

Use Promo Code NEWPROT50 to Save 50% On New Sample Size Recombinant Proteins

Product	Cat No	Size	Webpage
Recombinant ADAR1 (1-295) protein	82140	100 µg	https://www.activemotif.com/catalog/details/82140
Recombinant ADAR1 protein	82117	50 µg	https://www.activemotif.com/catalog/details/82117
Recombinant AHR / HIF1B Complex	82137	10 µg	https://www.activemotif.com/catalog/details/82137
Recombinant beta-Catenin protein	82067	20 µg	https://www.activemotif.com/catalog/details/82067
Recombinant Calmodulin-1 protein	82061	50 µg	https://www.activemotif.com/catalog/details/82061
Recombinant CDK1 / Cyclin B1 Complex	82134	20 µg	https://www.activemotif.com/catalog/details/82134
Recombinant CDK1 / Cyclin E1 Complex	82135	20 µg	https://www.activemotif.com/catalog/details/82135
Recombinant CDK1 protein	82082	20 µg	https://www.activemotif.com/catalog/details/82082
Recombinant CDK2 / Cyclin A2 Complex	82089	20 µg	https://www.activemotif.com/catalog/details/82089
Recombinant CDK2 / Cyclin B1 Complex	82136	20 µg	https://www.activemotif.com/catalog/details/82136
Recombinant CDK2 / Cyclin E1 Complex	82088	20 µg	https://www.activemotif.com/catalog/details/82088
Recombinant CDK2 protein	82083	20 µg	https://www.activemotif.com/catalog/details/82083
Recombinant CDK4 / Cyclin D1 Complex	82084	20 µg	https://www.activemotif.com/catalog/details/82084
Recombinant CDK4 / Cyclin D2 Complex	82085	20 µg	https://www.activemotif.com/catalog/details/82085
Recombinant CDK4 / Cyclin D3 Complex	82086	20 µg	https://www.activemotif.com/catalog/details/82086
Recombinant CDK6 / Cyclin D2 complex	82096	20 µg	https://www.activemotif.com/catalog/details/82096
Recombinant CDK6 / Cyclin D3 Complex	82087	20 µg	https://www.activemotif.com/catalog/details/82087
Recombinant CHD1L protein	82071	20 µg	https://www.activemotif.com/catalog/details/82071
Recombinant CHD2 (1-947) protein	82114	20 µg	https://www.activemotif.com/catalog/details/82114
Recombinant CHD6 (288-959) protein	82072	20 µg	https://www.activemotif.com/catalog/details/82072
Recombinant CHD7 protein	82065	10 µg	https://www.activemotif.com/catalog/details/82065
Recombinant CHD8 protein	82066	20 µg	https://www.activemotif.com/catalog/details/82066
Recombinant CLK1 protein	82109	20 µg	https://www.activemotif.com/catalog/details/82109
Recombinant CLK2 protein	82110	20 µg	https://www.activemotif.com/catalog/details/82110
Recombinant CLK3 protein	82111	20 µg	https://www.activemotif.com/catalog/details/82111
Recombinant CREB1 protein	82073	10 µg	https://www.activemotif.com/catalog/details/82073
Recombinant CRT1 protein	82075	10 µg	https://www.activemotif.com/catalog/details/82075
Recombinant CRT2 protein	82074	10 µg	https://www.activemotif.com/catalog/details/82074
Recombinant CTCF protein	82119	10 µg	https://www.activemotif.com/catalog/details/82119
Recombinant CUL5 protein	82103	50 µg	https://www.activemotif.com/catalog/details/82103
Recombinant DDX3X protein	82141	10 µg	https://www.activemotif.com/catalog/details/82141
Recombinant DDX5 protein	82125	20 µg	https://www.activemotif.com/catalog/details/82125

Recombinant DDX54 protein	82144	20 µg	https://www.activemotif.com/catalog/details/82144
Recombinant DNMT2 protein	82095	20 µg	https://www.activemotif.com/catalog/details/82095
Recombinant E2F1 protein	82076	20 µg	https://www.activemotif.com/catalog/details/82076
Recombinant EEF2 protein	82118	20 µg	https://www.activemotif.com/catalog/details/82118
Recombinant ELAVL1 protein	82152	50 µg	https://www.activemotif.com/catalog/details/82152
Recombinant FACT complex	82064	20 µg	https://www.activemotif.com/catalog/details/82064
Recombinant GLI1 protein	82129	10 µg	https://www.activemotif.com/catalog/details/82129
Recombinant HMGA2 protein	82151	50 µg	https://www.activemotif.com/catalog/details/82151
Recombinant IRAK4 (154-460) protein	82100	20 µg	https://www.activemotif.com/catalog/details/82100
Recombinant KRAS (1-186, A146T) protein	82060	100 µg	https://www.activemotif.com/catalog/details/82060
Recombinant KRAS (1-186, G12A) protein	82051	100 µg	https://www.activemotif.com/catalog/details/82051
Recombinant KRAS (1-186, G12C) protein	82052	100 µg	https://www.activemotif.com/catalog/details/82052
Recombinant KRAS (1-186, G12D) protein	82053	100 µg	https://www.activemotif.com/catalog/details/82053
Recombinant KRAS (1-186, G12S) protein	82054	100 µg	https://www.activemotif.com/catalog/details/82054
Recombinant KRAS (1-186, G12V) protein	82055	100 µg	https://www.activemotif.com/catalog/details/82055
Recombinant KRAS (1-186, G13C) protein	82056	100 µg	https://www.activemotif.com/catalog/details/82056
Recombinant KRAS (1-186, G13D) protein	82057	100 µg	https://www.activemotif.com/catalog/details/82057
Recombinant KRAS (1-186, G61R) protein	82059	100 µg	https://www.activemotif.com/catalog/details/82059
Recombinant KRAS (1-186, Q61H) protein	82058	100 µg	https://www.activemotif.com/catalog/details/82058
Recombinant MAPK1 (ERK2) protein	82081	20 µg	https://www.activemotif.com/catalog/details/82081
Recombinant MAPK8 (JNK1) protein	82062	100 µg	https://www.activemotif.com/catalog/details/82062
Recombinant MAPK9 (JNK2) protein	82063	100 µg	https://www.activemotif.com/catalog/details/82063
Recombinant MDM2 protein	82044	50 µg	https://www.activemotif.com/catalog/details/82044
Recombinant NEDD8 protein	82094	100 µg	https://www.activemotif.com/catalog/details/82094
Recombinant NFATC3 protein	82097	20 µg	https://www.activemotif.com/catalog/details/82097
Recombinant PADI4 protein	82098	20 µg	https://www.activemotif.com/catalog/details/82098
Recombinant PARP11 protein	82150	10 µg	https://www.activemotif.com/catalog/details/82150
Recombinant PARP7 protein	82149	10 µg	https://www.activemotif.com/catalog/details/82149
Recombinant PDPK1 protein, GST-Tag	82167	20 µg	https://www.activemotif.com/catalog/details/82167
Recombinant PIAS3 protein	82092	50 µg	https://www.activemotif.com/catalog/details/82092
Recombinant PPARα protein	82078	20 µg	https://www.activemotif.com/catalog/details/82078
Recombinant PRC2 Hexameric Complex	82180	10 mg	https://www.activemotif.com/catalog/details/82180
Recombinant PRDM16 protein	82040	20 µg	https://www.activemotif.com/catalog/details/82040
Recombinant PRKCA protein	82122	20 µg	https://www.activemotif.com/catalog/details/82122
Recombinant RANBP2 (2553-2838) protein	82101	100 µg	https://www.activemotif.com/catalog/details/82101

Recombinant RIPK1 (1-294) protein	82099	20 µg	https://www.activemotif.com/catalog/details/82099
Recombinant RNF4 protein	82003	100 µg	https://www.activemotif.com/catalog/details/82003
Recombinant SMAD1 protein	82069	20 µg	https://www.activemotif.com/catalog/details/82069
Recombinant SMAD2 protein	82070	20 µg	https://www.activemotif.com/catalog/details/82070
Recombinant SMARCA4 protein	82090	20 µg	https://www.activemotif.com/catalog/details/82090
Recombinant SOX9 protein	82079	20 µg	https://www.activemotif.com/catalog/details/82079
Recombinant SREBF1 (1-530) protein	82123	20 µg	https://www.activemotif.com/catalog/details/82123
Recombinant STAT1 protein	82080	20 µg	https://www.activemotif.com/catalog/details/82080
Recombinant STAT4 protein	82112	20 µg	https://www.activemotif.com/catalog/details/82112
Recombinant STAT5B protein	82113	20 µg	https://www.activemotif.com/catalog/details/82113
Recombinant STUB1 / CHIP protein	82008	100 µg	https://www.activemotif.com/catalog/details/82008
Recombinant SUMO1 protein - biotinylated	82105	100 µg	https://www.activemotif.com/catalog/details/82105
Recombinant SUMO2 protein - biotinylated	82106	100 µg	https://www.activemotif.com/catalog/details/82106
Recombinant SUMO3 protein - biotinylated	82107	100 µg	https://www.activemotif.com/catalog/details/82107
Recombinant TDG protein	82093	50 µg	https://www.activemotif.com/catalog/details/82093
Recombinant TOP2B protein	82115	20 µg	https://www.activemotif.com/catalog/details/82115
Recombinant UBA7 (UBE1L) protein	82077	20 µg	https://www.activemotif.com/catalog/details/82077
Recombinant UBE2W protein	82091	100 µg	https://www.activemotif.com/catalog/details/82091
Recombinant YAP1 protein	82068	20 µg	https://www.activemotif.com/catalog/details/82068
Recombinant YWHAB protein	82045	100 µg	https://www.activemotif.com/catalog/details/82045
Recombinant YWHA E protein	82046	100 µg	https://www.activemotif.com/catalog/details/82046
Recombinant YWHAG protein	82047	100 µg	https://www.activemotif.com/catalog/details/82047
Recombinant YWHAH protein	82048	100 µg	https://www.activemotif.com/catalog/details/82048
Recombinant YWHAQ protein	82049	100 µg	https://www.activemotif.com/catalog/details/82049
Recombinant YWHAZ protein	82050	100 µg	https://www.activemotif.com/catalog/details/82050