 49 | | | | 530 |
| PDK1/PDPK1 | 94 | 95 | 94 | 90 | 43 | 37 | 38 | 39 | | | 530 |
| BRAF | 99 | 87 | 93 | 89 | 45 | 40 | 33 | 42 | | | 528 |
| AMPK α 1/PRKAA1 | 96 | 96 | 98 | 83 | | 58 | 50 | 46 | | | 527 |
| JNK3/MAPK10 | 98 | 99 | 98 | 84 | 78 | 70 | | | | | 527 |
| AKT2 | 98 | 97 | 97 | | 61 | | 40 | 39 | 49 | 45 | 526 |
| NLK | 99 | 99 | 99 | 98 | 76 | 54 | | | | | 525 |
| CaMK2 δ | 92 | 92 | 91 | 68 | 74 | 67 | 39 | | | | 523 |

