

Peptide array: Histone H3(acetyl K4+K9 + K14 + K18 + K23 + K27) (ab300641)

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

Cell Ref	AUC	Peptide
A1	11.18710313	ab10112 - H4 K91ac
A2	8.993729167	ab10121 - H2B K85ac
A3	14.97223854	ab10127 - H2A K95me2
A4	10.78184375	ab10138 - H1.4 T17p
A5	12.97386563	ab10139 - H1.4 T146p
A6	13.16373229	ab113997 - H2A.X Y142p
A7	12.09142083	ab11477 - H3 S10p
A8	15.73031458	ab12946 - H2B K5me3
A9	13.04058854	ab12949 - H2B K23me2
A10	9.6636625	ab12952 - H2A K127me1
A11	13.31834479	ab13211 - H2B K5me1
A12	156.100075	ab1340 - H3 K4me1
A13	15.51400104	ab13570 - H2B K43me1
A14	10.22958854	ab13835 - HA
A15	11.88071042	ab13837 - Myc
A16	10.62824167	ab140408 - ScH2B T129p
A17	399.1489969	ab14103 - H3 T6p
A18	15.20344792	ab14724 - H4 S1p
A19	28.15424375	ab14799 - H3 T32p
A20	10.88517188	ab14943 - His
A21	10.76787396	ab14952 - H2A R77me2s
A22	11.02625521	ab14964 - H4 K20me2
A23	10.62834896	ab14997 - ScH2A K21ac
A24	22.34912396	ab154782 - H3 R26me2s
A25	242.1068625	ab15591 - H3 K14ac
B1	12.89348438	ab15644 - H3 K9me3 + S10p
B2	10.8415	ab15645 - H2A.X S139p
B3	28.42298958	ab15824 - H4 K8ac
B4	11.6489125	ab15829 - V5
B5	1040.467835	ab16635 - H3 K9ac
B6	10.82960208	ab166685 - H2B S112GlcNAc
B7	10.87623229	ab166688 - H2B S36GlcNAc
B8	13.82718021	ab16935 - H3 R17me2a
B9	12.19808333	ab17043 - H4 K20me1
B10	46.29039583	ab17162 - H3 P16hyd + P30hyd
B11	10.30188229	ab17567 - H4 K20me3
B12	10.74049583	ab17587 - H1.4 K25me3
B13	12.24903854	ab175905 - H2B K11ac
B14	13.59112188	ab176211 - H2B K23ac
B15	10.44190208	ab17632 - H4 K5bio
B16	17.07587917	ab1771 - H3 K9me1
B17	19.25034688	ab1773 - H3 K9me3
B18	11.81754792	ab17770 - H4 R3me1
B19	17.46767396	ab1781 - H3 K27me2
B20	47.21062917	ab1783 - H3 K36me1
B21	16.7482125	ab1784 - H3 K36me2
B22	15.71014479	ab1785 - H3 K36me3
B23	13.75028021	ab178667 - ScH2A unmod(1-30)
B24	12.8146125	ab178668 - ScH2A unmod(102-132)
B25	15.38465208	ab178670 - H2A.Z unmod(1-30)
C1	9.281926042	ab178833 - Flag
C2	17.82327708	ab178835 - H2B R99mutO
C3	11.13282188	ab178836 - H4 R23mutO
C4	12.98500521	ab178837 - H4 G48mutD
C5	76.03140208	ab178987 - H1.4 unmod(1-30)
C6	10.88413229	ab178988 - H1.4 S1p
C7	10.80067813	ab178989 - H1.4 T3p
C8	9.564069792	ab178990 - H1.4 S1p + T3p
C9	10.34132917	ab178991 - H1.4 K25me1
C10	9.35703125	ab178992 - H1.4 K25me3 + S26p
C11	16.478275	ab178993 - H1.4 unmod(130-160)
C12	12.87463646	ab178994 - H2A unmod(1-30)
C13	11.83657604	ab178995 - H2A S1p
C14	14.15744583	ab178996 - H2A S1p + K5ac
C15	12.423675	ab178997 - H2A S1p + R3me2a + K5ac
C16	14.84399167	ab178998 - H2A S1p + R3me2s + K5ac
C17	13.94973438	ab178999 - H2A S1p + R3cit + K5ac
C18	13.47310625	ab179000 - H2A S1ac + S1p + K5ac + K9ac + K13ac + K15ac
C19	13.48055104	ab179001 - H2A S1p + K5ac + K9ac + K13ac + K15ac
C20	10.05383125	ab179002 - H2A R3cit

C21	14.09923542	ab179003 - H2A R3cit + K5ac
C22	11.6207125	ab179004 - H2A R3me2a + K5ac
C23	15.78830833	ab179005 - H2A R3me2s + K5ac
C24	13.33140521	ab179006 - H2A S1ac + K5ac + K9ac + K13ac + K15ac
C25	10.74708438	ab179007 - H2A K5ac + K9ac + K13ac + K15ac
D1	11.47299792	ab179008 - H2A K9me2
D2	14.29038854	ab179009 - H2A K9me3
D3	11.09223021	ab179010 - H2A R11me1
D4	12.27807188	ab179011 - H2A R11me2s
D5	12.00346979	ab179012 - H2A R11me2a
D6	10.73464688	ab179013 - H2A K13ac
D7	12.00633646	ab179014 - H2A K15ac
D8	8.346301042	ab179015 - H2A R17me1
D9	11.16615417	ab179016 - H2A R17me2s
D10	11.75742292	ab179017 - H2A R17me2a
D11	13.40892604	ab179018 - H2A unmod(20-50)
D12	14.0002625	ab179019 - H2A P26hyd
D13	12.48378542	ab179020 - H2A R29me1
D14	13.2016625	ab179021 - H2A R29me2s
D15	12.26999896	ab179022 - H2A R29me2a
D16	172.4686594	ab179023 - H2A K36ac
D17	14.56407917	ab179024 - H2A K36cr
D18	11.46519688	ab179025 - H2A Y39hyd
D19	12.05008958	ab179026 - H2A R42me1
D20	11.75397396	ab179027 - H2A R42me2s
D21	12.53454792	ab179028 - H2A R42me2a
D22	11.912	ab179029 - H2A unmod(71-103)
D23	14.85869583	ab179030 - H2A R77me1
D24	13.06458438	ab179031 - H2A R77me2a
D25	11.34759792	ab179032 - H2A R88me1
E1	13.98682708	ab179033 - H2A R88me2s
E2	10.13706458	ab179034 - H2A R88me2a
E3	16.72746979	ab179035 - H2A K95me1
E4	15.86474375	ab179036 - H2A K95me3
E5	8.712966667	ab179037 - H2A unmod(92-123)
E6	12.05588958	ab179038 - H2A T101GlcNAc
E7	13.17268542	ab179039 - H2A Q104me1
E8	10.84533438	ab179040 - H2A unmod(112-129)
E9	9.895383333	ab179041 - H2A K118me1
E10	11.06931875	ab179042 - H2A K118me2
E11	10.51893229	ab179043 - H2A K118me3
E12	12.42504479	ab179044 - H2A K118fo
E13	11.30603021	ab179045 - H2A K118cr
E14	10.96610208	ab179046 - H2A K119cr
E15	10.82106667	ab179047 - H2A K118ubiq
E16	10.2883875	ab179048 - H2A K119ubiq
E17	12.85738958	ab179049 - H2A T120p
E18	12.36996667	ab179050 - H2A K125me1
E19	10.75279583	ab179051 - H2A K125me2
E20	13.03120938	ab179052 - H2A K125cr
E21	14.42345625	ab179053 - H2A.X unmod(1-30)
E22	18.60149063	ab179054 - H2A.X K5ac
E23	13.1274375	ab179055 - H2A.X unmod(114-142)
E24	11.818775	ab179056 - H2A.X K118ubiq
E25	13.12014896	ab179057 - H2A.X K119ubiq
F1	10.8435625	ab179058 - H2A.X T120p
F2	15.09074688	ab179059 - H2A.X K127ubiq
F3	13.29520729	ab179060 - H2A.X S139p + Y142p
F4	19.78765938	ab179061 - H2A.Z K4ac
F5	16.91978229	ab179062 - H2A.Z K7ac
F6	14.26270833	ab179063 - H2A.Z K11ac
F7	36.94903542	ab179064 - H2A.Z K4ac + K7ac + K11ac
F8	13.42890313	ab179065 - H2B unmod(1-30)
F9	35.62222813	ab179066 - H2B K5ac
F10	16.33150833	ab179067 - H2B K5cr
F11	14.77875313	ab179068 - H2B K5fo
F12	12.69065521	ab179069 - H2B P10hyd
F13	16.78931771	ab179070 - H2B K11me1
F14	12.14295417	ab179071 - H2B K11me2
F15	11.34690729	ab179072 - H2B K11me3
F16	12.58793333	ab179073 - H2B K11cr
F17	15.50881563	ab179074 - H2B K12ac
F18	13.01205938	ab179075 - H2B K12cr
F19	12.36452813	ab179076 - H2B K12me1

F20	10.98345104	ab179077 - H2B K12me2
F21	8.233360417	ab179078 - H2B K12me3
F22	106.1154823	ab179079 - H2B K12ac + K15ac
F23	9.282607292	ab179080 - H2B S14p
F24	16.9603875	ab179081 - H2B K15ac
F25	13.45611458	ab179082 - H2B K15cr
G1	16.73335625	ab179083 - H2B K15me1
G2	12.17547188	ab179084 - H2B K15me2
G3	17.08803646	ab179085 - H2B K15me3
G4	11.28061458	ab179086 - H2B K16ac
G5	10.00826146	ab179087 - H2B K16cr
G6	10.68018229	ab179088 - H2B K20ac
G7	11.48883854	ab179089 - H2B K20cr
G8	13.86694896	ab179090 - H2B K20me1
G9	9.038844792	ab179091 - H2B K20me2
G10	13.41062708	ab179092 - H2B K20me3
G11	14.84604792	ab179093 - H2B K23cr
G12	12.43895104	ab179094 - H2B K23me1
G13	12.48104583	ab179095 - H2B K23me3
G14	12.02188125	ab179096 - H2B K24ac
G15	17.00719271	ab179097 - H2B unmod(26-56)
G16	12.018575	ab179098 - H2B K34cr
G17	13.73743958	ab179099 - H2B K34fo
G18	14.77135104	ab179100 - H2B Y37hyd
G19	12.06646354	ab179101 - H2B Y37p
G20	10.35658438	ab179102 - H2B K43me2
G21	14.44126771	ab179103 - H2B K43me3
G22	10.25552396	ab179104 - H2B K46fo
G23	11.81611771	ab179105 - H2B K46ac
G24	14.37674271	ab179106 - H2B K46me1
G25	10.36891979	ab179107 - H2B K46me2
H1	13.24309896	ab179108 - H2B K46me3
H2	16.78726979	ab179109 - H2B unmod(51-81)
H3	16.45062813	ab179110 - H2B K57me1
H4	8.461516667	ab179111 - H2B K57me2
H5	16.01986563	ab179112 - H2B K57me3
H6	14.09730104	ab179113 - H2B unmod(73-103)
H7	12.63896042	ab179114 - H2B R79me1
H8	14.14486771	ab179115 - H2B R79me2s
H9	10.41549167	ab179116 - H2B R79me2a
H10	16.49645521	ab179117 - H2B Y83hyd
H11	13.06502396	ab179118 - H2B K85me1
H12	10.58656875	ab179119 - H2B K85me2
H13	15.99754271	ab179120 - H2B K85me3
H14	16.90067292	ab179121 - H2B unmod(94-125)
H15	15.42900833	ab179122 - H2B R99me1
H16	12.79955208	ab179123 - H2B R99me2s
H17	14.80774271	ab179124 - H2B R99me2a
H18	14.20433438	ab179125 - H2B K108fo
H19	11.54826979	ab179126 - H2B K116fo
H20	10.19769583	ab179127 - H2B K116me1
H21	11.31553542	ab179128 - H2B K116me2
H22	11.82798438	ab179129 - H2B K116me3
H23	15.28690417	ab179130 - H2B K116ubiq
H24	10.90050833	ab179131 - H2B K120ac
H25	13.09406458	ab179132 - H2B K120fo
I1	12.48190625	ab179133 - H2B K120ubiq
I2	12.12757292	ab179134 - H2B K125ubiq
I3	17.94421771	ab179135 - Sch2B unmod(102-132)
I4	311.9567229	ab179136 - H3 unmod(1-33)
I5	10.96178125	ab179137 - H3 R2cit
I6	28.39462292	ab179138 - H3 R2cit + K4me2
I7	10.10101354	ab179139 - H3 R2cit + K4me3
I8	1691.461384	ab179140 - H3 R2cit + K4me3 + K9ac + K14ac + K18ac
I9	2549.198054	ab179141 - H3 R2cit + K4ac + K9ac + K14ac + K18ac
I10	358.7207552	ab179142 - H3 R2me1
I11	19.89036458	ab179143 - H3 R2me1 + K4me2
I12	43.52552604	ab179144 - H3 R2me1 + K4me3
I13	1364.924968	ab179145 - H3 R2me1 + K4me3 + K9ac + K14ac + K18ac
I14	20.19716875	ab179146 - H3 R2me2s
I15	183.7168104	ab179147 - H3 R2me2s + K4me2
I16	149.3892813	ab179148 - H3 R2me2s + K4me3
I17	2244.849724	ab179149 - H3 R2me2s + K4me3 + K9ac + K14ac + K18ac
I18	18.28171979	ab179150 - H3 R2me2a

I19	468.691974	ab179151 - H3 R2me2a + K4me2
I20	16.1572125	ab179152 - H3 R2me2a + K4me3
I21	82.46433854	ab179153 - H3 R2me2a + K4me3 + S10p
I22	13.98962083	ab179154 - H3 R2me2a + T3p + K4me3
I23	2470.008223	ab179155 - H3 R2me2a + K4ac + K9ac + K14ac + K18ac
I24	2132.409966	ab179156 - H3 R2me2a + K4me3 + K9ac + K14ac + K18ac
I25	2195.242056	ab179157 - H3 R2me2a + T3p + K4me3 + K9ac + K14ac + K18ac
J1	1853.014913	ab179158 - H3 R2me2a + K4me3 + K9ac + S10p + K14ac + K18ac
J2	123.1132031	ab179159 - H3 T3p
J3	225.4048146	ab179160 - H3 T3p + K4me1
J4	36.75748854	ab179161 - H3 T3p + K4me2
J5	2485.444419	ab179162 - H3 T3p + K4me3 + K9ac + K14ac + K18ac
J6	25.75958229	ab179163 - H3 K4ac
J7	194.4330104	ab179164 - H3 K4ac + K9ac
J8	1203.465297	ab179165 - H3 K4ac + K14ac
J9	1046.213422	ab179166 - H3 K4ac + K18ac
J10	36.48634896	ab179167 - H3 K4ac + K9me3
J11	1037.784747	ab179168 - H3 K4ac + K9ac + K14ac
J12	2032.703177	ab179169 - H3 K4ac + K9ac + K18ac
J13	1255.827929	ab179170 - H3 K4ac + K14ac + K18ac
J14	2036.039916	ab179171 - H3 K4ac + K9ac + K14ac + K18ac
J15	2070.849323	ab179172 - H3 K4ac + K9me1 + K14ac + K18ac
J16	1804.392722	ab179173 - H3 K4ac + K9me2 + K14ac + K18ac
J17	2303.913847	ab179174 - H3 K4ac + K9me3 + K14ac + K18ac
J18	1987.429463	ab179175 - H3 K4ac + K9ac + S10p + K14ac + K18ac
J19	483.4170344	ab179176 - H3 K4cr
J20	53.37439896	ab179177 - H3 K4me1 + K9me2
J21	1800.869208	ab179178 - H3 K4me1 + K9ac + K18ac
J22	2213.330817	ab179179 - H3 K4me1 + K9ac + K14ac + K18ac
J23	369.0637719	ab179180 - H3 K4me2 + T6p
J24	45.973625	ab179181 - H3 K4me2 + K9me2
J25	2044.360093	ab179182 - H3 K4me2 + K9ac + K18ac
K1	2080.69707	ab179183 - H3 K4me2 + K9ac + K14ac + K18ac
K2	310.9157125	ab179184 - H3 K4me3 + T6p
K3	738.2448156	ab179185 - H3 K4me3 + K9ac
K4	43.67075	ab179186 - H3 K4me3 + K9me2
K5	75.43874479	ab179187 - H3 K4me3 + K9me3
K6	45.34421563	ab179188 - H3 K4me3 + S10p
K7	586.7489333	ab179189 - H3 K4me3 + K14ac
K8	1382.185565	ab179190 - H3 K4me3 + K18ac
K9	17.70669688	ab179191 - H3 K4me3 + R8me2s + K9me3
K10	20.74088958	ab179192 - H3 K4me3 + R8me2a + K9me3
K11	944.6769271	ab179193 - H3 K4me3 + K9ac + S10p
K12	1343.336435	ab179194 - H3 K4me3 + K9ac + K14ac
K13	2104.764865	ab179195 - H3 K4me3 + K9ac + K18ac
K14	1560.39723	ab179196 - H3 K4me3 + K14ac + K18ac
K15	2373.5127	ab179197 - H3 K4me3 + K9ac + K14ac + K18ac
K16	2092.145504	ab179198 - H3 K4me3 + K9ac + S10p + K14ac + K18ac
K17	2285.176241	ab179199 - H3 K4me3 + T6p + K9ac + K14ac + K18ac
K18	2096.281132	ab179200 - H3 K4me2 + T6p + K9ac + K14ac + K18ac
K19	73.35574688	ab179201 - H3 K4bio + K9bio + K18bio
K20	32.23819063	ab179202 - H3 T6p + K9me3
K21	30.98515521	ab179203 - H3 T6p + R8me2a + K9me3
K22	1646.78336	ab179204 - H3 T6p + K9ac + K14ac + K18ac
K23	316.8608438	ab179205 - H3 R8me1
K24	40.78090833	ab179206 - H3 R8me1 + K9me1
K25	96.42164792	ab179207 - H3 R8me1 + K9me2
L1	9.149023958	ab179208 - H3 R8me1 + K9me3
L2	28.89422604	ab179209 - H3 R8me2s
L3	13.96096146	ab179210 - H3 R8me2s + K9me1
L4	12.88545313	ab179211 - H3 R8me2s + K9me2
L5	6.5502125	ab179212 - H3 R8me2s + K9me3
L6	40.03115	ab179213 - H3 R8me2a
L7	17.242925	ab179214 - H3 R8me2a + K9me1
L8	14.74839167	ab179215 - H3 R8me2a + K9me2
L9	8.407121875	ab179216 - H3 R8me2a + K9me3
L10	14.40535833	ab179217 - H3 R8cit
L11	832.9548063	ab179218 - H3 K9ac + K14ac
L12	901.2676656	ab179219 - H3 K9ac + K18ac
L13	811.4525594	ab179220 - H3 K9ac + K14ac + K18ac
L14	2250.822435	ab179221 - H3 K9ac + K14ac + K18ac + K23ac + K27ac
L15	61.87360729	ab179222 - H3 K9ac + S10p
L16	533.9649344	ab179223 - H3 K9cr
L17	38.34328125	ab179224 - H3 K9me2

L18	15.33186146	ab179225 - H3 K9me2 + S10p
L19	31.56319375	ab179226 - H3 K9me2 + K27me2
L20	164.3253531	ab179227 - H3 S10p + S28p
L21	1482.559158	ab179228 - H3 S10p K14ac
L22	12.64852188	ab179229 - H3 K14me1
L23	13.45857292	ab179230 - H3 K14me2
L24	11.43642813	ab179231 - H3 K14me3
L25	726.8768552	ab179232 - H3 K14ac + K18ac
M1	15.77135208	ab179233 - H3 R17cit
M2	12.73162292	ab179234 - H3 R17me1
M3	13.86788854	ab179235 - H3 R17me2s
M4	14.01421979	ab179236 - H3 R17me2a + K18ac
M5	11.11081146	ab179237 - H3 K18cr
M6	13.46848438	ab179238 - H3 K18me1
M7	10.32561771	ab179239 - H3 K18me2
M8	276.6769219	ab179240 - H3 K18me3 + K36me3
M9	13.70000313	ab179241 - H3 K23cr
M10	17.66411042	ab179242 - H3 K23me1
M11	17.96160625	ab179243 - H3 K23me2
M12	14.41621146	ab179244 - H3 K23me3
M13	13.95967396	ab179245 - H3 T22p + K23me3
M14	21.60643229	ab179246 - H3 K23pr
M15	347.6539719	ab179247 - H3 unmod(20-51)
M16	82.00937708	ab179248 - H3 R26me1
M17	48.39518958	ab179249 - H3 R26me2a
M18	14.58606146	ab179250 - H3 R26me2a + K27me1
M19	11.77869063	ab179251 - H3 R26me2a + K27me3
M20	119.4183323	ab179252 - H3 K27ac + S28p
M21	135.4540604	ab179253 - H3 K27cr
M22	19.21315938	ab179254 - H3 K27me1
M23	16.72106979	ab179255 - H3 K27me2 + S28p
M24	10.15190313	ab179256 - H3 K27me3
M25	15.72930208	ab179257 - H3 K27me3 + S28p
N1	13.28227813	ab179258 - H3 P30hpro
N2	16.36246667	ab179259 - H3 P30(5,5me2)
N3	12.63936979	ab179260 - H3 P30
N4	14.30426042	ab179261 - H3 P30mutA
N5	11.90856458	ab179262 - H3 T32GlcNAc
N6	14.35502708	ab179263 - H3 K36ac
N7	13.67203542	ab179264 - H3 P38hpro
N8	11.69113542	ab179265 - H3 P38(5,5me2)
N9	18.98526563	ab179266 - H3 P38
N10	13.24436042	ab179267 - H3 P38mutA
N11	9.336334375	ab179268 - H3 Y41p
N12	17.07092813	ab179269 - H3 unmod(50-80)
N13	11.77589688	ab179270 - H3 K56ac
N14	13.2637375	ab179271 - H3 K56cr
N15	15.14293333	ab179272 - H3 K56fo
N16	10.42103125	ab179273 - H3 K56me1
N17	14.62286563	ab179274 - H3 K56me2
N18	11.71477813	ab179275 - H3 K56me3
N19	13.51948438	ab179276 - H3 R63me1
N20	10.14937604	ab179277 - H3 R63me2s
N21	13.46349063	ab179278 - H3 R63me2a
N22	13.56465938	ab179279 - H3 K64ac
N23	11.609225	ab179280 - H3 K64me2
N24	12.61525104	ab179281 - H3 K64me3
N25	11.91457813	ab179282 - H3 unmod(73-103)
O1	14.95742292	ab179283 - H3 K79ac
O2	13.063375	ab179284 - H3 K79fo
O3	6.6510625	ab179285 - H3 T80p
O4	13.77568333	ab179286 - H3 K79me3 + T80p
O5	12.34469271	ab179287 - H3 unmod(99-135)
O6	14.98413021	ab179288 - H3 T118O-ac
O7	13.26870208	ab179289 - H3 K122fo
O8	14.28571042	ab179290 - H3 K122me1
O9	12.47092188	ab179291 - H3 R128me1
O10	11.92946042	ab179292 - H3 R128me2s
O11	170.1829396	ab179293 - H3.3 unmod(10-40)
O12	11.23004063	ab179294 - H3.3 S28p
O13	15.97933021	ab179295 - H3.3 S28p + S31p
O14	12.50975521	ab179296 - H3.3 S31p
O15	10.68209583	ab179297 - H4 unmod(1-30)
O16	11.84501042	ab179298 - H4 S1p + R3me1

O17	14.7259	ab179299 - H4 S1p + R3me2s
O18	12.48291667	ab179300 - H4 S1p + R3me2a
O19	27.099125	ab179301 - H4 S1p + K5ac + K8ac + K12ac + K16ac
O20	11.29130625	ab179302 - H4 R3cit
O21	56.68974063	ab179303 - H4 R3me1 + K5ac + K8ac + K12ac + K16ac + K20ac
O22	11.23519792	ab179304 - H4 R3me2s
O23	49.07621042	ab179305 - H4 R3me2s + K5ac + K8ac + K12ac + K16ac + K20ac
O24	11.95474271	ab179306 - H4 R3me2a
O25	11.16461979	ab179307 - H4 R3me2a + K5ac
P1	177.7826656	ab179308 - H4 R3me2a + K5ac + K8ac + K12ac + K16ac + K20ac
P2	9.069522917	ab179309 - H4 K5ac
P3	16.37148125	ab179310 - H4 K5ac + K8ac
P4	23.39422292	ab179311 - H4 K5ac + K12ac
P5	14.33553229	ab179312 - H4 K5ac + K16ac
P6	31.28786979	ab179313 - H4 K5ac + K8ac + K12ac
P7	14.77748021	ab179314 - H4 K5ac + K8me1 + K12ac
P8	23.71702708	ab179315 - H4 K5ac + K8ac + K12ac + K16ac
P9	12.88634792	ab179316 - H4 K5ac + K8me1 + K12ac + K16ac
P10	19.23759583	ab179317 - H4 K5ac + K8ac + K12me1 + K16ac
P11	11.57502083	ab179318 - H4 K5me1
P12	11.07408958	ab179319 - H4 K5me1 + K8ac + K12me1
P13	12.32349167	ab179320 - H4 K5me1 + K8me1 + K12me1
P14	11.57995	ab179321 - H4 K5me1 + K8ac + K12ac + K16ac
P15	11.54096354	ab179322 - H4 K5me2
P16	12.65625	ab179323 - H4 K5me3
P17	12.13660625	ab179324 - H4 K5cr
P18	10.35699688	ab179325 - H4 K8ac + K12ac
P19	17.07515625	ab179326 - H4 K8ac + K16ac
P20	12.246575	ab179327 - H4 K8cr
P21	14.93159792	ab179328 - H4 K8me1
P22	10.29945	ab179329 - H4 K12ac
P23	16.41090521	ab179330 - H4 K12ac + K16ac
P24	15.7811125	ab179331 - H4 K12ac + K16ac + K20me2
P25	12.68583333	ab179332 - H4 K12ac + K16ac + K20me3
Q1	14.91530938	ab179333 - H4 K5mutQ + K8mutQ + K12ac + K16mutQ + K20mutQ
Q2	13.52202917	ab179334 - H4 K12cr
Q3	10.33032083	ab179335 - H4 K12fo
Q4	10.8080125	ab179336 - H4 K12me1
Q5	12.45114063	ab179337 - H4 K12me2
Q6	10.58747917	ab179338 - H4 K12me3
Q7	13.01085313	ab179339 - H4 K16ac
Q8	19.87755729	ab179340 - H4 K16ac + K20ac
Q9	11.04650833	ab179341 - H4 K16cr
Q10	12.71127604	ab179342 - H4 K16me1
Q11	15.41658021	ab179343 - H4 K16me3
Q12	11.25717396	ab179344 - H4 R17me1
Q13	12.97173854	ab179345 - H4 R17me2s
Q14	10.09855313	ab179346 - H4 R17me2a
Q15	9.918598958	ab179347 - H4 R19me1
Q16	10.24422396	ab179348 - H4 R19me2s
Q17	11.47681979	ab179349 - H4 R19me2a
Q18	16.47335625	ab179350 - H4 K20ac
Q19	11.77346771	ab179351 - H4 R23me1
Q20	13.41866667	ab179352 - H4 R23me2s
Q21	10.10810417	ab179353 - H4 R23me2a
Q22	14.13610938	ab179354 - H4 unmod(24-54)
Q23	13.74354792	ab179355 - H4 K31fo
Q24	10.16730729	ab179356 - H4 K31me1
Q25	13.87634792	ab179357 - H4 K31me2
R1	11.4281875	ab179358 - H4 K31me3
R2	11.05415208	ab179359 - H4 P32hyd
R3	11.47173438	ab179360 - H4 R35me1
R4	11.65874167	ab179361 - H4 R35me2s
R5	11.19678438	ab179362 - H4 R35me2a
R6	11.74259375	ab179363 - H4 S47O-ac
R7	10.68036563	ab179364 - H4 S47GlcNAc
R8	13.54817292	ab179365 - H4 S47p
R9	15.21596458	ab179366 - H4 unmod(45-75)
R10	10.59581875	ab179367 - H4 Y51hyd
R11	9.240975	ab179368 - H4 R55me1
R12	13.80213229	ab179369 - H4 R55me2s
R13	11.66437917	ab179370 - H4 R55me2a
R14	14.11928438	ab179371 - H4 K59fo
R15	14.21061146	ab179372 - H4 K59me1

R16	12.81602708	ab179373 - H4 K59me2
R17	10.02515938	ab179374 - H4 R67me1
R18	10.78974167	ab179375 - H4 R67me2s
R19	9.954559375	ab179376 - H4 R67me2a
R20	11.69679896	ab179377 - H4 unmod(71-102)
R21	10.19285313	ab179378 - H4 K77me1
R22	10.07767188	ab179379 - H4 K77me2
R23	12.34331563	ab179380 - H4 K77me3
R24	10.57933021	ab179381 - H4 K79fo
R25	13.09633958	ab179382 - H4 K79me1
S1	12.01548229	ab179383 - H4 K79me3
S2	12.2113	ab179384 - H4 T82O-ac
S3	11.55495417	ab179385 - H4 Y88hyd
S4	9.65055	ab179386 - H4 Y88nitro
S5	11.0263125	ab179387 - H4 K91fo
S6	10.48135208	ab179403 - H2A K5ac
S7	9.681767708	ab179404 - H2A K9me1
S8	13.84061146	ab18504 - H2B S32p
S9	10.01730625	ab19828 - ScH2A S129p
S10	77.45755625	ab20631 - H3 R26cit
S11	34.80182708	ab20635 - H3 T3p + K4me3
S12	13.06341563	ab21043 - H3 K37me1
S13	12.29365521	ab21998 - H1.4 K25me2
S14	12.70539792	ab22079 - H4 K59me3
S15	14.17463229	ab22080 - H3 K122me2
S16	12.07818021	ab22082 - H3 K122me3
S17	15.1313625	ab22214 - H3 K37me2
S18	11.19628229	ab22398 - H2A R3me2a
S19	8.3955875	ab22399 - H2A R3me2s
S20	15.97704271	ab23539 - H3 T22p
S21	550.6801875	ab24003 - H3 K18ac
S22	1661.016443	ab24404 - H3 K27ac
S23	11.72442188	ab24417 - H3 K37me3
S24	37.11206875	ab24444 - H3 T11p
S25	10.47555833	ab24659 - H2A K127me3
T1	13.10339375	ab24660 - H2A K125me3
T2	13.11503333	ab26307 - H3 T45p
T3	12.45419063	ab26342 - H2B K5me2
T4	9.043570833	ab26739 - H4 T30p
T5	10.35261458	ab27896 - H4 T96p
T6	10.97562813	ab27898 - H4 T80p
T7	12.41002292	ab28756 - H2A R3me1
T8	21.73934167	ab32876 - H3 R2cit + R8cit + R17cit
T9	13.47353229	ab33047 - H2A K127me2
T10	11.32346458	ab33505 - H3 T118p
T11	8.079015625	ab33594 - H3 R128me2a
T12	15.1255375	ab34466 - H3 K122ac
T13	11.94939167	ab36927 - H2A S122p
T14	10.88293958	ab4555 - H3 K79me1
T15	11.48314063	ab4556 - H3 K79me2
T16	11.11485104	ab4557 - H3 K79me3
T17	10.92731771	ab4560 - H4 K79me2
T18	11.85452813	ab46854 - H3 K18me3
T19	257.134351	ab48359 - H3 K23ac
T20	12.50791875	ab53244 - H3 K64me1
T21	15.16100313	ab54016 - H2A K9ac
T22	156.7091125	ab5499 - H3 S28p
T23	586.2948281	ab7768 - H3 K4me2
T24	16.42844479	ab89018 - H1.4 S26p
T25	474.5042354	ab92374 - H3 K4me3 (HEX1)
U1	9.397917708	ab95444 - H4 K16me2