

Peptide array: Histone H3(acetyl K14) (ab52946)

Area under curve (AUC) data for all peptides on the pad. Antigen-containing peptides are highlighted in green.

Cell Ref	AUC	Peptide
A1	7.743763542	ab10112 - H4 K91ac
A2	4.990241667	ab10121 - H2B K85ac
A3	4.6146	ab10127 - H2A K95me2
A4	5.780186458	ab10138 - H1.4 T17p
A5	7.69729375	ab10139 - H1.4 T146p
A6	8.563495833	ab113997 - H2A.X Y142p
A7	6.846858333	ab11477 - H3 S10p
A8	6.389183333	ab12946 - H2B K5me3
A9	6.322772917	ab12949 - H2B K23me2
A10	6.01686875	ab12952 - H2A K127me1
A11	5.701210417	ab13211 - H2B K5me1
A12	360.1935531	ab1340 - H3 K4me1
A13	4.189308333	ab13570 - H2B K43me1
A14	8.848139583	ab13835 - HA
A15	6.725952083	ab13837 - Myc
A16	7.208692708	ab140408 - ScH2B T129p
A17	6.309985417	ab14103 - H3 T6p
A18	7.398358333	ab14724 - H4 S1p
A19	7.579836458	ab14799 - H3 T32p
A20	6.482447917	ab14943 - His
A21	4.879867708	ab14952 - H2A R77me2s
A22	7.462257292	ab14964 - H4 K20me2
A23	6.428129167	ab14997 - ScH2A K21ac
A24	6.758172917	ab154782 - H3 R26me2s
A25	698.533426	ab15591 - H3 K14ac
B1	8.845992708	ab15644 - H3 K9me3 + S10p
B2	4.23189375	ab15645 - H2A.X S139p
B3	24.82171771	ab15824 - H4 K8ac
B4	6.3583	ab15829 - V5
B5	3.513202083	ab16635 - H3 K9ac
B6	4.968539583	ab166685 - H2B S112GlcNAc
B7	2.325280208	ab166688 - H2B S36GlcNAc
B8	5.599296875	ab16935 - H3 R17me2a
B9	6.515152083	ab17043 - H4 K20me1
B10	7.1308375	ab17162 - H3 P16hyd + P30hyd
B11	7.531175	ab17567 - H4 K20me3
B12	3.753532292	ab17587 - H1.4 K25me3
B13	7.097652083	ab175905 - H2B K11ac
B14	5.108103125	ab176211 - H2B K23ac
B15	6.829225	ab17632 - H4 K5bio
B16	6.573328125	ab1771 - H3 K9me1
B17	6.316905208	ab1773 - H3 K9me3
B18	7.167990625	ab17770 - H4 R3me1
B19	4.959027083	ab1781 - H3 K27me2
B20	8.0753375	ab1783 - H3 K36me1
B21	4.55063125	ab1784 - H3 K36me2
B22	6.133953125	ab1785 - H3 K36me3
B23	7.978701042	ab178667 - ScH2A unmod(1-30)
B24	6.350145833	ab178668 - ScH2A unmod(102-132)
B25	8.864592708	ab178670 - H2A.Z unmod(1-30)
C1	6.499021875	ab178833 - Flag
C2	5.5456	ab178835 - H2B R99mutO
C3	6.783492708	ab178836 - H4 R23mutO
C4	6.728545833	ab178837 - H4 G48mutD
C5	7.335214583	ab178987 - H1.4 unmod(1-30)
C6	3.0970125	ab178988 - H1.4 S1p
C7	6.919339583	ab178989 - H1.4 T3p
C8	2.846777083	ab178990 - H1.4 S1p + T3p
C9	3.2592375	ab178991 - H1.4 K25me1
C10	6.065013542	ab178992 - H1.4 K25me3 + S26p
C11	5.149640625	ab178993 - H1.4 unmod(130-160)
C12	9.6427125	ab178994 - H2A unmod(1-30)
C13	6.927201042	ab178995 - H2A S1p
C14	8.185180208	ab178996 - H2A S1p + K5ac
C15	5.480855208	ab178997 - H2A S1p + R3me2a + K5ac
C16	5.349136458	ab178998 - H2A S1p + R3me2s + K5ac
C17	6.63090625	ab178999 - H2A S1p + R3cit + K5ac
C18	8.436915625	ab179000 - H2A S1ac + S1p + K5ac + K9ac + K13ac + K15ac
C19	6.803928125	ab179001 - H2A S1p + K5ac + K9ac + K13ac + K15ac
C20	5.8021125	ab179002 - H2A R3cit

C21	7.3753	ab179003 - H2A R3cit + K5ac
C22	6.306530208	ab179004 - H2A R3me2a + K5ac
C23	6.16694375	ab179005 - H2A R3me2s + K5ac
C24	7.52635625	ab179006 - H2A S1ac + K5ac + K9ac + K13ac + K15ac
C25	6.103007292	ab179007 - H2A K5ac + K9ac + K13ac + K15ac
D1	5.58989375	ab179008 - H2A K9me2
D2	4.106441667	ab179009 - H2A K9me3
D3	6.19504375	ab179010 - H2A R11me1
D4	3.260871875	ab179011 - H2A R11me2s
D5	6.542054167	ab179012 - H2A R11me2a
D6	3.474975	ab179013 - H2A K13ac
D7	7.545619792	ab179014 - H2A K15ac
D8	2.750285417	ab179015 - H2A R17me1
D9	4.675188542	ab179016 - H2A R17me2s
D10	7.525526042	ab179017 - H2A R17me2a
D11	12.28766042	ab179018 - H2A unmod(20-50)
D12	6.878404167	ab179019 - H2A P26hyd
D13	15.41500208	ab179020 - H2A R29me1
D14	5.986944792	ab179021 - H2A R29me2s
D15	5.222060417	ab179022 - H2A R29me2a
D16	6.970026042	ab179023 - H2A K36ac
D17	9.44415625	ab179024 - H2A K36cr
D18	6.104194792	ab179025 - H2A Y39hyd
D19	5.609553125	ab179026 - H2A R42me1
D20	4.76350625	ab179027 - H2A R42me2s
D21	6.613227083	ab179028 - H2A R42me2a
D22	5.189961458	ab179029 - H2A unmod(71-103)
D23	6.179613542	ab179030 - H2A R77me1
D24	8.769578125	ab179031 - H2A R77me2a
D25	6.462244792	ab179032 - H2A R88me1
E1	7.797084375	ab179033 - H2A R88me2s
E2	6.7044	ab179034 - H2A R88me2a
E3	5.501382292	ab179035 - H2A K95me1
E4	6.597532292	ab179036 - H2A K95me3
E5	35.69598854	ab179037 - H2A unmod(92-123)
E6	4.112965625	ab179038 - H2A T101GlcNAc
E7	4.710296875	ab179039 - H2A Q104me1
E8	6.4266	ab179040 - H2A unmod(112-129)
E9	3.231097917	ab179041 - H2A K118me1
E10	6.151711458	ab179042 - H2A K118me2
E11	3.845205208	ab179043 - H2A K118me3
E12	7.523727083	ab179044 - H2A K118fo
E13	4.388907292	ab179045 - H2A K118cr
E14	6.2688125	ab179046 - H2A K119cr
E15	6.526213542	ab179047 - H2A K118ubiq
E16	5.689870833	ab179048 - H2A K119ubiq
E17	5.210302083	ab179049 - H2A T120p
E18	7.629775	ab179050 - H2A K125me1
E19	3.993861458	ab179051 - H2A K125me2
E20	5.936077083	ab179052 - H2A K125cr
E21	20.81092083	ab179053 - H2A.X unmod(1-30)
E22	7.644496875	ab179054 - H2A.X K5ac
E23	8.05985	ab179055 - H2A.X unmod(114-142)
E24	4.340346875	ab179056 - H2A.X K118ubiq
E25	7.291536458	ab179057 - H2A.X K119ubiq
F1	4.65959375	ab179058 - H2A.X T120p
F2	6.406410417	ab179059 - H2A.X K127ubiq
F3	6.491573958	ab179060 - H2A.X S139p + Y142p
F4	6.828633333	ab179061 - H2A.Z K4ac
F5	7.249373958	ab179062 - H2A.Z K7ac
F6	5.148471875	ab179063 - H2A.Z K11ac
F7	7.939677083	ab179064 - H2A.Z K4ac + K7ac + K11ac
F8	4.150967708	ab179065 - H2B unmod(1-30)
F9	5.76335	ab179066 - H2B K5ac
F10	5.588785417	ab179067 - H2B K5cr
F11	5.005975	ab179068 - H2B K5fo
F12	4.489783333	ab179069 - H2B P10hyd
F13	4.347561458	ab179070 - H2B K11me1
F14	5.791745833	ab179071 - H2B K11me2
F15	4.775257292	ab179072 - H2B K11me3
F16	5.378907292	ab179073 - H2B K11cr
F17	7.38670625	ab179074 - H2B K12ac
F18	6.199004167	ab179075 - H2B K12cr
F19	5.190120833	ab179076 - H2B K12me1

F20	4.402935417	ab179077 - H2B K12me2
F21	6.648232292	ab179078 - H2B K12me3
F22	6.961238542	ab179079 - H2B K12ac + K15ac
F23	13.60656667	ab179080 - H2B S14p
F24	5.44624375	ab179081 - H2B K15ac
F25	7.189465625	ab179082 - H2B K15cr
G1	7.422947917	ab179083 - H2B K15me1
G2	3.615960417	ab179084 - H2B K15me2
G3	6.069071875	ab179085 - H2B K15me3
G4	5.140815625	ab179086 - H2B K16ac
G5	5.753258333	ab179087 - H2B K16cr
G6	7.955497917	ab179088 - H2B K20ac
G7	5.285376042	ab179089 - H2B K20cr
G8	5.827298958	ab179090 - H2B K20me1
G9	4.08235625	ab179091 - H2B K20me2
G10	6.862752083	ab179092 - H2B K20me3
G11	4.780314583	ab179093 - H2B K23cr
G12	7.513966667	ab179094 - H2B K23me1
G13	6.755660417	ab179095 - H2B K23me3
G14	4.080357292	ab179096 - H2B K24ac
G15	7.239741667	ab179097 - H2B unmod(26-56)
G16	8.548920833	ab179098 - H2B K34cr
G17	6.124084375	ab179099 - H2B K34fo
G18	6.501751042	ab179100 - H2B Y37hyd
G19	5.390021875	ab179101 - H2B Y37p
G20	5.096840625	ab179102 - H2B K43me2
G21	7.075672917	ab179103 - H2B K43me3
G22	4.967373958	ab179104 - H2B K46fo
G23	4.564527083	ab179105 - H2B K46ac
G24	6.38430625	ab179106 - H2B K46me1
G25	7.484759375	ab179107 - H2B K46me2
H1	5.287025	ab179108 - H2B K46me3
H2	4.796059375	ab179109 - H2B unmod(51-81)
H3	6.751653125	ab179110 - H2B K57me1
H4	6.208516667	ab179111 - H2B K57me2
H5	6.609208333	ab179112 - H2B K57me3
H6	7.346307292	ab179113 - H2B unmod(73-103)
H7	5.163580208	ab179114 - H2B R79me1
H8	5.916489583	ab179115 - H2B R79me2s
H9	4.085957292	ab179116 - H2B R79me2a
H10	6.301396875	ab179117 - H2B Y83hyd
H11	5.69556875	ab179118 - H2B K85me1
H12	6.152984375	ab179119 - H2B K85me2
H13	8.359847917	ab179120 - H2B K85me3
H14	4.594782292	ab179121 - H2B unmod(94-125)
H15	7.702971875	ab179122 - H2B R99me1
H16	7.523261458	ab179123 - H2B R99me2s
H17	6.102269792	ab179124 - H2B R99me2a
H18	6.526528125	ab179125 - H2B K108fo
H19	5.189123958	ab179126 - H2B K116fo
H20	4.260232292	ab179127 - H2B K116me1
H21	4.435948958	ab179128 - H2B K116me2
H22	3.159158333	ab179129 - H2B K116me3
H23	7.205473958	ab179130 - H2B K116ubiq
H24	4.506109375	ab179131 - H2B K120ac
H25	5.112513542	ab179132 - H2B K120fo
I1	4.26215625	ab179133 - H2B K120ubiq
I2	7.584079167	ab179134 - H2B K125ubiq
I3	4.793533333	ab179135 - Sch2B unmod(102-132)
I4	5.903226042	ab179136 - H3 unmod(1-33)
I5	4.868833333	ab179137 - H3 R2cit
I6	6.679069792	ab179138 - H3 R2cit + K4me2
I7	3.612551042	ab179139 - H3 R2cit + K4me3
I8	236.0596635	ab179140 - H3 R2cit + K4me3 + K9ac + K14ac + K18ac
I9	38.25292917	ab179141 - H3 R2cit + K4ac + K9ac + K14ac + K18ac
I10	11.23532083	ab179142 - H3 R2me1
I11	6.413879167	ab179143 - H3 R2me1 + K4me2
I12	7.508363542	ab179144 - H3 R2me1 + K4me3
I13	25.86424479	ab179145 - H3 R2me1 + K4me3 + K9ac + K14ac + K18ac
I14	5.889690625	ab179146 - H3 R2me2s
I15	6.386034375	ab179147 - H3 R2me2s + K4me2
I16	56.39580104	ab179148 - H3 R2me2s + K4me3
I17	19.24466771	ab179149 - H3 R2me2s + K4me3 + K9ac + K14ac + K18ac
I18	9.023678125	ab179150 - H3 R2me2a

I19	7.21933125	ab179151 - H3 R2me2a + K4me2
I20	9.888982292	ab179152 - H3 R2me2a + K4me3
I21	6.604139583	ab179153 - H3 R2me2a + K4me3 + S10p
I22	6.47856875	ab179154 - H3 R2me2a + T3p + K4me3
I23	19.71184583	ab179155 - H3 R2me2a + K4ac + K9ac + K14ac + K18ac
I24	16.71485521	ab179156 - H3 R2me2a + K4me3 + K9ac + K14ac + K18ac
I25	31.38820208	ab179157 - H3 R2me2a + T3p + K4me3 + K9ac + K14ac + K18ac
J1	1089.318415	ab179158 - H3 R2me2a + K4me3 + K9ac + S10p + K14ac + K18ac
J2	5.511961458	ab179159 - H3 T3p
J3	17.16456875	ab179160 - H3 T3p + K4me1
J4	4.399109375	ab179161 - H3 T3p + K4me2
J5	37.28294167	ab179162 - H3 T3p + K4me3 + K9ac + K14ac + K18ac
J6	6.574690625	ab179163 - H3 K4ac
J7	5.664348958	ab179164 - H3 K4ac + K9ac
J8	895.9098615	ab179165 - H3 K4ac + K14ac
J9	8.532223958	ab179166 - H3 K4ac + K18ac
J10	10.34022813	ab179167 - H3 K4ac + K9me3
J11	831.2563771	ab179168 - H3 K4ac + K9ac + K14ac
J12	5.369501042	ab179169 - H3 K4ac + K9ac + K18ac
J13	100.9547635	ab179170 - H3 K4ac + K14ac + K18ac
J14	52.59948646	ab179171 - H3 K4ac + K9ac + K14ac + K18ac
J15	21.23186979	ab179172 - H3 K4ac + K9me1 + K14ac + K18ac
J16	66.78364271	ab179173 - H3 K4ac + K9me2 + K14ac + K18ac
J17	20.60543021	ab179174 - H3 K4ac + K9me3 + K14ac + K18ac
J18	5431.327783	ab179175 - H3 K4ac + K9ac + S10p + K14ac + K18ac
J19	7.611339583	ab179176 - H3 K4cr
J20	5.663682292	ab179177 - H3 K4me1 + K9me2
J21	7.724995833	ab179178 - H3 K4me1 + K9ac + K18ac
J22	60.38076771	ab179179 - H3 K4me1 + K9ac + K14ac + K18ac
J23	6.805134375	ab179180 - H3 K4me2 + T6p
J24	6.321903125	ab179181 - H3 K4me2 + K9me2
J25	5.534060417	ab179182 - H3 K4me2 + K9ac + K18ac
K1	47.24845417	ab179183 - H3 K4me2 + K9ac + K14ac + K18ac
K2	6.211821875	ab179184 - H3 K4me3 + T6p
K3	8.374830208	ab179185 - H3 K4me3 + K9ac
K4	6.23066875	ab179186 - H3 K4me3 + K9me2
K5	5.224580208	ab179187 - H3 K4me3 + K9me3
K6	8.008953125	ab179188 - H3 K4me3 + S10p
K7	1458.867905	ab179189 - H3 K4me3 + K14ac
K8	5.755329167	ab179190 - H3 K4me3 + K18ac
K9	13.74012292	ab179191 - H3 K4me3 + R8me2s + K9me3
K10	8.495579167	ab179192 - H3 K4me3 + R8me2a + K9me3
K11	5.213619792	ab179193 - H3 K4me3 + K9ac + S10p
K12	920.8027167	ab179194 - H3 K4me3 + K9ac + K14ac
K13	4.98001875	ab179195 - H3 K4me3 + K9ac + K18ac
K14	668.4255083	ab179196 - H3 K4me3 + K14ac + K18ac
K15	36.61162604	ab179197 - H3 K4me3 + K9ac + K14ac + K18ac
K16	3881.592292	ab179198 - H3 K4me3 + K9ac + S10p + K14ac + K18ac
K17	26.77483229	ab179199 - H3 K4me3 + T6p + K9ac + K14ac + K18ac
K18	39.91894063	ab179200 - H3 K4me2 + T6p + K9ac + K14ac + K18ac
K19	9.295995833	ab179201 - H3 K4bio + K9bio + K18bio
K20	5.0856125	ab179202 - H3 T6p + K9me3
K21	6.089442708	ab179203 - H3 T6p + R8me2a + K9me3
K22	41.59995625	ab179204 - H3 T6p + K9ac + K14ac + K18ac
K23	5.534572917	ab179205 - H3 R8me1
K24	6.437760417	ab179206 - H3 R8me1 + K9me1
K25	137.6149969	ab179207 - H3 R8me1 + K9me2
L1	4.21708125	ab179208 - H3 R8me1 + K9me3
L2	10.81601354	ab179209 - H3 R8me2s
L3	4.116930208	ab179210 - H3 R8me2s + K9me1
L4	7.881691667	ab179211 - H3 R8me2s + K9me2
L5	3.235985417	ab179212 - H3 R8me2s + K9me3
L6	5.983214583	ab179213 - H3 R8me2a
L7	6.156619792	ab179214 - H3 R8me2a + K9me1
L8	7.103171875	ab179215 - H3 R8me2a + K9me2
L9	4.6491	ab179216 - H3 R8me2a + K9me3
L10	6.947008333	ab179217 - H3 R8cit
L11	2126.850757	ab179218 - H3 K9ac + K14ac
L12	779.5163427	ab179219 - H3 K9ac + K18ac
L13	87.92009271	ab179220 - H3 K9ac + K14ac + K18ac
L14	20.82482604	ab179221 - H3 K9ac + K14ac + K18ac + K23ac + K27ac
L15	8.234063542	ab179222 - H3 K9ac + S10p
L16	10.56236875	ab179223 - H3 K9cr
L17	6.880666667	ab179224 - H3 K9me2

L18	6.757514583	ab179225 - H3 K9me2 + S10p
L19	3.714235417	ab179226 - H3 K9me2 + K27me2
L20	5.472560417	ab179227 - H3 S10p + S28p
L21	3995.382013	ab179228 - H3 S10p K14ac
L22	7.786691667	ab179229 - H3 K14me1
L23	5.749334375	ab179230 - H3 K14me2
L24	5.737888542	ab179231 - H3 K14me3
L25	203.2873573	ab179232 - H3 K14ac + K18ac
M1	7.791397917	ab179233 - H3 R17cit
M2	2.961941667	ab179234 - H3 R17me1
M3	4.420439583	ab179235 - H3 R17me2s
M4	6.456592708	ab179236 - H3 R17me2a + K18ac
M5	4.986939583	ab179237 - H3 K18cr
M6	6.254508333	ab179238 - H3 K18me1
M7	5.036339583	ab179239 - H3 K18me2
M8	94.76198542	ab179240 - H3 K18me3 + K36me3
M9	4.980545833	ab179241 - H3 K23cr
M10	7.195925	ab179242 - H3 K23me1
M11	5.504	ab179243 - H3 K23me2
M12	7.610982292	ab179244 - H3 K23me3
M13	6.335970833	ab179245 - H3 T22p + K23me3
M14	9.414534375	ab179246 - H3 K23pr
M15	9.168563542	ab179247 - H3 unmod(20-51)
M16	6.208252083	ab179248 - H3 R26me1
M17	5.457695833	ab179249 - H3 R26me2a
M18	7.040951042	ab179250 - H3 R26me2a + K27me1
M19	4.18954375	ab179251 - H3 R26me2a + K27me3
M20	7.575408333	ab179252 - H3 K27ac + S28p
M21	3.888333333	ab179253 - H3 K27cr
M22	4.909767708	ab179254 - H3 K27me1
M23	4.427460417	ab179255 - H3 K27me2 + S28p
M24	4.275	ab179256 - H3 K27me3
M25	5.355116667	ab179257 - H3 K27me3 + S28p
N1	4.271148958	ab179258 - H3 P30hpro
N2	6.1539375	ab179259 - H3 P30(5,5me2)
N3	7.830505208	ab179260 - H3 P30
N4	6.441470833	ab179261 - H3 P30mutA
N5	6.33308125	ab179262 - H3 T32GlcNAc
N6	4.918283333	ab179263 - H3 K36ac
N7	5.750034375	ab179264 - H3 P38hpro
N8	4.738259375	ab179265 - H3 P38(5,5me2)
N9	3.548710417	ab179266 - H3 P38
N10	6.173884375	ab179267 - H3 P38mutA
N11	7.052810417	ab179268 - H3 Y41p
N12	9.546390625	ab179269 - H3 unmod(50-80)
N13	2.857921875	ab179270 - H3 K56ac
N14	6.759707292	ab179271 - H3 K56cr
N15	5.430234375	ab179272 - H3 K56fo
N16	9.234478125	ab179273 - H3 K56me1
N17	8.152410417	ab179274 - H3 K56me2
N18	9.690975	ab179275 - H3 K56me3
N19	9.375766667	ab179276 - H3 R63me1
N20	7.264438542	ab179277 - H3 R63me2s
N21	6.7136875	ab179278 - H3 R63me2a
N22	8.334422917	ab179279 - H3 K64ac
N23	6.6142875	ab179280 - H3 K64me2
N24	7.638523958	ab179281 - H3 K64me3
N25	4.948423958	ab179282 - H3 unmod(73-103)
O1	6.220811458	ab179283 - H3 K79ac
O2	6.829944792	ab179284 - H3 K79fo
O3	5.915734375	ab179285 - H3 T80p
O4	8.317807292	ab179286 - H3 K79me3 + T80p
O5	3.290721875	ab179287 - H3 unmod(99-135)
O6	6.654433333	ab179288 - H3 T118O-ac
O7	4.794264583	ab179289 - H3 K122fo
O8	6.301116667	ab179290 - H3 K122me1
O9	6.133410417	ab179291 - H3 R128me1
O10	6.6349	ab179292 - H3 R128me2s
O11	8.6963625	ab179293 - H3.3 unmod(10-40)
O12	6.022191667	ab179294 - H3.3 S28p
O13	6.290358333	ab179295 - H3.3 S28p + S31p
O14	5.821161458	ab179296 - H3.3 S31p
O15	20.98545208	ab179297 - H4 unmod(1-30)
O16	8.083117708	ab179298 - H4 S1p + R3me1

O17	9.705535417	ab179299 - H4 S1p + R3me2s
O18	9.386715625	ab179300 - H4 S1p + R3me2a
O19	36.41172813	ab179301 - H4 S1p + K5ac + K8ac + K12ac + K16ac
O20	8.309907292	ab179302 - H4 R3cit
O21	20.34015	ab179303 - H4 R3me1 + K5ac + K8ac + K12ac + K16ac + K20ac
O22	7.170676042	ab179304 - H4 R3me2s
O23	14.45241875	ab179305 - H4 R3me2s + K5ac + K8ac + K12ac + K16ac + K20ac
O24	6.656564583	ab179306 - H4 R3me2a
O25	8.871502083	ab179307 - H4 R3me2a + K5ac
P1	115.507101	ab179308 - H4 R3me2a + K5ac + K8ac + K12ac + K16ac + K20ac
P2	12.33923542	ab179309 - H4 K5ac
P3	43.82641875	ab179310 - H4 K5ac + K8ac
P4	39.32337083	ab179311 - H4 K5ac + K12ac
P5	36.56816458	ab179312 - H4 K5ac + K16ac
P6	43.38765625	ab179313 - H4 K5ac + K8ac + K12ac
P7	58.28206563	ab179314 - H4 K5ac + K8me1 + K12ac
P8	46.40600938	ab179315 - H4 K5ac + K8ac + K12ac + K16ac
P9	46.3168625	ab179316 - H4 K5ac + K8me1 + K12ac + K16ac
P10	45.210175	ab179317 - H4 K5ac + K8ac + K12me1 + K16ac
P11	7.959383333	ab179318 - H4 K5me1
P12	6.936455208	ab179319 - H4 K5me1 + K8ac + K12me1
P13	8.812310417	ab179320 - H4 K5me1 + K8me1 + K12me1
P14	7.141033333	ab179321 - H4 K5me1 + K8ac + K12ac + K16ac
P15	6.104022917	ab179322 - H4 K5me2
P16	6.218875	ab179323 - H4 K5me3
P17	7.051861458	ab179324 - H4 K5cr
P18	7.816776042	ab179325 - H4 K8ac + K12ac
P19	24.24913229	ab179326 - H4 K8ac + K16ac
P20	6.160719792	ab179327 - H4 K8cr
P21	10.19455833	ab179328 - H4 K8me1
P22	6.9296625	ab179329 - H4 K12ac
P23	10.32596875	ab179330 - H4 K12ac + K16ac
P24	9.696511458	ab179331 - H4 K12ac + K16ac + K20me2
P25	6.941190625	ab179332 - H4 K12ac + K16ac + K20me3
Q1	6.88050625	ab179333 - H4 K5mutQ + K8mutQ + K12ac + K16mutQ + K20mutQ
Q2	6.342891667	ab179334 - H4 K12cr
Q3	6.000338542	ab179335 - H4 K12fo
Q4	6.762253125	ab179336 - H4 K12me1
Q5	7.852092708	ab179337 - H4 K12me2
Q6	6.644042708	ab179338 - H4 K12me3
Q7	7.811623958	ab179339 - H4 K16ac
Q8	12.45385625	ab179340 - H4 K16ac + K20ac
Q9	6.834254167	ab179341 - H4 K16cr
Q10	23.91655208	ab179342 - H4 K16me1
Q11	29.68835417	ab179343 - H4 K16me3
Q12	7.579177083	ab179344 - H4 R17me1
Q13	10.75233438	ab179345 - H4 R17me2s
Q14	9.476148958	ab179346 - H4 R17me2a
Q15	8.844923958	ab179347 - H4 R19me1
Q16	6.935330208	ab179348 - H4 R19me2s
Q17	7.313461458	ab179349 - H4 R19me2a
Q18	8.232247917	ab179350 - H4 K20ac
Q19	8.22119375	ab179351 - H4 R23me1
Q20	7.322741667	ab179352 - H4 R23me2s
Q21	6.676034375	ab179353 - H4 R23me2a
Q22	7.250242708	ab179354 - H4 unmod(24-54)
Q23	5.340871875	ab179355 - H4 K31fo
Q24	7.923048958	ab179356 - H4 K31me1
Q25	5.788823958	ab179357 - H4 K31me2
R1	5.867261458	ab179358 - H4 K31me3
R2	10.23785729	ab179359 - H4 P32hyd
R3	6.617745833	ab179360 - H4 R35me1
R4	8.054755208	ab179361 - H4 R35me2s
R5	4.888417708	ab179362 - H4 R35me2a
R6	7.511554167	ab179363 - H4 S47O-ac
R7	5.935198958	ab179364 - H4 S47GlcNAc
R8	8.721244792	ab179365 - H4 S47p
R9	7.671072917	ab179366 - H4 unmod(45-75)
R10	7.916197917	ab179367 - H4 Y51hyd
R11	6.423180208	ab179368 - H4 R55me1
R12	5.40114375	ab179369 - H4 R55me2s
R13	8.146101042	ab179370 - H4 R55me2a
R14	6.220884375	ab179371 - H4 K59fo
R15	7.237086458	ab179372 - H4 K59me1

R16	5.290448958	ab179373 - H4 K59me2
R17	6.583902083	ab179374 - H4 R67me1
R18	6.51024375	ab179375 - H4 R67me2s
R19	6.017290625	ab179376 - H4 R67me2a
R20	9.538041667	ab179377 - H4 unmod(71-102)
R21	6.315671875	ab179378 - H4 K77me1
R22	7.529320833	ab179379 - H4 K77me2
R23	7.409002083	ab179380 - H4 K77me3
R24	7.858634375	ab179381 - H4 K79fo
R25	5.135179167	ab179382 - H4 K79me1
S1	6.99115	ab179383 - H4 K79me3
S2	6.559146875	ab179384 - H4 T82O-ac
S3	6.936936458	ab179385 - H4 Y88hyd
S4	8.372185417	ab179386 - H4 Y88nitro
S5	6.948528125	ab179387 - H4 K91fo
S6	6.610402083	ab179403 - H2A K5ac
S7	3.668679167	ab179404 - H2A K9me1
S8	5.88945	ab18504 - H2B S32p
S9	4.508220833	ab19828 - ScH2A S129p
S10	7.874505208	ab20631 - H3 R26cit
S11	5.037026042	ab20635 - H3 T3p + K4me3
S12	3.144766667	ab21043 - H3 K37me1
S13	5.29621875	ab21998 - H1.4 K25me2
S14	6.816302083	ab22079 - H4 K59me3
S15	8.619019792	ab22080 - H3 K122me2
S16	8.155416667	ab22082 - H3 K122me3
S17	9.530441667	ab22214 - H3 K37me2
S18	6.881955208	ab22398 - H2A R3me2a
S19	328.2839177	ab22399 - H2A R3me2s
S20	5.53118125	ab23539 - H3 T22p
S21	4.064209375	ab24003 - H3 K18ac
S22	5.990451042	ab24404 - H3 K27ac
S23	3.436411458	ab24417 - H3 K37me3
S24	5.760919792	ab24444 - H3 T11p
S25	4.7964	ab24659 - H2A K127me3
T1	8.459740625	ab24660 - H2A K125me3
T2	5.354745833	ab26307 - H3 T45p
T3	4.504613542	ab26342 - H2B K5me2
T4	6.902394792	ab26739 - H4 T30p
T5	7.45275625	ab27896 - H4 T96p
T6	25.88189792	ab27898 - H4 T80p
T7	6.944921875	ab28756 - H2A R3me1
T8	6.829336458	ab32876 - H3 R2cit + R8cit + R17cit
T9	7.201973958	ab33047 - H2A K127me2
T10	4.455041667	ab33505 - H3 T118p
T11	7.5769375	ab33594 - H3 R128me2a
T12	8.911010417	ab34466 - H3 K122ac
T13	5.600446875	ab36927 - H2A S122p
T14	7.490321875	ab4555 - H3 K79me1
T15	8.074707292	ab4556 - H3 K79me2
T16	5.993363542	ab4557 - H3 K79me3
T17	7.340716667	ab4560 - H4 K79me2
T18	4.471402083	ab46854 - H3 K18me3
T19	11.06325208	ab48359 - H3 K23ac
T20	7.192303125	ab53244 - H3 K64me1
T21	7.616183333	ab54016 - H2A K9ac
T22	5.290917708	ab5499 - H3 S28p
T23	6.219342708	ab7768 - H3 K4me2
T24	7.697273958	ab89018 - H1.4 S26p
T25	6.7290875	ab92374 - H3 K4me3 (HEX1)
U1	8.212561458	ab95444 - H4 K16me2