

Peptide array: Histone H2A.X (ab322655)

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

Cell Ref	AUC	Peptide
A1	22.55538125	ab10112 - H4 K91ac
A2	21.79534063	ab10121 - H2B K85ac
A3	19.03461458	ab10127 - H2A K95me2
A4	13.55506667	ab10138 - H1.4 T17p
A5	20.08466354	ab10139 - H1.4 T146p
A6	19.3918875	ab113997 - H2A.X Y142p
A7	23.52391667	ab11477 - H3 S10p
A8	22.77532083	ab12946 - H2B K5me3
A9	19.04850729	ab12949 - H2B K23me2
A10	18.8702875	ab12952 - H2A K127me1
A11	18.57007708	ab13211 - H2B K5me1
A12	23.96734896	ab1340 - H3 K4me1
A13	21.87417813	ab13570 - H2B K43me1
A14	20.00605833	ab13835 - HA
A15	19.32398854	ab13837 - Myc
A16	23.45226771	ab140408 - ScH2B T129p
A17	23.86217083	ab14103 - H3 T6p
A18	20.93570521	ab14724 - H4 S1p
A19	22.33772083	ab14799 - H3 T32p
A20	21.99256875	ab14943 - His
A21	21.51212917	ab14952 - H2A R77me2s
A22	17.40593229	ab14964 - H4 K20me2
A23	20.94584688	ab14997 - ScH2A K21ac
A24	23.02836875	ab154782 - H3 R26me2s
A25	21.57946667	ab15591 - H3 K14ac
B1	21.15477188	ab15644 - H3 K9me3 + S10p
B2	18.99417292	ab15645 - H2A.X S139p
B3	17.77525417	ab15824 - H4 K8ac
B4	22.08767396	ab15829 - V5
B5	14.19980938	ab16635 - H3 K9ac
B6	30.165425	ab166685 - H2B S112GlcNAc
B7	20.38993854	ab166688 - H2B S36GlcNAc
B8	22.80417396	ab16935 - H3 R17me2a
B9	23.84860208	ab17043 - H4 K20me1
B10	24.28316667	ab17162 - H3 P16hyd + P30hyd
B11	25.38872188	ab17567 - H4 K20me3
B12	13.72211563	ab17587 - H1.4 K25me3
B13	23.10321354	ab175905 - H2B K11ac
B14	23.58484792	ab176211 - H2B K23ac
B15	15.54274167	ab17632 - H4 K5bio
B16	14.20837188	ab1771 - H3 K9me1
B17	17.28237813	ab1773 - H3 K9me3
B18	23.9489	ab17770 - H4 R3me1
B19	19.34204063	ab1781 - H3 K27me2
B20	24.45703229	ab1783 - H3 K36me1
B21	23.93389375	ab1784 - H3 K36me2
B22	23.71125625	ab1785 - H3 K36me3
B23	4191.450966	ab178667 - ScH2A unmod(1-30)
B24	24.41288854	ab178668 - ScH2A unmod(102-132)
B25	11253.57599	ab178670 - H2A.Z unmod(1-30)
C1	23.8405125	ab178833 - Flag
C2	44.11266771	ab178835 - H2B R99mutO
C3	17.70614375	ab178836 - H4 R23mutO
C4	22.28637604	ab178837 - H4 G48mutD
C5	18.81718958	ab178987 - H1.4 unmod(1-30)
C6	14.47755938	ab178988 - H1.4 S1p
C7	15.73953333	ab178989 - H1.4 T3p
C8	12.12859375	ab178990 - H1.4 S1p + T3p
C9	20.73142917	ab178991 - H1.4 K25me1
C10	19.42057083	ab178992 - H1.4 K25me3 + S26p
C11	21.41261458	ab178993 - H1.4 unmod(130-160)
C12	6095.295558	ab178994 - H2A unmod(1-30)
C13	24.14687917	ab178995 - H2A S1p
C14	27.83837083	ab178996 - H2A S1p + K5ac
C15	24.26410625	ab178997 - H2A S1p + R3me2a + K5ac
C16	23.20030313	ab178998 - H2A S1p + R3me2s + K5ac
C17	19.45662396	ab178999 - H2A S1p + R3cit + K5ac
C18	22.89499167	ab179000 - H2A S1ac + S1p + K5ac + K9ac + K13ac + K15ac
C19	22.13930521	ab179001 - H2A S1p + K5ac + K9ac + K13ac + K15ac
C20	27.32406458	ab179002 - H2A R3cit

C21	25.34471042	ab179003 - H2A R3cit + K5ac
C22	18.83823229	ab179004 - H2A R3me2a + K5ac
C23	24.03601979	ab179005 - H2A R3me2s + K5ac
C24	16.61799896	ab179006 - H2A S1ac + K5ac + K9ac + K13ac + K15ac
C25	30.14791667	ab179007 - H2A K5ac + K9ac + K13ac + K15ac
D1	14.65014167	ab179008 - H2A K9me2
D2	12.83240938	ab179009 - H2A K9me3
D3	18.14742917	ab179010 - H2A R11me1
D4	15.05689375	ab179011 - H2A R11me2s
D5	15.27001771	ab179012 - H2A R11me2a
D6	17.20824792	ab179013 - H2A K13ac
D7	14.61275938	ab179014 - H2A K15ac
D8	13.97922188	ab179015 - H2A R17me1
D9	22.94463542	ab179016 - H2A R17me2s
D10	23.66669375	ab179017 - H2A R17me2a
D11	1678.697316	ab179018 - H2A unmod(20-50)
D12	167.9324438	ab179019 - H2A P26hyd
D13	40.80934792	ab179020 - H2A R29me1
D14	21.58300521	ab179021 - H2A R29me2s
D15	26.81659792	ab179022 - H2A R29me2a
D16	21.71382604	ab179023 - H2A K36ac
D17	17.93251667	ab179024 - H2A K36cr
D18	20.49853542	ab179025 - H2A Y39hyd
D19	22.57213333	ab179026 - H2A R42me1
D20	27.45506979	ab179027 - H2A R42me2s
D21	21.319525	ab179028 - H2A R42me2a
D22	17.16615833	ab179029 - H2A unmod(71-103)
D23	19.86455417	ab179030 - H2A R77me1
D24	29.496075	ab179031 - H2A R77me2a
D25	19.37990729	ab179032 - H2A R88me1
E1	21.63183438	ab179033 - H2A R88me2s
E2	15.48324688	ab179034 - H2A R88me2a
E3	21.34722188	ab179035 - H2A K95me1
E4	18.13448125	ab179036 - H2A K95me3
E5	19.41393854	ab179037 - H2A unmod(92-123)
E6	17.95551875	ab179038 - H2A T101GlcNAc
E7	17.14022188	ab179039 - H2A Q104me1
E8	15.42045625	ab179040 - H2A unmod(112-129)
E9	13.67662292	ab179041 - H2A K118me1
E10	11.31290104	ab179042 - H2A K118me2
E11	12.3083875	ab179043 - H2A K118me3
E12	19.67561146	ab179044 - H2A K118fo
E13	16.594875	ab179045 - H2A K118cr
E14	34.71939167	ab179046 - H2A K119cr
E15	19.53444167	ab179047 - H2A K118ubiq
E16	24.93713958	ab179048 - H2A K119ubiq
E17	29.69475208	ab179049 - H2A T120p
E18	21.00784896	ab179050 - H2A K125me1
E19	20.16045833	ab179051 - H2A K125me2
E20	19.69655	ab179052 - H2A K125cr
E21	9700.932507	ab179053 - H2A.X unmod(1-30)
E22	18.3830875	ab179054 - H2A.X K5ac
E23	5510.277839	ab179055 - H2A.X unmod(114-142)
E24	65.14174375	ab179056 - H2A.X K118ubiq
E25	26.92486354	ab179057 - H2A.X K119ubiq
F1	18.57047708	ab179058 - H2A.X T120p
F2	24.61268542	ab179059 - H2A.X K127ubiq
F3	33.59055208	ab179060 - H2A.X S139p + Y142p
F4	44.61505417	ab179061 - H2A.Z K4ac
F5	963.1713188	ab179062 - H2A.Z K7ac
F6	22.65119375	ab179063 - H2A.Z K11ac
F7	66.90206354	ab179064 - H2A.Z K4ac + K7ac + K11ac
F8	31.18192917	ab179065 - H2B unmod(1-30)
F9	18.44508021	ab179066 - H2B K5ac
F10	22.04320521	ab179067 - H2B K5cr
F11	21.20252813	ab179068 - H2B K5fo
F12	17.45343646	ab179069 - H2B P10hyd
F13	17.16373438	ab179070 - H2B K11me1
F14	21.93355833	ab179071 - H2B K11me2
F15	17.88971979	ab179072 - H2B K11me3
F16	22.2310125	ab179073 - H2B K11cr
F17	22.2401875	ab179074 - H2B K12ac
F18	20.05875104	ab179075 - H2B K12cr
F19	19.4500125	ab179076 - H2B K12me1

F20	21.66570833	ab179077 - H2B K12me2
F21	17.6441125	ab179078 - H2B K12me3
F22	15.67066354	ab179079 - H2B K12ac + K15ac
F23	20.44781042	ab179080 - H2B S14p
F24	14.98320521	ab179081 - H2B K15ac
F25	21.68759479	ab179082 - H2B K15cr
G1	14.87599479	ab179083 - H2B K15me1
G2	21.89492292	ab179084 - H2B K15me2
G3	15.43666875	ab179085 - H2B K15me3
G4	88.58109583	ab179086 - H2B K16ac
G5	20.76439375	ab179087 - H2B K16cr
G6	30.18949167	ab179088 - H2B K20ac
G7	22.79969063	ab179089 - H2B K20cr
G8	22.93352083	ab179090 - H2B K20me1
G9	20.97591146	ab179091 - H2B K20me2
G10	16.30543438	ab179092 - H2B K20me3
G11	22.72911563	ab179093 - H2B K23cr
G12	20.40527813	ab179094 - H2B K23me1
G13	18.32069792	ab179095 - H2B K23me3
G14	21.78557813	ab179096 - H2B K24ac
G15	24.77156458	ab179097 - H2B unmod(26-56)
G16	18.49678646	ab179098 - H2B K34cr
G17	17.22861979	ab179099 - H2B K34fo
G18	19.15594271	ab179100 - H2B Y37hyd
G19	21.24399583	ab179101 - H2B Y37p
G20	20.16503438	ab179102 - H2B K43me2
G21	14.44822188	ab179103 - H2B K43me3
G22	19.88642917	ab179104 - H2B K46fo
G23	23.30131042	ab179105 - H2B K46ac
G24	14.69356354	ab179106 - H2B K46me1
G25	23.11974167	ab179107 - H2B K46me2
H1	13.7644375	ab179108 - H2B K46me3
H2	23.58986771	ab179109 - H2B unmod(51-81)
H3	16.08195313	ab179110 - H2B K57me1
H4	20.21003021	ab179111 - H2B K57me2
H5	12.22254896	ab179112 - H2B K57me3
H6	58.52016354	ab179113 - H2B unmod(73-103)
H7	22.30412292	ab179114 - H2B R79me1
H8	25.48233229	ab179115 - H2B R79me2s
H9	23.804125	ab179116 - H2B R79me2a
H10	24.76345417	ab179117 - H2B Y83hyd
H11	26.23825521	ab179118 - H2B K85me1
H12	19.10093021	ab179119 - H2B K85me2
H13	21.03142604	ab179120 - H2B K85me3
H14	25.31198333	ab179121 - H2B unmod(94-125)
H15	23.44727396	ab179122 - H2B R99me1
H16	20.24659271	ab179123 - H2B R99me2s
H17	19.94977604	ab179124 - H2B R99me2a
H18	20.561925	ab179125 - H2B K108fo
H19	16.00215833	ab179126 - H2B K116fo
H20	20.62533542	ab179127 - H2B K116me1
H21	19.901925	ab179128 - H2B K116me2
H22	17.96467604	ab179129 - H2B K116me3
H23	21.40525938	ab179130 - H2B K116ubiq
H24	18.18166979	ab179131 - H2B K120ac
H25	19.89108333	ab179132 - H2B K120fo
I1	21.63581667	ab179133 - H2B K120ubiq
I2	27.5494375	ab179134 - H2B K125ubiq
I3	16.26115104	ab179135 - Sch2B unmod(102-132)
I4	12.21049375	ab179136 - H3 unmod(1-33)
I5	21.71149479	ab179137 - H3 R2cit
I6	11.95586771	ab179138 - H3 R2cit + K4me2
I7	21.94112708	ab179139 - H3 R2cit + K4me3
I8	17.92560313	ab179140 - H3 R2cit + K4me3 + K9ac + K14ac + K18ac
I9	20.69242708	ab179141 - H3 R2cit + K4ac + K9ac + K14ac + K18ac
I10	22.95676563	ab179142 - H3 R2me1
I11	15.74956667	ab179143 - H3 R2me1 + K4me2
I12	22.34333958	ab179144 - H3 R2me1 + K4me3
I13	14.66141458	ab179145 - H3 R2me1 + K4me3 + K9ac + K14ac + K18ac
I14	20.62123021	ab179146 - H3 R2me2s
I15	14.24036979	ab179147 - H3 R2me2s + K4me2
I16	20.68260521	ab179148 - H3 R2me2s + K4me3
I17	22.5797625	ab179149 - H3 R2me2s + K4me3 + K9ac + K14ac + K18ac
I18	22.18243646	ab179150 - H3 R2me2a

I19	24.23904063	ab179151 - H3 R2me2a + K4me2
I20	19.60400208	ab179152 - H3 R2me2a + K4me3
I21	19.32960313	ab179153 - H3 R2me2a + K4me3 + S10p
I22	19.63469167	ab179154 - H3 R2me2a + T3p + K4me3
I23	22.59705625	ab179155 - H3 R2me2a + K4ac + K9ac + K14ac + K18ac
I24	19.52028229	ab179156 - H3 R2me2a + K4me3 + K9ac + K14ac + K18ac
I25	21.23980313	ab179157 - H3 R2me2a + T3p + K4me3 + K9ac + K14ac + K18ac
J1	23.77865938	ab179158 - H3 R2me2a + K4me3 + K9ac + S10p + K14ac + K18ac
J2	21.88838021	ab179159 - H3 T3p
J3	24.08360729	ab179160 - H3 T3p + K4me1
J4	17.40138229	ab179161 - H3 T3p + K4me2
J5	21.63223958	ab179162 - H3 T3p + K4me3 + K9ac + K14ac + K18ac
J6	21.52135	ab179163 - H3 K4ac
J7	15.03416875	ab179164 - H3 K4ac + K9ac
J8	21.30426667	ab179165 - H3 K4ac + K14ac
J9	17.39169688	ab179166 - H3 K4ac + K18ac
J10	22.41149271	ab179167 - H3 K4ac + K9me3
J11	15.56444896	ab179168 - H3 K4ac + K9ac + K14ac
J12	23.10595833	ab179169 - H3 K4ac + K9ac + K18ac
J13	16.55671042	ab179170 - H3 K4ac + K14ac + K18ac
J14	24.03567604	ab179171 - H3 K4ac + K9ac + K14ac + K18ac
J15	22.62582813	ab179172 - H3 K4ac + K9me1 + K14ac + K18ac
J16	20.78339583	ab179173 - H3 K4ac + K9me2 + K14ac + K18ac
J17	24.02072917	ab179174 - H3 K4ac + K9me3 + K14ac + K18ac
J18	21.61254896	ab179175 - H3 K4ac + K9ac + S10p + K14ac + K18ac
J19	22.69331458	ab179176 - H3 K4cr
J20	21.16137708	ab179177 - H3 K4me1 + K9me2
J21	22.30119167	ab179178 - H3 K4me1 + K9ac + K18ac
J22	20.92328229	ab179179 - H3 K4me1 + K9ac + K14ac + K18ac
J23	21.92633333	ab179180 - H3 K4me2 + T6p
J24	17.97806667	ab179181 - H3 K4me2 + K9me2
J25	22.49289167	ab179182 - H3 K4me2 + K9ac + K18ac
K1	23.4750125	ab179183 - H3 K4me2 + K9ac + K14ac + K18ac
K2	20.83341146	ab179184 - H3 K4me3 + T6p
K3	13.66670521	ab179185 - H3 K4me3 + K9ac
K4	23.04695208	ab179186 - H3 K4me3 + K9me2
K5	15.38309167	ab179187 - H3 K4me3 + K9me3
K6	20.22227604	ab179188 - H3 K4me3 + S10p
K7	11.78175938	ab179189 - H3 K4me3 + K14ac
K8	22.10609583	ab179190 - H3 K4me3 + K18ac
K9	13.23211875	ab179191 - H3 K4me3 + R8me2s + K9me3
K10	19.93580104	ab179192 - H3 K4me3 + R8me2a + K9me3
K11	23.66523333	ab179193 - H3 K4me3 + K9ac + S10p
K12	21.95130833	ab179194 - H3 K4me3 + K9ac + K14ac
K13	20.18547083	ab179195 - H3 K4me3 + K9ac + K18ac
K14	21.5922125	ab179196 - H3 K4me3 + K14ac + K18ac
K15	20.10439063	ab179197 - H3 K4me3 + K9ac + K14ac + K18ac
K16	23.93625208	ab179198 - H3 K4me3 + K9ac + S10p + K14ac + K18ac
K17	21.26623646	ab179199 - H3 K4me3 + T6p + K9ac + K14ac + K18ac
K18	23.06343021	ab179200 - H3 K4me2 + T6p + K9ac + K14ac + K18ac
K19	25.10415313	ab179201 - H3 K4bio + K9bio + K18bio
K20	18.59623646	ab179202 - H3 T6p + K9me3
K21	24.13093646	ab179203 - H3 T6p + R8me2a + K9me3
K22	16.78120729	ab179204 - H3 T6p + K9ac + K14ac + K18ac
K23	24.84570313	ab179205 - H3 R8me1
K24	16.08210625	ab179206 - H3 R8me1 + K9me1
K25	17.43438646	ab179207 - H3 R8me1 + K9me2
L1	15.34978021	ab179208 - H3 R8me1 + K9me3
L2	22.7602625	ab179209 - H3 R8me2s
L3	18.62063125	ab179210 - H3 R8me2s + K9me1
L4	18.01485104	ab179211 - H3 R8me2s + K9me2
L5	15.68254688	ab179212 - H3 R8me2s + K9me3
L6	20.68341875	ab179213 - H3 R8me2a
L7	19.21414688	ab179214 - H3 R8me2a + K9me1
L8	20.5304875	ab179215 - H3 R8me2a + K9me2
L9	17.85310833	ab179216 - H3 R8me2a + K9me3
L10	21.57136979	ab179217 - H3 R8cit
L11	18.39131042	ab179218 - H3 K9ac + K14ac
L12	27.15737396	ab179219 - H3 K9ac + K18ac
L13	21.19519896	ab179220 - H3 K9ac + K14ac + K18ac
L14	23.46261563	ab179221 - H3 K9ac + K14ac + K18ac + K23ac + K27ac
L15	19.43644063	ab179222 - H3 K9ac + S10p
L16	22.92394479	ab179223 - H3 K9cr
L17	25.74629792	ab179224 - H3 K9me2

L18	21.94281979	ab179225 - H3 K9me2 + S10p
L19	16.62948854	ab179226 - H3 K9me2 + K27me2
L20	16.84029688	ab179227 - H3 S10p + S28p
L21	21.86462604	ab179228 - H3 S10p K14ac
L22	24.20408333	ab179229 - H3 K14me1
L23	21.34772083	ab179230 - H3 K14me2
L24	21.77162188	ab179231 - H3 K14me3
L25	23.4500625	ab179232 - H3 K14ac + K18ac
M1	23.78845313	ab179233 - H3 R17cit
M2	21.10541042	ab179234 - H3 R17me1
M3	20.31862813	ab179235 - H3 R17me2s
M4	20.33910625	ab179236 - H3 R17me2a + K18ac
M5	20.42442292	ab179237 - H3 K18cr
M6	21.53888021	ab179238 - H3 K18me1
M7	22.8217875	ab179239 - H3 K18me2
M8	23.98682396	ab179240 - H3 K18me3 + K36me3
M9	18.66393333	ab179241 - H3 K23cr
M10	21.38115938	ab179242 - H3 K23me1
M11	16.59094896	ab179243 - H3 K23me2
M12	23.79643542	ab179244 - H3 K23me3
M13	17.67318229	ab179245 - H3 T22p + K23me3
M14	22.38725417	ab179246 - H3 K23pr
M15	25.01815104	ab179247 - H3 unmod(20-51)
M16	24.06726042	ab179248 - H3 R26me1
M17	24.32213542	ab179249 - H3 R26me2a
M18	22.55473438	ab179250 - H3 R26me2a + K27me1
M19	23.45807396	ab179251 - H3 R26me2a + K27me3
M20	17.85703125	ab179252 - H3 K27ac + S28p
M21	21.28173333	ab179253 - H3 K27cr
M22	19.86440104	ab179254 - H3 K27me1
M23	16.87241042	ab179255 - H3 K27me2 + S28p
M24	23.54940313	ab179256 - H3 K27me3
M25	16.34602292	ab179257 - H3 K27me3 + S28p
N1	17.73480625	ab179258 - H3 P30hpro
N2	23.87812396	ab179259 - H3 P30(5,5me2)
N3	15.42591875	ab179260 - H3 P30
N4	20.86145208	ab179261 - H3 P30mutA
N5	19.08029896	ab179262 - H3 T32GlcNAc
N6	17.83803125	ab179263 - H3 K36ac
N7	21.51929479	ab179264 - H3 P38hpro
N8	22.98844271	ab179265 - H3 P38(5,5me2)
N9	18.16458854	ab179266 - H3 P38
N10	21.64345208	ab179267 - H3 P38mutA
N11	17.58194375	ab179268 - H3 Y41p
N12	23.07804792	ab179269 - H3 unmod(50-80)
N13	20.43357188	ab179270 - H3 K56ac
N14	20.60759271	ab179271 - H3 K56cr
N15	16.72868542	ab179272 - H3 K56fo
N16	20.27511563	ab179273 - H3 K56me1
N17	22.74347083	ab179274 - H3 K56me2
N18	20.89127188	ab179275 - H3 K56me3
N19	20.95127708	ab179276 - H3 R63me1
N20	21.43346563	ab179277 - H3 R63me2s
N21	20.31160208	ab179278 - H3 R63me2a
N22	17.51502813	ab179279 - H3 K64ac
N23	21.46669583	ab179280 - H3 K64me2
N24	25.94469375	ab179281 - H3 K64me3
N25	21.62182917	ab179282 - H3 unmod(73-103)
O1	25.63773021	ab179283 - H3 K79ac
O2	20.58256979	ab179284 - H3 K79fo
O3	20.35145833	ab179285 - H3 T80p
O4	23.14337292	ab179286 - H3 K79me3 + T80p
O5	22.36840521	ab179287 - H3 unmod(99-135)
O6	20.93843438	ab179288 - H3 T118O-ac
O7	20.5249125	ab179289 - H3 K122fo
O8	20.74728229	ab179290 - H3 K122me1
O9	16.43781875	ab179291 - H3 R128me1
O10	24.4928	ab179292 - H3 R128me2s
O11	25.44083854	ab179293 - H3.3 unmod(10-40)
O12	17.27880938	ab179294 - H3.3 S28p
O13	23.54630208	ab179295 - H3.3 S28p + S31p
O14	20.05665625	ab179296 - H3.3 S31p
O15	27.698325	ab179297 - H4 unmod(1-30)
O16	17.09020208	ab179298 - H4 S1p + R3me1

O17	19.71231458	ab179299 - H4 S1p + R3me2s
O18	21.00000208	ab179300 - H4 S1p + R3me2a
O19	21.10940833	ab179301 - H4 S1p + K5ac + K8ac + K12ac + K16ac
O20	22.94110208	ab179302 - H4 R3cit
O21	22.37950104	ab179303 - H4 R3me1 + K5ac + K8ac + K12ac + K16ac + K20ac
O22	23.98458438	ab179304 - H4 R3me2s
O23	24.34690625	ab179305 - H4 R3me2s + K5ac + K8ac + K12ac + K16ac + K20ac
O24	22.67284271	ab179306 - H4 R3me2a
O25	17.17799375	ab179307 - H4 R3me2a + K5ac
P1	25.80795521	ab179308 - H4 R3me2a + K5ac + K8ac + K12ac + K16ac + K20ac
P2	19.53279271	ab179309 - H4 K5ac
P3	18.55567188	ab179310 - H4 K5ac + K8ac
P4	21.58321979	ab179311 - H4 K5ac + K12ac
P5	23.35828646	ab179312 - H4 K5ac + K16ac
P6	20.99471979	ab179313 - H4 K5ac + K8ac + K12ac
P7	37.09807813	ab179314 - H4 K5ac + K8me1 + K12ac
P8	22.22771875	ab179315 - H4 K5ac + K8ac + K12ac + K16ac
P9	23.34018958	ab179316 - H4 K5ac + K8me1 + K12ac + K16ac
P10	22.92100521	ab179317 - H4 K5ac + K8ac + K12me1 + K16ac
P11	25.71188542	ab179318 - H4 K5me1
P12	18.20826458	ab179319 - H4 K5me1 + K8ac + K12me1
P13	22.73557604	ab179320 - H4 K5me1 + K8me1 + K12me1
P14	18.70150208	ab179321 - H4 K5me1 + K8ac + K12ac + K16ac
P15	22.37864375	ab179322 - H4 K5me2
P16	17.17638021	ab179323 - H4 K5me3
P17	26.86973958	ab179324 - H4 K5cr
P18	19.01409896	ab179325 - H4 K8ac + K12ac
P19	21.04330521	ab179326 - H4 K8ac + K16ac
P20	20.28694688	ab179327 - H4 K8cr
P21	19.960625	ab179328 - H4 K8me1
P22	16.42760313	ab179329 - H4 K12ac
P23	20.23080729	ab179330 - H4 K12ac + K16ac
P24	15.70531042	ab179331 - H4 K12ac + K16ac + K20me2
P25	19.69405625	ab179332 - H4 K12ac + K16ac + K20me3
Q1	25.39811563	ab179333 - H4 K5mutQ + K8mutQ + K12ac + K16mutQ + K20mutQ
Q2	21.35831875	ab179334 - H4 K12cr
Q3	23.65003229	ab179335 - H4 K12fo
Q4	17.69408333	ab179336 - H4 K12me1
Q5	24.20096458	ab179337 - H4 K12me2
Q6	17.64537396	ab179338 - H4 K12me3
Q7	24.28805938	ab179339 - H4 K16ac
Q8	18.59897396	ab179340 - H4 K16ac + K20ac
Q9	19.84249271	ab179341 - H4 K16cr
Q10	22.20268854	ab179342 - H4 K16me1
Q11	21.68894688	ab179343 - H4 K16me3
Q12	19.899725	ab179344 - H4 R17me1
Q13	23.27349375	ab179345 - H4 R17me2s
Q14	22.50801771	ab179346 - H4 R17me2a
Q15	22.93510208	ab179347 - H4 R19me1
Q16	15.96533438	ab179348 - H4 R19me2s
Q17	23.8179125	ab179349 - H4 R19me2a
Q18	27.50577188	ab179350 - H4 K20ac
Q19	17.87926146	ab179351 - H4 R23me1
Q20	20.56637708	ab179352 - H4 R23me2s
Q21	21.20298854	ab179353 - H4 R23me2a
Q22	25.48605417	ab179354 - H4 unmod(24-54)
Q23	19.87522917	ab179355 - H4 K31fo
Q24	22.09689583	ab179356 - H4 K31me1
Q25	22.07509688	ab179357 - H4 K31me2
R1	19.37432813	ab179358 - H4 K31me3
R2	24.99150729	ab179359 - H4 P32hyd
R3	18.48465417	ab179360 - H4 R35me1
R4	24.62415625	ab179361 - H4 R35me2s
R5	19.76418958	ab179362 - H4 R35me2a
R6	23.57124583	ab179363 - H4 S47O-ac
R7	20.40471979	ab179364 - H4 S47GlcNAc
R8	25.18546771	ab179365 - H4 S47p
R9	20.00749167	ab179366 - H4 unmod(45-75)
R10	23.1349625	ab179367 - H4 Y51hyd
R11	21.19698438	ab179368 - H4 R55me1
R12	20.83674063	ab179369 - H4 R55me2s
R13	17.06951979	ab179370 - H4 R55me2a
R14	20.86822083	ab179371 - H4 K59fo
R15	19.16418229	ab179372 - H4 K59me1

R16	20.24960729	ab179373 - H4 K59me2
R17	18.08558333	ab179374 - H4 R67me1
R18	20.50170833	ab179375 - H4 R67me2s
R19	19.23172083	ab179376 - H4 R67me2a
R20	23.43757292	ab179377 - H4 unmod(71-102)
R21	19.97121771	ab179378 - H4 K77me1
R22	22.24527083	ab179379 - H4 K77me2
R23	23.89836563	ab179380 - H4 K77me3
R24	23.22982083	ab179381 - H4 K79fo
R25	23.18687396	ab179382 - H4 K79me1
S1	21.12198854	ab179383 - H4 K79me3
S2	20.99866146	ab179384 - H4 T82O-ac
S3	25.92142083	ab179385 - H4 Y88hyd
S4	17.6643375	ab179386 - H4 Y88nitro
S5	19.9966375	ab179387 - H4 K91fo
S6	19.34257708	ab179403 - H2A K5ac
S7	26.07656667	ab179404 - H2A K9me1
S8	22.29798125	ab18504 - H2B S32p
S9	24.83296667	ab19828 - ScH2A S129p
S10	20.67409063	ab20631 - H3 R26cit
S11	19.54740313	ab20635 - H3 T3p + K4me3
S12	23.050875	ab21043 - H3 K37me1
S13	17.94925104	ab21998 - H1.4 K25me2
S14	20.58927188	ab22079 - H4 K59me3
S15	17.96213958	ab22080 - H3 K122me2
S16	24.18852813	ab22082 - H3 K122me3
S17	24.96752396	ab22214 - H3 K37me2
S18	24.77961146	ab22398 - H2A R3me2a
S19	43.80041146	ab22399 - H2A R3me2s
S20	20.42065417	ab23539 - H3 T22p
S21	18.19386458	ab24003 - H3 K18ac
S22	21.70531771	ab24404 - H3 K27ac
S23	21.84129375	ab24417 - H3 K37me3
S24	16.67240417	ab24444 - H3 T11p
S25	67.06116354	ab24659 - H2A K127me3
T1	21.99744896	ab24660 - H2A K125me3
T2	18.91555521	ab26307 - H3 T45p
T3	18.79984688	ab26342 - H2B K5me2
T4	20.86020521	ab26739 - H4 T30p
T5	23.80985417	ab27896 - H4 T96p
T6	21.99999896	ab27898 - H4 T80p
T7	20.44760833	ab28756 - H2A R3me1
T8	15.12164271	ab32876 - H3 R2cit + R8cit + R17cit
T9	20.75006563	ab33047 - H2A K127me2
T10	23.49098542	ab33505 - H3 T118p
T11	20.79751979	ab33594 - H3 R128me2a
T12	25.76139479	ab34466 - H3 K122ac
T13	20.13029167	ab36927 - H2A S122p
T14	26.9520125	ab4555 - H3 K79me1
T15	21.39550208	ab4556 - H3 K79me2
T16	24.45127708	ab4557 - H3 K79me3
T17	21.68890729	ab4560 - H4 K79me2
T18	21.13693021	ab46854 - H3 K18me3
T19	23.72018542	ab48359 - H3 K23ac
T20	22.68086458	ab53244 - H3 K64me1
T21	23.95346771	ab54016 - H2A K9ac
T22	22.122375	ab5499 - H3 S28p
T23	24.81909479	ab7768 - H3 K4me2
T24	22.4583125	ab89018 - H1.4 S26p
T25	20.09603021	ab92374 - H3 K4me3 (HEX1)
U1	18.86790625	ab95444 - H4 K16me2