

Peptide array: gamma H2A.X(phospho S139) (ab303656)

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

Cell Ref	AUC	Peptide
A1	2.403164583	ab10112 - H4 K91ac
A2	2.906578125	ab10121 - H2B K85ac
A3	1.3734625	ab10127 - H2A K95me2
A4	2.5791875	ab10138 - H1.4 T17p
A5	1.937275	ab10139 - H1.4 T146p
A6	14.65022292	ab113997 - H2A.X Y142p
A7	1.781889583	ab11477 - H3 S10p
A8	2.073698958	ab12946 - H2B K5me3
A9	2.030945833	ab12949 - H2B K23me2
A10	1.178484375	ab12952 - H2A K127me1
A11	1.861811458	ab13211 - H2B K5me1
A12	1.871772917	ab1340 - H3 K4me1
A13	2.217269792	ab13570 - H2B K43me1
A14	2.456889583	ab13835 - HA
A15	1.947769792	ab13837 - Myc
A16	2.1946875	ab140408 - ScH2B T129p
A17	1.712453125	ab14103 - H3 T6p
A18	1.899857292	ab14724 - H4 S1p
A19	1.637078125	ab14799 - H3 T32p
A20	2.316101042	ab14943 - His
A21	2.160414583	ab14952 - H2A R77me2s
A22	1.681153125	ab14964 - H4 K20me2
A23	2.664503125	ab14997 - ScH2A K21ac
A24	1.898346875	ab154782 - H3 R26me2s
A25	2.902771875	ab15591 - H3 K14ac
B1	1.527729167	ab15644 - H3 K9me3 + S10p
B2	104.7812948	ab15645 - H2A.X S139p
B3	2.138653125	ab15824 - H4 K8ac
B4	1.959954167	ab15829 - V5
B5	2.145694792	ab16635 - H3 K9ac
B6	1.851185417	ab166685 - H2B S112GlcNAc
B7	2.91523125	ab166688 - H2B S36GlcNAc
B8	3.30663125	ab16935 - H3 R17me2a
B9	1.952952083	ab17043 - H4 K20me1
B10	2.790594792	ab17162 - H3 P16hyd + P30hyd
B11	2.258472917	ab17567 - H4 K20me3
B12	2.305141667	ab17587 - H1.4 K25me3
B13	2.477985417	ab175905 - H2B K11ac
B14	2.452147917	ab176211 - H2B K23ac
B15	1.945626042	ab17632 - H4 K5bio
B16	2.271376042	ab1771 - H3 K9me1
B17	2.374486458	ab1773 - H3 K9me3
B18	2.107970833	ab17770 - H4 R3me1
B19	2.410909375	ab1781 - H3 K27me2
B20	2.380051042	ab1783 - H3 K36me1
B21	2.178013542	ab1784 - H3 K36me2
B22	2.10194375	ab1785 - H3 K36me3
B23	2.117236458	ab178667 - ScH2A unmod(1-30)
B24	1.957471875	ab178668 - ScH2A unmod(102-132)
B25	1.48671875	ab178670 - H2A.Z unmod(1-30)
C1	2.814160417	ab178833 - Flag
C2	1.742504167	ab178835 - H2B R99mutO
C3	1.598966667	ab178836 - H4 R23mutO
C4	2.0719125	ab178837 - H4 G48mutD
C5	1.5600875	ab178987 - H1.4 unmod(1-30)
C6	2.766360417	ab178988 - H1.4 S1p
C7	1.528721875	ab178989 - H1.4 T3p
C8	3.136391667	ab178990 - H1.4 S1p + T3p
C9	2.178185417	ab178991 - H1.4 K25me1
C10	2.003173958	ab178992 - H1.4 K25me3 + S26p
C11	1.758847917	ab178993 - H1.4 unmod(130-160)
C12	1.3242625	ab178994 - H2A unmod(1-30)
C13	2.299622917	ab178995 - H2A S1p
C14	2.195591667	ab178996 - H2A S1p + K5ac
C15	2.562684375	ab178997 - H2A S1p + R3me2a + K5ac
C16	2.075126042	ab178998 - H2A S1p + R3me2s + K5ac
C17	2.028239583	ab178999 - H2A S1p + R3cit + K5ac
C18	1.363890625	ab179000 - H2A S1ac + S1p + K5ac + K9ac + K13ac + K15ac
C19	2.003102083	ab179001 - H2A S1p + K5ac + K9ac + K13ac + K15ac
C20	2.17503125	ab179002 - H2A R3cit

C21	2.267752083	ab179003 - H2A R3cit + K5ac
C22	1.306807292	ab179004 - H2A R3me2a + K5ac
C23	1.688830208	ab179005 - H2A R3me2s + K5ac
C24	1.881139583	ab179006 - H2A S1ac + K5ac + K9ac + K13ac + K15ac
C25	3.704830208	ab179007 - H2A K5ac + K9ac + K13ac + K15ac
D1	1.628290625	ab179008 - H2A K9me2
D2	2.109442708	ab179009 - H2A K9me3
D3	1.454603125	ab179010 - H2A R11me1
D4	2.464865625	ab179011 - H2A R11me2s
D5	1.490236458	ab179012 - H2A R11me2a
D6	2.917553125	ab179013 - H2A K13ac
D7	1.593739583	ab179014 - H2A K15ac
D8	3.173255208	ab179015 - H2A R17me1
D9	1.529340625	ab179016 - H2A R17me2s
D10	2.467763542	ab179017 - H2A R17me2a
D11	2.310582292	ab179018 - H2A unmod(20-50)
D12	1.821273958	ab179019 - H2A P26hyd
D13	2.021602083	ab179020 - H2A R29me1
D14	1.740622917	ab179021 - H2A R29me2s
D15	1.859952083	ab179022 - H2A R29me2a
D16	2.130877083	ab179023 - H2A K36ac
D17	1.583730208	ab179024 - H2A K36cr
D18	2.205763542	ab179025 - H2A Y39hyd
D19	1.768223958	ab179026 - H2A R42me1
D20	1.772434375	ab179027 - H2A R42me2s
D21	2.338440625	ab179028 - H2A R42me2a
D22	2.17009375	ab179029 - H2A unmod(71-103)
D23	2.022034375	ab179030 - H2A R77me1
D24	1.852945833	ab179031 - H2A R77me2a
D25	2.329592708	ab179032 - H2A R88me1
E1	2.27014375	ab179033 - H2A R88me2s
E2	1.026796875	ab179034 - H2A R88me2a
E3	1.392096875	ab179035 - H2A K95me1
E4	1.524364583	ab179036 - H2A K95me3
E5	0.992102083	ab179037 - H2A unmod(92-123)
E6	2.461019792	ab179038 - H2A T101GlcNAc
E7	2.896013542	ab179039 - H2A Q104me1
E8	1.505802083	ab179040 - H2A unmod(112-129)
E9	3.115054167	ab179041 - H2A K118me1
E10	1.301053125	ab179042 - H2A K118me2
E11	2.700267708	ab179043 - H2A K118me3
E12	2.125173958	ab179044 - H2A K118fo
E13	2.996369792	ab179045 - H2A K118cr
E14	1.750469792	ab179046 - H2A K119cr
E15	1.822661458	ab179047 - H2A K118ubiq
E16	1.724023958	ab179048 - H2A K119ubiq
E17	2.018959375	ab179049 - H2A T120p
E18	2.160752083	ab179050 - H2A K125me1
E19	2.070203125	ab179051 - H2A K125me2
E20	1.729209375	ab179052 - H2A K125cr
E21	1.92379375	ab179053 - H2A.X unmod(1-30)
E22	1.90849375	ab179054 - H2A.X K5ac
E23	1.567016667	ab179055 - H2A.X unmod(114-142)
E24	2.055163542	ab179056 - H2A.X K118ubiq
E25	2.121211458	ab179057 - H2A.X K119ubiq
F1	1.883432292	ab179058 - H2A.X T120p
F2	1.379889583	ab179059 - H2A.X K127ubiq
F3	13.78661875	ab179060 - H2A.X S139p + Y142p
F4	2.522244792	ab179061 - H2A.Z K4ac
F5	1.87289375	ab179062 - H2A.Z K7ac
F6	2.471917708	ab179063 - H2A.Z K11ac
F7	2.068907292	ab179064 - H2A.Z K4ac + K7ac + K11ac
F8	1.977113542	ab179065 - H2B unmod(1-30)
F9	2.63453125	ab179066 - H2B K5ac
F10	1.964869792	ab179067 - H2B K5cr
F11	1.830304167	ab179068 - H2B K5fo
F12	1.679247917	ab179069 - H2B P10hyd
F13	2.711084375	ab179070 - H2B K11me1
F14	2.237632292	ab179071 - H2B K11me2
F15	3.0525625	ab179072 - H2B K11me3
F16	2.835332292	ab179073 - H2B K11cr
F17	3.028458333	ab179074 - H2B K12ac
F18	2.652407292	ab179075 - H2B K12cr
F19	3.046013542	ab179076 - H2B K12me1

F20	3.938242708	ab179077 - H2B K12me2
F21	1.850733333	ab179078 - H2B K12me3
F22	1.863848958	ab179079 - H2B K12ac + K15ac
F23	1.108010417	ab179080 - H2B S14p
F24	1.630382292	ab179081 - H2B K15ac
F25	1.36756875	ab179082 - H2B K15cr
G1	1.950345833	ab179083 - H2B K15me1
G2	2.240604167	ab179084 - H2B K15me2
G3	2.607657292	ab179085 - H2B K15me3
G4	1.708448958	ab179086 - H2B K16ac
G5	1.929755208	ab179087 - H2B K16cr
G6	1.85174375	ab179088 - H2B K20ac
G7	1.84184375	ab179089 - H2B K20cr
G8	1.637073958	ab179090 - H2B K20me1
G9	2.04905625	ab179091 - H2B K20me2
G10	1.718630208	ab179092 - H2B K20me3
G11	2.4442875	ab179093 - H2B K23cr
G12	2.184103125	ab179094 - H2B K23me1
G13	2.238514583	ab179095 - H2B K23me3
G14	2.222782292	ab179096 - H2B K24ac
G15	2.642334375	ab179097 - H2B unmod(26-56)
G16	2.267525	ab179098 - H2B K34cr
G17	3.303782292	ab179099 - H2B K34fo
G18	3.630766667	ab179100 - H2B Y37hyd
G19	2.646391667	ab179101 - H2B Y37p
G20	2.991132292	ab179102 - H2B K43me2
G21	2.670017708	ab179103 - H2B K43me3
G22	2.554025	ab179104 - H2B K46fo
G23	1.433876042	ab179105 - H2B K46ac
G24	2.567022917	ab179106 - H2B K46me1
G25	0.8424125	ab179107 - H2B K46me2
H1	2.538194792	ab179108 - H2B K46me3
H2	2.525444792	ab179109 - H2B unmod(51-81)
H3	2.350659375	ab179110 - H2B K57me1
H4	1.667719792	ab179111 - H2B K57me2
H5	2.508954167	ab179112 - H2B K57me3
H6	1.667965625	ab179113 - H2B unmod(73-103)
H7	2.353669792	ab179114 - H2B R79me1
H8	1.282667708	ab179115 - H2B R79me2s
H9	1.976401042	ab179116 - H2B R79me2a
H10	1.966815625	ab179117 - H2B Y83hyd
H11	2.049911458	ab179118 - H2B K85me1
H12	2.197738542	ab179119 - H2B K85me2
H13	1.69878125	ab179120 - H2B K85me3
H14	1.755786458	ab179121 - H2B unmod(94-125)
H15	1.818128125	ab179122 - H2B R99me1
H16	2.004397917	ab179123 - H2B R99me2s
H17	1.626975	ab179124 - H2B R99me2a
H18	2.106639583	ab179125 - H2B K108fo
H19	3.182697917	ab179126 - H2B K116fo
H20	2.462157292	ab179127 - H2B K116me1
H21	3.624053125	ab179128 - H2B K116me2
H22	3.011992708	ab179129 - H2B K116me3
H23	3.450611458	ab179130 - H2B K116ubiq
H24	2.231819792	ab179131 - H2B K120ac
H25	3.252869792	ab179132 - H2B K120fo
I1	2.522946875	ab179133 - H2B K120ubiq
I2	1.596797917	ab179134 - H2B K125ubiq
I3	2.108241667	ab179135 - Sch2B unmod(102-132)
I4	1.93123125	ab179136 - H3 unmod(1-33)
I5	1.447485417	ab179137 - H3 R2cit
I6	2.875521875	ab179138 - H3 R2cit + K4me2
I7	1.606480208	ab179139 - H3 R2cit + K4me3
I8	3.138116667	ab179140 - H3 R2cit + K4me3 + K9ac + K14ac + K18ac
I9	1.553128125	ab179141 - H3 R2cit + K4ac + K9ac + K14ac + K18ac
I10	1.750621875	ab179142 - H3 R2me1
I11	2.620927083	ab179143 - H3 R2me1 + K4me2
I12	2.005688542	ab179144 - H3 R2me1 + K4me3
I13	2.924589583	ab179145 - H3 R2me1 + K4me3 + K9ac + K14ac + K18ac
I14	1.804439583	ab179146 - H3 R2me2s
I15	1.698913542	ab179147 - H3 R2me2s + K4me2
I16	2.241855208	ab179148 - H3 R2me2s + K4me3
I17	2.501279167	ab179149 - H3 R2me2s + K4me3 + K9ac + K14ac + K18ac
I18	2.242852083	ab179150 - H3 R2me2a

I19	2.265261458	ab179151 - H3 R2me2a + K4me2
I20	1.332121875	ab179152 - H3 R2me2a + K4me3
I21	1.804517708	ab179153 - H3 R2me2a + K4me3 + S10p
I22	1.396353125	ab179154 - H3 R2me2a + T3p + K4me3
I23	1.888884375	ab179155 - H3 R2me2a + K4ac + K9ac + K14ac + K18ac
I24	1.597538542	ab179156 - H3 R2me2a + K4me3 + K9ac + K14ac + K18ac
I25	2.082433333	ab179157 - H3 R2me2a + T3p + K4me3 + K9ac + K14ac + K18ac
J1	1.631139583	ab179158 - H3 R2me2a + K4me3 + K9ac + S10p + K14ac + K18ac
J2	2.22005	ab179159 - H3 T3p
J3	1.208922917	ab179160 - H3 T3p + K4me1
J4	1.8065	ab179161 - H3 T3p + K4me2
J5	2.067755208	ab179162 - H3 T3p + K4me3 + K9ac + K14ac + K18ac
J6	2.002858333	ab179163 - H3 K4ac
J7	1.71160625	ab179164 - H3 K4ac + K9ac
J8	1.604142708	ab179165 - H3 K4ac + K14ac
J9	2.156588542	ab179166 - H3 K4ac + K18ac
J10	1.964809375	ab179167 - H3 K4ac + K9me3
J11	1.928283333	ab179168 - H3 K4ac + K9ac + K14ac
J12	2.137333333	ab179169 - H3 K4ac + K9ac + K18ac
J13	2.783835417	ab179170 - H3 K4ac + K14ac + K18ac
J14	1.969457292	ab179171 - H3 K4ac + K9ac + K14ac + K18ac
J15	2.354672917	ab179172 - H3 K4ac + K9me1 + K14ac + K18ac
J16	1.837146875	ab179173 - H3 K4ac + K9me2 + K14ac + K18ac
J17	11.99686354	ab179174 - H3 K4ac + K9me3 + K14ac + K18ac
J18	2.273371875	ab179175 - H3 K4ac + K9ac + S10p + K14ac + K18ac
J19	2.446909375	ab179176 - H3 K4cr
J20	2.007455208	ab179177 - H3 K4me1 + K9me2
J21	1.906513542	ab179178 - H3 K4me1 + K9ac + K18ac
J22	2.294948958	ab179179 - H3 K4me1 + K9ac + K14ac + K18ac
J23	1.526229167	ab179180 - H3 K4me2 + T6p
J24	1.776448958	ab179181 - H3 K4me2 + K9me2
J25	2.245792708	ab179182 - H3 K4me2 + K9ac + K18ac
K1	2.036488542	ab179183 - H3 K4me2 + K9ac + K14ac + K18ac
K2	1.525707292	ab179184 - H3 K4me3 + T6p
K3	1.96545625	ab179185 - H3 K4me3 + K9ac
K4	2.25844375	ab179186 - H3 K4me3 + K9me2
K5	2.562815625	ab179187 - H3 K4me3 + K9me3
K6	1.916535417	ab179188 - H3 K4me3 + S10p
K7	1.861680208	ab179189 - H3 K4me3 + K14ac
K8	2.356184375	ab179190 - H3 K4me3 + K18ac
K9	2.040602083	ab179191 - H3 K4me3 + R8me2s + K9me3
K10	2.51933125	ab179192 - H3 K4me3 + R8me2a + K9me3
K11	2.10088125	ab179193 - H3 K4me3 + K9ac + S10p
K12	1.622215625	ab179194 - H3 K4me3 + K9ac + K14ac
K13	2.56449375	ab179195 - H3 K4me3 + K9ac + K18ac
K14	1.897985417	ab179196 - H3 K4me3 + K14ac + K18ac
K15	2.066914583	ab179197 - H3 K4me3 + K9ac + K14ac + K18ac
K16	1.584588542	ab179198 - H3 K4me3 + K9ac + S10p + K14ac + K18ac
K17	2.46168125	ab179199 - H3 K4me3 + T6p + K9ac + K14ac + K18ac
K18	1.328533333	ab179200 - H3 K4me2 + T6p + K9ac + K14ac + K18ac
K19	2.090315625	ab179201 - H3 K4bio + K9bio + K18bio
K20	1.685476042	ab179202 - H3 T6p + K9me3
K21	1.490117708	ab179203 - H3 T6p + R8me2a + K9me3
K22	1.892451042	ab179204 - H3 T6p + K9ac + K14ac + K18ac
K23	1.741209375	ab179205 - H3 R8me1
K24	1.777290625	ab179206 - H3 R8me1 + K9me1
K25	2.11359375	ab179207 - H3 R8me1 + K9me2
L1	2.399778125	ab179208 - H3 R8me1 + K9me3
L2	1.91269375	ab179209 - H3 R8me2s
L3	2.610347917	ab179210 - H3 R8me2s + K9me1
L4	2.043484375	ab179211 - H3 R8me2s + K9me2
L5	3.74250625	ab179212 - H3 R8me2s + K9me3
L6	2.067532292	ab179213 - H3 R8me2a
L7	1.986371875	ab179214 - H3 R8me2a + K9me1
L8	1.956938542	ab179215 - H3 R8me2a + K9me2
L9	2.186445833	ab179216 - H3 R8me2a + K9me3
L10	1.901090625	ab179217 - H3 R8cit
L11	2.46781875	ab179218 - H3 K9ac + K14ac
L12	2.183316667	ab179219 - H3 K9ac + K18ac
L13	1.923915625	ab179220 - H3 K9ac + K14ac + K18ac
L14	2.023344792	ab179221 - H3 K9ac + K14ac + K18ac + K23ac + K27ac
L15	2.227673958	ab179222 - H3 K9ac + S10p
L16	2.743310417	ab179223 - H3 K9cr
L17	2.590576042	ab179224 - H3 K9me2

L18	1.717090625	ab179225 - H3 K9me2 + S10p
L19	2.051794792	ab179226 - H3 K9me2 + K27me2
L20	1.789608333	ab179227 - H3 S10p + S28p
L21	2.124834375	ab179228 - H3 S10p K14ac
L22	2.356601042	ab179229 - H3 K14me1
L23	2.3992125	ab179230 - H3 K14me2
L24	2.025151042	ab179231 - H3 K14me3
L25	2.166760417	ab179232 - H3 K14ac + K18ac
M1	2.423167708	ab179233 - H3 R17cit
M2	2.865645833	ab179234 - H3 R17me1
M3	2.700271875	ab179235 - H3 R17me2s
M4	3.662708333	ab179236 - H3 R17me2a + K18ac
M5	3.97533125	ab179237 - H3 K18cr
M6	3.209196875	ab179238 - H3 K18me1
M7	1.716914583	ab179239 - H3 K18me2
M8	2.054385417	ab179240 - H3 K18me3 + K36me3
M9	1.386636458	ab179241 - H3 K23cr
M10	1.529657292	ab179242 - H3 K23me1
M11	2.057630208	ab179243 - H3 K23me2
M12	1.926017708	ab179244 - H3 K23me3
M13	2.3304875	ab179245 - H3 T22p + K23me3
M14	2.422315625	ab179246 - H3 K23pr
M15	2.759355208	ab179247 - H3 unmod(20-51)
M16	2.15891875	ab179248 - H3 R26me1
M17	3.154834375	ab179249 - H3 R26me2a
M18	2.397973958	ab179250 - H3 R26me2a + K27me1
M19	3.135010417	ab179251 - H3 R26me2a + K27me3
M20	3.584288542	ab179252 - H3 K27ac + S28p
M21	1.496471875	ab179253 - H3 K27cr
M22	3.69704375	ab179254 - H3 K27me1
M23	3.084375	ab179255 - H3 K27me2 + S28p
M24	1.684454167	ab179256 - H3 K27me3
M25	3.3540125	ab179257 - H3 K27me3 + S28p
N1	2.418859375	ab179258 - H3 P30hpro
N2	1.941530208	ab179259 - H3 P30(5,5me2)
N3	1.533245833	ab179260 - H3 P30
N4	2.274192708	ab179261 - H3 P30mutA
N5	2.14550625	ab179262 - H3 T32GlcNAc
N6	2.035142708	ab179263 - H3 K36ac
N7	2.257089583	ab179264 - H3 P38hpro
N8	2.176923958	ab179265 - H3 P38(5,5me2)
N9	1.95835625	ab179266 - H3 P38
N10	3.338334375	ab179267 - H3 P38mutA
N11	2.780673958	ab179268 - H3 Y41p
N12	2.723032292	ab179269 - H3 unmod(50-80)
N13	2.878317708	ab179270 - H3 K56ac
N14	2.827744792	ab179271 - H3 K56cr
N15	3.151133333	ab179272 - H3 K56fo
N16	1.860380208	ab179273 - H3 K56me1
N17	1.583971875	ab179274 - H3 K56me2
N18	1.748153125	ab179275 - H3 K56me3
N19	1.829638542	ab179276 - H3 R63me1
N20	2.253048958	ab179277 - H3 R63me2s
N21	1.643117708	ab179278 - H3 R63me2a
N22	1.616808333	ab179279 - H3 K64ac
N23	2.449703125	ab179280 - H3 K64me2
N24	2.779117708	ab179281 - H3 K64me3
N25	1.607975	ab179282 - H3 unmod(73-103)
O1	3.517475	ab179283 - H3 K79ac
O2	2.240701042	ab179284 - H3 K79fo
O3	2.180655208	ab179285 - H3 T80p
O4	3.764229167	ab179286 - H3 K79me3 + T80p
O5	3.086875	ab179287 - H3 unmod(99-135)
O6	3.028460417	ab179288 - H3 T118O-ac
O7	2.601252083	ab179289 - H3 K122fo
O8	3.522491667	ab179290 - H3 K122me1
O9	2.15095	ab179291 - H3 R128me1
O10	2.06085	ab179292 - H3 R128me2s
O11	2.040125	ab179293 - H3.3 unmod(10-40)
O12	2.370739583	ab179294 - H3.3 S28p
O13	1.896372917	ab179295 - H3.3 S28p + S31p
O14	1.938721875	ab179296 - H3.3 S31p
O15	1.546339583	ab179297 - H4 unmod(1-30)
O16	1.764989583	ab179298 - H4 S1p + R3me1

O17	1.976098958	ab179299 - H4 S1p + R3me2s
O18	1.838551042	ab179300 - H4 S1p + R3me2a
O19	2.047134375	ab179301 - H4 S1p + K5ac + K8ac + K12ac + K16ac
O20	2.18123125	ab179302 - H4 R3cit
O21	2.854855208	ab179303 - H4 R3me1 + K5ac + K8ac + K12ac + K16ac + K20ac
O22	2.099221875	ab179304 - H4 R3me2s
O23	2.726571875	ab179305 - H4 R3me2s + K5ac + K8ac + K12ac + K16ac + K20ac
O24	1.76918125	ab179306 - H4 R3me2a
O25	2.485678125	ab179307 - H4 R3me2a + K5ac
P1	2.669957292	ab179308 - H4 R3me2a + K5ac + K8ac + K12ac + K16ac + K20ac
P2	2.600435417	ab179309 - H4 K5ac
P3	2.496847917	ab179310 - H4 K5ac + K8ac
P4	2.10533125	ab179311 - H4 K5ac + K12ac
P5	2.337605208	ab179312 - H4 K5ac + K16ac
P6	2.240929167	ab179313 - H4 K5ac + K8ac + K12ac
P7	1.838789583	ab179314 - H4 K5ac + K8me1 + K12ac
P8	3.108951042	ab179315 - H4 K5ac + K8ac + K12ac + K16ac
P9	2.399401042	ab179316 - H4 K5ac + K8me1 + K12ac + K16ac
P10	1.764867708	ab179317 - H4 K5ac + K8ac + K12me1 + K16ac
P11	2.210248958	ab179318 - H4 K5me1
P12	1.979766667	ab179319 - H4 K5me1 + K8ac + K12me1
P13	2.460635417	ab179320 - H4 K5me1 + K8me1 + K12me1
P14	2.282497917	ab179321 - H4 K5me1 + K8ac + K12ac + K16ac
P15	1.849186458	ab179322 - H4 K5me2
P16	2.33596875	ab179323 - H4 K5me3
P17	2.26660625	ab179324 - H4 K5cr
P18	1.757228125	ab179325 - H4 K8ac + K12ac
P19	2.359778125	ab179326 - H4 K8ac + K16ac
P20	1.824401042	ab179327 - H4 K8cr
P21	1.9172125	ab179328 - H4 K8me1
P22	1.814091667	ab179329 - H4 K12ac
P23	1.918078125	ab179330 - H4 K12ac + K16ac
P24	2.337555208	ab179331 - H4 K12ac + K16ac + K20me2
P25	2.082225	ab179332 - H4 K12ac + K16ac + K20me3
Q1	2.389679167	ab179333 - H4 K5mutQ + K8mutQ + K12ac + K16mutQ + K20mutQ
Q2	1.708408333	ab179334 - H4 K12cr
Q3	2.543702083	ab179335 - H4 K12fo
Q4	2.0021625	ab179336 - H4 K12me1
Q5	2.573747917	ab179337 - H4 K12me2
Q6	1.977644792	ab179338 - H4 K12me3
Q7	3.067117708	ab179339 - H4 K16ac
Q8	1.644044792	ab179340 - H4 K16ac + K20ac
Q9	3.142095833	ab179341 - H4 K16cr
Q10	1.962809375	ab179342 - H4 K16me1
Q11	1.885189583	ab179343 - H4 K16me3
Q12	2.241472917	ab179344 - H4 R17me1
Q13	1.949620833	ab179345 - H4 R17me2s
Q14	1.666889583	ab179346 - H4 R17me2a
Q15	2.099095833	ab179347 - H4 R19me1
Q16	1.540913542	ab179348 - H4 R19me2s
Q17	2.4821625	ab179349 - H4 R19me2a
Q18	1.791645833	ab179350 - H4 K20ac
Q19	2.519889583	ab179351 - H4 R23me1
Q20	1.871838542	ab179352 - H4 R23me2s
Q21	2.08735625	ab179353 - H4 R23me2a
Q22	2.413040625	ab179354 - H4 unmod(24-54)
Q23	1.61115	ab179355 - H4 K31fo
Q24	2.045827083	ab179356 - H4 K31me1
Q25	1.594813542	ab179357 - H4 K31me2
R1	2.181760417	ab179358 - H4 K31me3
R2	2.583303125	ab179359 - H4 P32hyd
R3	2.005103125	ab179360 - H4 R35me1
R4	2.350820833	ab179361 - H4 R35me2s
R5	2.373841667	ab179362 - H4 R35me2a
R6	2.670292708	ab179363 - H4 S47O-ac
R7	2.000008333	ab179364 - H4 S47GlcNAc
R8	2.803738542	ab179365 - H4 S47p
R9	2.389666667	ab179366 - H4 unmod(45-75)
R10	2.690629167	ab179367 - H4 Y51hyd
R11	1.884253125	ab179368 - H4 R55me1
R12	2.412121875	ab179369 - H4 R55me2s
R13	2.761508333	ab179370 - H4 R55me2a
R14	1.695129167	ab179371 - H4 K59fo
R15	2.056147917	ab179372 - H4 K59me1

R16	2.429054167	ab179373 - H4 K59me2
R17	2.3732625	ab179374 - H4 R67me1
R18	2.851017708	ab179375 - H4 R67me2s
R19	1.935140625	ab179376 - H4 R67me2a
R20	2.435182292	ab179377 - H4 unmod(71-102)
R21	1.860978125	ab179378 - H4 K77me1
R22	1.658884375	ab179379 - H4 K77me2
R23	1.365257292	ab179380 - H4 K77me3
R24	2.080257292	ab179381 - H4 K79fo
R25	1.8927875	ab179382 - H4 K79me1
S1	2.2606875	ab179383 - H4 K79me3
S2	2.577778125	ab179384 - H4 T82O-ac
S3	2.451188542	ab179385 - H4 Y88hyd
S4	1.672505208	ab179386 - H4 Y88nitro
S5	2.3235375	ab179387 - H4 K91fo
S6	1.477883333	ab179403 - H2A K5ac
S7	1.824152083	ab179404 - H2A K9me1
S8	4.521478125	ab18504 - H2B S32p
S9	1.929371875	ab19828 - ScH2A S129p
S10	2.229075	ab20631 - H3 R26cit
S11	1.859577083	ab20635 - H3 T3p + K4me3
S12	2.020701042	ab21043 - H3 K37me1
S13	2.183486458	ab21998 - H1.4 K25me2
S14	2.146840625	ab22079 - H4 K59me3
S15	2.581704167	ab22080 - H3 K122me2
S16	2.513359375	ab22082 - H3 K122me3
S17	2.019053125	ab22214 - H3 K37me2
S18	1.496786458	ab22398 - H2A R3me2a
S19	1.922555208	ab22399 - H2A R3me2s
S20	1.635582292	ab23539 - H3 T22p
S21	3.332015625	ab24003 - H3 K18ac
S22	2.2701625	ab24404 - H3 K27ac
S23	2.537696875	ab24417 - H3 K37me3
S24	2.338459375	ab24444 - H3 T11p
S25	1.429032292	ab24659 - H2A K127me3
T1	1.694569792	ab24660 - H2A K125me3
T2	3.203235417	ab26307 - H3 T45p
T3	2.149566667	ab26342 - H2B K5me2
T4	2.644775	ab26739 - H4 T30p
T5	2.539854167	ab27896 - H4 T96p
T6	2.468447917	ab27898 - H4 T80p
T7	1.661251042	ab28756 - H2A R3me1
T8	2.593736458	ab32876 - H3 R2cit + R8cit + R17cit
T9	1.491244792	ab33047 - H2A K127me2
T10	2.392328125	ab33505 - H3 T118p
T11	1.860070833	ab33594 - H3 R128me2a
T12	4.066158333	ab34466 - H3 K122ac
T13	1.631448958	ab36927 - H2A S122p
T14	1.91096875	ab4555 - H3 K79me1
T15	2.024275	ab4556 - H3 K79me2
T16	2.8303125	ab4557 - H3 K79me3
T17	2.644610417	ab4560 - H4 K79me2
T18	2.664478125	ab46854 - H3 K18me3
T19	1.515486458	ab48359 - H3 K23ac
T20	1.924902083	ab53244 - H3 K64me1
T21	1.366633333	ab54016 - H2A K9ac
T22	1.369008333	ab5499 - H3 S28p
T23	1.568944792	ab7768 - H3 K4me2
T24	1.498538542	ab89018 - H1.4 S26p
T25	2.24079375	ab92374 - H3 K4me3 (HEX1)
U1	1.988851042	ab95444 - H4 K16me2