

Peptide array: Histone H2A(symmetrical di methyl R3) + Histone H4 (symmetrical di methyl R3)(ab309354)
 Area under curve (AUC) data for all peptides on the pad. Antigen-containing peptides are highlighted in green.

Cell Ref	AUC	Peptide
A1	19.14833542	ab10112 - H4 K91ac
A2	18.58471771	ab10121 - H2B K85ac
A3	15.60381979	ab10127 - H2A K95me2
A4	12.99278021	ab10138 - H1.4 T17p
A5	12.03073125	ab10139 - H1.4 T146p
A6	20.2084	ab113997 - H2A.X Y142p
A7	16.53706979	ab11477 - H3 S10p
A8	18.38590313	ab12946 - H2B K5me3
A9	13.17216875	ab12949 - H2B K23me2
A10	18.74283854	ab12952 - H2A K127me1
A11	17.20389792	ab13211 - H2B K5me1
A12	17.27482188	ab1340 - H3 K4me1
A13	16.96113021	ab13570 - H2B K43me1
A14	17.48687917	ab13835 - HA
A15	17.68335938	ab13837 - Myc
A16	20.25134896	ab140408 - ScH2B T129p
A17	19.80237917	ab14103 - H3 T6p
A18	18.34832292	ab14724 - H4 S1p
A19	18.61479479	ab14799 - H3 T32p
A20	18.62996875	ab14943 - His
A21	21.08499688	ab14952 - H2A R77me2s
A22	17.94293021	ab14964 - H4 K20me2
A23	16.93868854	ab14997 - ScH2A K21ac
A24	21.16318125	ab154782 - H3 R26me2s
A25	19.59250417	ab15591 - H3 K14ac
B1	16.00229896	ab15644 - H3 K9me3 + S10p
B2	20.97079271	ab15645 - H2A.X S139p
B3	18.9955125	ab15824 - H4 K8ac
B4	25.6502375	ab15829 - V5
B5	18.45204792	ab16635 - H3 K9ac
B6	18.73279063	ab166685 - H2B S112GlcNAc
B7	23.94018229	ab166688 - H2B S36GlcNAc
B8	21.90192917	ab16935 - H3 R17me2a
B9	22.26805521	ab17043 - H4 K20me1
B10	23.96479688	ab17162 - H3 P16hyd + P30hyd
B11	20.76886146	ab17567 - H4 K20me3
B12	29.06597188	ab17587 - H1.4 K25me3
B13	29.37441458	ab175905 - H2B K11ac
B14	24.07836979	ab176211 - H2B K23ac
B15	17.44623958	ab17632 - H4 K5bio
B16	19.79261771	ab1771 - H3 K9me1
B17	17.52661771	ab1773 - H3 K9me3
B18	18.59332083	ab17770 - H4 R3me1
B19	17.71945	ab1781 - H3 K27me2
B20	19.6439875	ab1783 - H3 K36me1
B21	21.07773854	ab1784 - H3 K36me2
B22	22.36654896	ab1785 - H3 K36me3
B23	28.2260625	ab178667 - ScH2A unmod(1-30)
B24	21.07616354	ab178668 - ScH2A unmod(102-132)
B25	28.82891771	ab178670 - H2A.Z unmod(1-30)
C1	20.47767292	ab178833 - Flag
C2	23.49654896	ab178835 - H2B R99mutO
C3	20.13448542	ab178836 - H4 R23mutO
C4	18.27360104	ab178837 - H4 G48mutD
C5	18.92154167	ab178987 - H1.4 unmod(1-30)
C6	12.59302292	ab178988 - H1.4 S1p
C7	11.83688021	ab178989 - H1.4 T3p
C8	12.81394792	ab178990 - H1.4 S1p + T3p
C9	17.56539583	ab178991 - H1.4 K25me1
C10	18.17163333	ab178992 - H1.4 K25me3 + S26p
C11	15.28861563	ab178993 - H1.4 unmod(130-160)
C12	24.913275	ab178994 - H2A unmod(1-30)
C13	20.24839688	ab178995 - H2A S1p
C14	19.28869479	ab178996 - H2A S1p + K5ac
C15	17.44912188	ab178997 - H2A S1p + R3me2a + K5ac
C16	5085.506793	ab178998 - H2A S1p + R3me2s + K5ac
C17	21.9875875	ab178999 - H2A S1p + R3cit + K5ac
C18	28.38675938	ab179000 - H2A S1ac + S1p + K5ac + K9ac + K13ac + K15ac
C19	15.55497813	ab179001 - H2A S1p + K5ac + K9ac + K13ac + K15ac
C20	193.6028635	ab179002 - H2A R3cit

C21	31.24868958	ab179003 - H2A R3cit + K5ac
C22	20.28532292	ab179004 - H2A R3me2a + K5ac
C23	3706.979975	ab179005 - H2A R3me2s + K5ac
C24	74.15755729	ab179006 - H2A S1ac + K5ac + K9ac + K13ac + K15ac
C25	21.52325	ab179007 - H2A K5ac + K9ac + K13ac + K15ac
D1	17.89614583	ab179008 - H2A K9me2
D2	13.25198229	ab179009 - H2A K9me3
D3	15.18059583	ab179010 - H2A R11me1
D4	13.55920521	ab179011 - H2A R11me2s
D5	16.56616146	ab179012 - H2A R11me2a
D6	12.89790833	ab179013 - H2A K13ac
D7	17.16364896	ab179014 - H2A K15ac
D8	12.00436667	ab179015 - H2A R17me1
D9	20.05049688	ab179016 - H2A R17me2s
D10	15.45187292	ab179017 - H2A R17me2a
D11	18.91807083	ab179018 - H2A unmod(20-50)
D12	13.26660625	ab179019 - H2A P26hyd
D13	19.1761125	ab179020 - H2A R29me1
D14	11.47291563	ab179021 - H2A R29me2s
D15	16.43443229	ab179022 - H2A R29me2a
D16	13.46191458	ab179023 - H2A K36ac
D17	17.23505208	ab179024 - H2A K36cr
D18	19.11358021	ab179025 - H2A Y39hyd
D19	22.96567396	ab179026 - H2A R42me1
D20	15.65575208	ab179027 - H2A R42me2s
D21	19.50449271	ab179028 - H2A R42me2a
D22	17.28951042	ab179029 - H2A unmod(71-103)
D23	22.37167813	ab179030 - H2A R77me1
D24	20.85254063	ab179031 - H2A R77me2a
D25	17.48977396	ab179032 - H2A R88me1
E1	22.48315625	ab179033 - H2A R88me2s
E2	16.10573542	ab179034 - H2A R88me2a
E3	22.65277188	ab179035 - H2A K95me1
E4	20.06847708	ab179036 - H2A K95me3
E5	25.25019271	ab179037 - H2A unmod(92-123)
E6	15.36690104	ab179038 - H2A T101GlcNAc
E7	13.42274167	ab179039 - H2A Q104me1
E8	17.40971146	ab179040 - H2A unmod(112-129)
E9	21.28991458	ab179041 - H2A K118me1
E10	13.08069479	ab179042 - H2A K118me2
E11	12.60525938	ab179043 - H2A K118me3
E12	17.21935938	ab179044 - H2A K118fo
E13	14.68953333	ab179045 - H2A K118cr
E14	18.64317188	ab179046 - H2A K119cr
E15	14.00838438	ab179047 - H2A K118ubiq
E16	16.74506042	ab179048 - H2A K119ubiq
E17	25.68165417	ab179049 - H2A T120p
E18	12.95170833	ab179050 - H2A K125me1
E19	18.98427083	ab179051 - H2A K125me2
E20	22.63179896	ab179052 - H2A K125cr
E21	50.21416875	ab179053 - H2A.X unmod(1-30)
E22	17.55716979	ab179054 - H2A.X K5ac
E23	21.54196875	ab179055 - H2A.X unmod(114-142)
E24	19.12811458	ab179056 - H2A.X K118ubiq
E25	20.50640521	ab179057 - H2A.X K119ubiq
F1	21.65314271	ab179058 - H2A.X T120p
F2	19.10886458	ab179059 - H2A.X K127ubiq
F3	20.95425417	ab179060 - H2A.X S139p + Y142p
F4	22.60064063	ab179061 - H2A.Z K4ac
F5	19.89098125	ab179062 - H2A.Z K7ac
F6	18.38343646	ab179063 - H2A.Z K11ac
F7	17.68425938	ab179064 - H2A.Z K4ac + K7ac + K11ac
F8	49.54609583	ab179065 - H2B unmod(1-30)
F9	16.88232813	ab179066 - H2B K5ac
F10	20.06078854	ab179067 - H2B K5cr
F11	15.86651146	ab179068 - H2B K5fo
F12	13.54638958	ab179069 - H2B P10hyd
F13	15.11983542	ab179070 - H2B K11me1
F14	24.03178958	ab179071 - H2B K11me2
F15	13.97935208	ab179072 - H2B K11me3
F16	21.77407917	ab179073 - H2B K11cr
F17	16.00972604	ab179074 - H2B K12ac
F18	20.96295208	ab179075 - H2B K12cr
F19	17.16448021	ab179076 - H2B K12me1

F20	21.0282875	ab179077 - H2B K12me2
F21	19.58930521	ab179078 - H2B K12me3
F22	18.47210521	ab179079 - H2B K12ac + K15ac
F23	18.08320938	ab179080 - H2B S14p
F24	15.57940625	ab179081 - H2B K15ac
F25	18.29511875	ab179082 - H2B K15cr
G1	15.66703438	ab179083 - H2B K15me1
G2	19.28458333	ab179084 - H2B K15me2
G3	16.31289583	ab179085 - H2B K15me3
G4	16.41775208	ab179086 - H2B K16ac
G5	14.57324792	ab179087 - H2B K16cr
G6	17.94902604	ab179088 - H2B K20ac
G7	22.78226563	ab179089 - H2B K20cr
G8	14.58227604	ab179090 - H2B K20me1
G9	21.76055208	ab179091 - H2B K20me2
G10	16.11456458	ab179092 - H2B K20me3
G11	19.56560729	ab179093 - H2B K23cr
G12	20.81884792	ab179094 - H2B K23me1
G13	17.52968438	ab179095 - H2B K23me3
G14	18.47743021	ab179096 - H2B K24ac
G15	18.35279896	ab179097 - H2B unmod(26-56)
G16	20.31671042	ab179098 - H2B K34cr
G17	15.83199375	ab179099 - H2B K34fo
G18	15.48701979	ab179100 - H2B Y37hyd
G19	22.21381979	ab179101 - H2B Y37p
G20	20.94360417	ab179102 - H2B K43me2
G21	15.31875104	ab179103 - H2B K43me3
G22	23.81300521	ab179104 - H2B K46fo
G23	19.9132125	ab179105 - H2B K46ac
G24	14.71702604	ab179106 - H2B K46me1
G25	17.93489271	ab179107 - H2B K46me2
H1	14.38998021	ab179108 - H2B K46me3
H2	14.24880521	ab179109 - H2B unmod(51-81)
H3	16.83020521	ab179110 - H2B K57me1
H4	25.38127708	ab179111 - H2B K57me2
H5	15.20426979	ab179112 - H2B K57me3
H6	18.92510833	ab179113 - H2B unmod(73-103)
H7	22.45565833	ab179114 - H2B R79me1
H8	17.94653021	ab179115 - H2B R79me2s
H9	18.23966875	ab179116 - H2B R79me2a
H10	15.36655	ab179117 - H2B Y83hyd
H11	16.529625	ab179118 - H2B K85me1
H12	23.64814688	ab179119 - H2B K85me2
H13	16.52762083	ab179120 - H2B K85me3
H14	23.35283542	ab179121 - H2B unmod(94-125)
H15	17.21220833	ab179122 - H2B R99me1
H16	20.01959688	ab179123 - H2B R99me2s
H17	16.03027396	ab179124 - H2B R99me2a
H18	15.35847917	ab179125 - H2B K108fo
H19	18.66660104	ab179126 - H2B K116fo
H20	22.64872708	ab179127 - H2B K116me1
H21	18.13225833	ab179128 - H2B K116me2
H22	21.3149125	ab179129 - H2B K116me3
H23	17.98523438	ab179130 - H2B K116ubiq
H24	23.84944896	ab179131 - H2B K120ac
H25	14.872975	ab179132 - H2B K120fo
I1	31.29217083	ab179133 - H2B K120ubiq
I2	29.10942396	ab179134 - H2B K125ubiq
I3	17.72070625	ab179135 - ScH2B unmod(102-132)
I4	19.85676146	ab179136 - H3 unmod(1-33)
I5	17.73381042	ab179137 - H3 R2cit
I6	16.59560729	ab179138 - H3 R2cit + K4me2
I7	20.30219271	ab179139 - H3 R2cit + K4me3
I8	26.33168958	ab179140 - H3 R2cit + K4me3 + K9ac + K14ac + K18ac
I9	16.2630125	ab179141 - H3 R2cit + K4ac + K9ac + K14ac + K18ac
I10	15.10838542	ab179142 - H3 R2me1
I11	17.49092917	ab179143 - H3 R2me1 + K4me2
I12	18.20468646	ab179144 - H3 R2me1 + K4me3
I13	17.34862292	ab179145 - H3 R2me1 + K4me3 + K9ac + K14ac + K18ac
I14	16.55312708	ab179146 - H3 R2me2s
I15	17.07773958	ab179147 - H3 R2me2s + K4me2
I16	17.42422292	ab179148 - H3 R2me2s + K4me3
I17	18.01255	ab179149 - H3 R2me2s + K4me3 + K9ac + K14ac + K18ac
I18	18.52859063	ab179150 - H3 R2me2a

I19	18.41384583	ab179151 - H3 R2me2a + K4me2
I20	16.31053333	ab179152 - H3 R2me2a + K4me3
I21	14.72359063	ab179153 - H3 R2me2a + K4me3 + S10p
I22	18.86106042	ab179154 - H3 R2me2a + T3p + K4me3
I23	16.54306042	ab179155 - H3 R2me2a + K4ac + K9ac + K14ac + K18ac
I24	21.39490833	ab179156 - H3 R2me2a + K4me3 + K9ac + K14ac + K18ac
I25	18.85509479	ab179157 - H3 R2me2a + T3p + K4me3 + K9ac + K14ac + K18ac
J1	16.05053958	ab179158 - H3 R2me2a + K4me3 + K9ac + S10p + K14ac + K18ac
J2	17.3900625	ab179159 - H3 T3p
J3	17.49144688	ab179160 - H3 T3p + K4me1
J4	18.20319896	ab179161 - H3 T3p + K4me2
J5	20.59641146	ab179162 - H3 T3p + K4me3 + K9ac + K14ac + K18ac
J6	15.26642292	ab179163 - H3 K4ac
J7	13.61720729	ab179164 - H3 K4ac + K9ac
J8	15.42639375	ab179165 - H3 K4ac + K14ac
J9	20.26171458	ab179166 - H3 K4ac + K18ac
J10	18.49195625	ab179167 - H3 K4ac + K9me3
J11	15.74201146	ab179168 - H3 K4ac + K9ac + K14ac
J12	15.74288542	ab179169 - H3 K4ac + K9ac + K18ac
J13	18.54308958	ab179170 - H3 K4ac + K14ac + K18ac
J14	16.3058	ab179171 - H3 K4ac + K9ac + K14ac + K18ac
J15	19.60668646	ab179172 - H3 K4ac + K9me1 + K14ac + K18ac
J16	16.39387604	ab179173 - H3 K4ac + K9me2 + K14ac + K18ac
J17	18.09627396	ab179174 - H3 K4ac + K9me3 + K14ac + K18ac
J18	16.03267292	ab179175 - H3 K4ac + K9ac + S10p + K14ac + K18ac
J19	16.01095521	ab179176 - H3 K4cr
J20	19.91685104	ab179177 - H3 K4me1 + K9me2
J21	20.27668646	ab179178 - H3 K4me1 + K9ac + K18ac
J22	18.03923021	ab179179 - H3 K4me1 + K9ac + K14ac + K18ac
J23	18.59904063	ab179180 - H3 K4me2 + T6p
J24	19.31750104	ab179181 - H3 K4me2 + K9me2
J25	18.93822292	ab179182 - H3 K4me2 + K9ac + K18ac
K1	23.97591979	ab179183 - H3 K4me2 + K9ac + K14ac + K18ac
K2	13.63919271	ab179184 - H3 K4me3 + T6p
K3	13.98460938	ab179185 - H3 K4me3 + K9ac
K4	16.78380625	ab179186 - H3 K4me3 + K9me2
K5	17.87696875	ab179187 - H3 K4me3 + K9me3
K6	17.55825208	ab179188 - H3 K4me3 + S10p
K7	14.11274896	ab179189 - H3 K4me3 + K14ac
K8	17.13756875	ab179190 - H3 K4me3 + K18ac
K9	13.68749896	ab179191 - H3 K4me3 + R8me2s + K9me3
K10	12.27592188	ab179192 - H3 K4me3 + R8me2a + K9me3
K11	20.02760729	ab179193 - H3 K4me3 + K9ac + S10p
K12	19.57037188	ab179194 - H3 K4me3 + K9ac + K14ac
K13	21.0369	ab179195 - H3 K4me3 + K9ac + K18ac
K14	18.53852292	ab179196 - H3 K4me3 + K14ac + K18ac
K15	20.96296354	ab179197 - H3 K4me3 + K9ac + K14ac + K18ac
K16	18.13135104	ab179198 - H3 K4me3 + K9ac + S10p + K14ac + K18ac
K17	17.37693438	ab179199 - H3 K4me3 + T6p + K9ac + K14ac + K18ac
K18	18.90281563	ab179200 - H3 K4me2 + T6p + K9ac + K14ac + K18ac
K19	22.34497083	ab179201 - H3 K4bio + K9bio + K18bio
K20	16.32998021	ab179202 - H3 T6p + K9me3
K21	23.24571563	ab179203 - H3 T6p + R8me2a + K9me3
K22	21.67982708	ab179204 - H3 T6p + K9ac + K14ac + K18ac
K23	20.30125417	ab179205 - H3 R8me1
K24	18.87066771	ab179206 - H3 R8me1 + K9me1
K25	17.16421354	ab179207 - H3 R8me1 + K9me2
L1	17.32643542	ab179208 - H3 R8me1 + K9me3
L2	13.83210938	ab179209 - H3 R8me2s
L3	14.3314375	ab179210 - H3 R8me2s + K9me1
L4	16.71499063	ab179211 - H3 R8me2s + K9me2
L5	14.60543958	ab179212 - H3 R8me2s + K9me3
L6	16.36668229	ab179213 - H3 R8me2a
L7	15.56736458	ab179214 - H3 R8me2a + K9me1
L8	20.12835729	ab179215 - H3 R8me2a + K9me2
L9	15.99788646	ab179216 - H3 R8me2a + K9me3
L10	16.04939792	ab179217 - H3 R8cit
L11	15.86540313	ab179218 - H3 K9ac + K14ac
L12	15.79500104	ab179219 - H3 K9ac + K18ac
L13	18.99531146	ab179220 - H3 K9ac + K14ac + K18ac
L14	24.53600938	ab179221 - H3 K9ac + K14ac + K18ac + K23ac + K27ac
L15	16.59391667	ab179222 - H3 K9ac + S10p
L16	19.69330521	ab179223 - H3 K9cr
L17	19.282425	ab179224 - H3 K9me2

L18	20.56752917	ab179225 - H3 K9me2 + S10p
L19	14.42434271	ab179226 - H3 K9me2 + K27me2
L20	16.05449688	ab179227 - H3 S10p + S28p
L21	17.99028438	ab179228 - H3 S10p K14ac
L22	19.60158229	ab179229 - H3 K14me1
L23	20.80174792	ab179230 - H3 K14me2
L24	23.67892083	ab179231 - H3 K14me3
L25	20.29754271	ab179232 - H3 K14ac + K18ac
M1	20.07599896	ab179233 - H3 R17cit
M2	21.02721979	ab179234 - H3 R17me1
M3	16.96809479	ab179235 - H3 R17me2s
M4	14.66701563	ab179236 - H3 R17me2a + K18ac
M5	14.07534583	ab179237 - H3 K18cr
M6	18.65737813	ab179238 - H3 K18me1
M7	17.44956354	ab179239 - H3 K18me2
M8	21.90243958	ab179240 - H3 K18me3 + K36me3
M9	20.01625938	ab179241 - H3 K23cr
M10	17.88737708	ab179242 - H3 K23me1
M11	16.11251042	ab179243 - H3 K23me2
M12	18.53469375	ab179244 - H3 K23me3
M13	17.36082083	ab179245 - H3 T22p + K23me3
M14	18.93992396	ab179246 - H3 K23pr
M15	26.91696563	ab179247 - H3 unmod(20-51)
M16	22.74227083	ab179248 - H3 R26me1
M17	23.99315208	ab179249 - H3 R26me2a
M18	19.22752396	ab179250 - H3 R26me2a + K27me1
M19	25.30957708	ab179251 - H3 R26me2a + K27me3
M20	16.06528021	ab179252 - H3 K27ac + S28p
M21	16.85911667	ab179253 - H3 K27cr
M22	16.37966979	ab179254 - H3 K27me1
M23	16.65673854	ab179255 - H3 K27me2 + S28p
M24	18.18669271	ab179256 - H3 K27me3
M25	13.64350313	ab179257 - H3 K27me3 + S28p
N1	15.54673854	ab179258 - H3 P30hpro
N2	16.28762396	ab179259 - H3 P30(5,5me2)
N3	14.82426146	ab179260 - H3 P30
N4	18.51536146	ab179261 - H3 P30mutA
N5	17.46228021	ab179262 - H3 T32GlcNAc
N6	18.82795417	ab179263 - H3 K36ac
N7	19.45495833	ab179264 - H3 P38hpro
N8	19.80533854	ab179265 - H3 P38(5,5me2)
N9	17.97726667	ab179266 - H3 P38
N10	15.1342	ab179267 - H3 P38mutA
N11	19.12122917	ab179268 - H3 Y41p
N12	17.73787813	ab179269 - H3 unmod(50-80)
N13	14.92329271	ab179270 - H3 K56ac
N14	18.30390729	ab179271 - H3 K56cr
N15	16.31622708	ab179272 - H3 K56fo
N16	16.04111875	ab179273 - H3 K56me1
N17	17.69330313	ab179274 - H3 K56me2
N18	13.88317604	ab179275 - H3 K56me3
N19	21.64844583	ab179276 - H3 R63me1
N20	13.96308229	ab179277 - H3 R63me2s
N21	20.82024896	ab179278 - H3 R63me2a
N22	23.04721146	ab179279 - H3 K64ac
N23	19.67146146	ab179280 - H3 K64me2
N24	26.28602292	ab179281 - H3 K64me3
N25	20.06194688	ab179282 - H3 unmod(73-103)
O1	25.85708333	ab179283 - H3 K79ac
O2	18.90994688	ab179284 - H3 K79fo
O3	16.36844896	ab179285 - H3 T80p
O4	17.67932188	ab179286 - H3 K79me3 + T80p
O5	18.15634896	ab179287 - H3 unmod(99-135)
O6	14.57028438	ab179288 - H3 T118O-ac
O7	17.74378646	ab179289 - H3 K122fo
O8	15.68422604	ab179290 - H3 K122me1
O9	16.87477396	ab179291 - H3 R128me1
O10	19.70149271	ab179292 - H3 R128me2s
O11	19.61390208	ab179293 - H3.3 unmod(10-40)
O12	21.31991875	ab179294 - H3.3 S28p
O13	19.99878958	ab179295 - H3.3 S28p + S31p
O14	19.06731667	ab179296 - H3.3 S31p
O15	19.26755104	ab179297 - H4 unmod(1-30)
O16	46.01401146	ab179298 - H4 S1p + R3me1

O17	4382.734455	ab179299 - H4 S1p + R3me2s
O18	17.36252917	ab179300 - H4 S1p + R3me2a
O19	24.46975833	ab179301 - H4 S1p + K5ac + K8ac + K12ac + K16ac
O20	18.01274167	ab179302 - H4 R3cit
O21	17.28693021	ab179303 - H4 R3me1 + K5ac + K8ac + K12ac + K16ac + K20ac
O22	4666.709179	ab179304 - H4 R3me2s
O23	2392.588605	ab179305 - H4 R3me2s + K5ac + K8ac + K12ac + K16ac + K20ac
O24	60.25896771	ab179306 - H4 R3me2a
O25	136.2211719	ab179307 - H4 R3me2a + K5ac
P1	22.85069271	ab179308 - H4 R3me2a + K5ac + K8ac + K12ac + K16ac + K20ac
P2	17.143925	ab179309 - H4 K5ac
P3	21.29738438	ab179310 - H4 K5ac + K8ac
P4	20.46697292	ab179311 - H4 K5ac + K12ac
P5	17.5580625	ab179312 - H4 K5ac + K16ac
P6	16.76675938	ab179313 - H4 K5ac + K8ac + K12ac
P7	27.88877396	ab179314 - H4 K5ac + K8me1 + K12ac
P8	15.37124479	ab179315 - H4 K5ac + K8ac + K12ac + K16ac
P9	18.36182396	ab179316 - H4 K5ac + K8me1 + K12ac + K16ac
P10	17.14907396	ab179317 - H4 K5ac + K8ac + K12me1 + K16ac
P11	16.49736979	ab179318 - H4 K5me1
P12	21.30636563	ab179319 - H4 K5me1 + K8ac + K12me1
P13	18.56831771	ab179320 - H4 K5me1 + K8me1 + K12me1
P14	18.14873021	ab179321 - H4 K5me1 + K8ac + K12ac + K16ac
P15	17.47469063	ab179322 - H4 K5me2
P16	17.06286458	ab179323 - H4 K5me3
P17	21.40404688	ab179324 - H4 K5cr
P18	18.95862917	ab179325 - H4 K8ac + K12ac
P19	21.14994063	ab179326 - H4 K8ac + K16ac
P20	15.45383542	ab179327 - H4 K8cr
P21	20.37429583	ab179328 - H4 K8me1
P22	20.65535313	ab179329 - H4 K12ac
P23	18.75367396	ab179330 - H4 K12ac + K16ac
P24	18.33188542	ab179331 - H4 K12ac + K16ac + K20me2
P25	20.63657813	ab179332 - H4 K12ac + K16ac + K20me3
Q1	21.45625521	ab179333 - H4 K5mutQ + K8mutQ + K12ac + K16mutQ + K20mutQ
Q2	18.90458021	ab179334 - H4 K12cr
Q3	20.84464063	ab179335 - H4 K12fo
Q4	17.58581771	ab179336 - H4 K12me1
Q5	20.54875938	ab179337 - H4 K12me2
Q6	21.86816146	ab179338 - H4 K12me3
Q7	21.53615104	ab179339 - H4 K16ac
Q8	20.71539688	ab179340 - H4 K16ac + K20ac
Q9	18.31369688	ab179341 - H4 K16cr
Q10	18.60761979	ab179342 - H4 K16me1
Q11	20.10333542	ab179343 - H4 K16me3
Q12	22.0609	ab179344 - H4 R17me1
Q13	24.11318438	ab179345 - H4 R17me2s
Q14	16.13557292	ab179346 - H4 R17me2a
Q15	22.09269271	ab179347 - H4 R19me1
Q16	19.74078438	ab179348 - H4 R19me2s
Q17	24.07239167	ab179349 - H4 R19me2a
Q18	19.96933125	ab179350 - H4 K20ac
Q19	15.7715375	ab179351 - H4 R23me1
Q20	17.76059792	ab179352 - H4 R23me2s
Q21	17.92873021	ab179353 - H4 R23me2a
Q22	19.3852375	ab179354 - H4 unmod(24-54)
Q23	17.64347396	ab179355 - H4 K31fo
Q24	17.90755104	ab179356 - H4 K31me1
Q25	31.73266354	ab179357 - H4 K31me2
R1	20.52299375	ab179358 - H4 K31me3
R2	21.36545833	ab179359 - H4 P32hyd
R3	17.76472604	ab179360 - H4 R35me1
R4	22.5312125	ab179361 - H4 R35me2s
R5	17.36339479	ab179362 - H4 R35me2a
R6	17.80718542	ab179363 - H4 S47O-ac
R7	19.91980313	ab179364 - H4 S47GlcNAc
R8	21.54968854	ab179365 - H4 S47p
R9	19.06484479	ab179366 - H4 unmod(45-75)
R10	21.41373854	ab179367 - H4 Y51hyd
R11	18.48428646	ab179368 - H4 R55me1
R12	15.35571875	ab179369 - H4 R55me2s
R13	15.54638958	ab179370 - H4 R55me2a
R14	22.34898333	ab179371 - H4 K59fo
R15	15.27156146	ab179372 - H4 K59me1

R16	23.91747604	ab179373 - H4 K59me2
R17	26.64506458	ab179374 - H4 R67me1
R18	22.94523542	ab179375 - H4 R67me2s
R19	24.79175	ab179376 - H4 R67me2a
R20	23.3903	ab179377 - H4 unmod(71-102)
R21	23.58100313	ab179378 - H4 K77me1
R22	19.96195104	ab179379 - H4 K77me2
R23	16.63541875	ab179380 - H4 K77me3
R24	18.16696146	ab179381 - H4 K79fo
R25	17.31134688	ab179382 - H4 K79me1
S1	20.91796771	ab179383 - H4 K79me3
S2	15.70463958	ab179384 - H4 T82O-ac
S3	16.61416771	ab179385 - H4 Y88hyd
S4	19.549175	ab179386 - H4 Y88nitro
S5	18.89159792	ab179387 - H4 K91fo
S6	18.84630417	ab179403 - H2A K5ac
S7	20.87327083	ab179404 - H2A K9me1
S8	24.40836667	ab18504 - H2B S32p
S9	16.87736875	ab19828 - ScH2A S129p
S10	20.78177604	ab20631 - H3 R26cit
S11	19.48839688	ab20635 - H3 T3p + K4me3
S12	19.36583021	ab21043 - H3 K37me1
S13	12.82169479	ab21998 - H1.4 K25me2
S14	23.249125	ab22079 - H4 K59me3
S15	20.23826458	ab22080 - H3 K122me2
S16	16.42662292	ab22082 - H3 K122me3
S17	20.55652604	ab22214 - H3 K37me2
S18	24.43159583	ab22398 - H2A R3me2a
S19	4045.365734	ab22399 - H2A R3me2s
S20	22.14098333	ab23539 - H3 T22p
S21	16.29154271	ab24003 - H3 K18ac
S22	20.49134375	ab24404 - H3 K27ac
S23	19.67653958	ab24417 - H3 K37me3
S24	20.11607917	ab24444 - H3 T11p
S25	14.83342292	ab24659 - H2A K127me3
T1	16.97743958	ab24660 - H2A K125me3
T2	16.91661667	ab26307 - H3 T45p
T3	14.30225729	ab26342 - H2B K5me2
T4	19.4265	ab26739 - H4 T30p
T5	22.23257917	ab27896 - H4 T96p
T6	16.79089167	ab27898 - H4 T80p
T7	278.1768531	ab28756 - H2A R3me1
T8	15.96031667	ab32876 - H3 R2cit + R8cit + R17cit
T9	18.86716146	ab33047 - H2A K127me2
T10	18.88565625	ab33505 - H3 T118p
T11	18.09547708	ab33594 - H3 R128me2a
T12	17.79883021	ab34466 - H3 K122ac
T13	18.53266979	ab36927 - H2A S122p
T14	25.28997604	ab4555 - H3 K79me1
T15	23.90932708	ab4556 - H3 K79me2
T16	23.03515208	ab4557 - H3 K79me3
T17	18.26293333	ab4560 - H4 K79me2
T18	15.28561875	ab46854 - H3 K18me3
T19	23.2974875	ab48359 - H3 K23ac
T20	22.97331563	ab53244 - H3 K64me1
T21	39.98533958	ab54016 - H2A K9ac
T22	18.74212604	ab5499 - H3 S28p
T23	20.44222813	ab7768 - H3 K4me2
T24	13.75871771	ab89018 - H1.4 S26p
T25	20.8405	ab92374 - H3 K4me3 (HEX1)
U1	21.81558854	ab95444 - H4 K16me2