

Peptide array: Histone H3 (ab322707)

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

Cell Ref	AUC	Peptide
A1	24.78918229	ab10112 - H4 K91ac
A2	19.360675	ab10121 - H2B K85ac
A3	25.63813646	ab10127 - H2A K95me2
A4	22.39342604	ab10138 - H1.4 T17p
A5	22.02428333	ab10139 - H1.4 T146p
A6	19.78778333	ab113997 - H2A.X Y142p
A7	27.03007396	ab11477 - H3 S10p
A8	23.64616042	ab12946 - H2B K5me3
A9	23.64844375	ab12949 - H2B K23me2
A10	23.41467708	ab12952 - H2A K127me1
A11	19.74412708	ab13211 - H2B K5me1
A12	32.58234271	ab1340 - H3 K4me1
A13	21.24715625	ab13570 - H2B K43me1
A14	26.95502813	ab13835 - HA
A15	23.81414375	ab13837 - Myc
A16	22.37403542	ab140408 - ScH2B T129p
A17	19.15310938	ab14103 - H3 T6p
A18	28.05788542	ab14724 - H4 S1p
A19	24.79400729	ab14799 - H3 T32p
A20	27.79130313	ab14943 - His
A21	23.66599167	ab14952 - H2A R77me2s
A22	24.55741979	ab14964 - H4 K20me2
A23	19.62827604	ab14997 - ScH2A K21ac
A24	24.38902917	ab154782 - H3 R26me2s
A25	18.23495625	ab15591 - H3 K14ac
B1	26.99634271	ab15644 - H3 K9me3 + S10p
B2	20.87027604	ab15645 - H2A.X S139p
B3	28.68214583	ab15824 - H4 K8ac
B4	26.55403438	ab15829 - V5
B5	21.36305208	ab16635 - H3 K9ac
B6	25.39332604	ab166685 - H2B S112GlcNAc
B7	23.24608542	ab166688 - H2B S36GlcNAc
B8	20.96801042	ab16935 - H3 R17me2a
B9	21.02521354	ab17043 - H4 K20me1
B10	24.08291354	ab17162 - H3 P16hyd + P30hyd
B11	22.52328229	ab17567 - H4 K20me3
B12	17.92755313	ab17587 - H1.4 K25me3
B13	19.21597604	ab175905 - H2B K11ac
B14	19.95535104	ab176211 - H2B K23ac
B15	19.67770938	ab17632 - H4 K5bio
B16	20.65973438	ab1771 - H3 K9me1
B17	23.45919271	ab1773 - H3 K9me3
B18	23.31455208	ab17770 - H4 R3me1
B19	23.16252396	ab1781 - H3 K27me2
B20	26.47257813	ab1783 - H3 K36me1
B21	49.62822708	ab1784 - H3 K36me2
B22	24.88423542	ab1785 - H3 K36me3
B23	31.36759896	ab178667 - ScH2A unmod(1-30)
B24	31.75581979	ab178668 - ScH2A unmod(102-132)
B25	35.67399792	ab178670 - H2A.Z unmod(1-30)
C1	27.29114688	ab178833 - Flag
C2	21.35411458	ab178835 - H2B R99mutO
C3	18.48047083	ab178836 - H4 R23mutO
C4	25.86626771	ab178837 - H4 G48mutD
C5	19.65697813	ab178987 - H1.4 unmod(1-30)
C6	20.08887083	ab178988 - H1.4 S1p
C7	19.55468125	ab178989 - H1.4 T3p
C8	20.81123125	ab178990 - H1.4 S1p + T3p
C9	17.83756667	ab178991 - H1.4 K25me1
C10	24.29948021	ab178992 - H1.4 K25me3 + S26p
C11	24.72384375	ab178993 - H1.4 unmod(130-160)
C12	30.02948333	ab178994 - H2A unmod(1-30)
C13	21.39154896	ab178995 - H2A S1p
C14	27.54289688	ab178996 - H2A S1p + K5ac
C15	18.77704167	ab178997 - H2A S1p + R3me2a + K5ac
C16	29.96734375	ab178998 - H2A S1p + R3me2s + K5ac
C17	21.64190521	ab178999 - H2A S1p + R3cit + K5ac
C18	27.46275208	ab179000 - H2A S1ac + S1p + K5ac + K9ac + K13ac + K15ac
C19	26.61559688	ab179001 - H2A S1p + K5ac + K9ac + K13ac + K15ac
C20	22.44910313	ab179002 - H2A R3cit

C21	26.32180729	ab179003 - H2A R3cit + K5ac
C22	21.76536458	ab179004 - H2A R3me2a + K5ac
C23	24.47249896	ab179005 - H2A R3me2s + K5ac
C24	28.80828229	ab179006 - H2A S1ac + K5ac + K9ac + K13ac + K15ac
C25	22.36893229	ab179007 - H2A K5ac + K9ac + K13ac + K15ac
D1	19.86817813	ab179008 - H2A K9me2
D2	17.91825417	ab179009 - H2A K9me3
D3	20.70647396	ab179010 - H2A R11me1
D4	20.08344792	ab179011 - H2A R11me2s
D5	16.40704063	ab179012 - H2A R11me2a
D6	19.41063125	ab179013 - H2A K13ac
D7	21.77128854	ab179014 - H2A K15ac
D8	16.51744688	ab179015 - H2A R17me1
D9	21.78657292	ab179016 - H2A R17me2s
D10	22.38338333	ab179017 - H2A R17me2a
D11	26.76084375	ab179018 - H2A unmod(20-50)
D12	23.99521667	ab179019 - H2A P26hyd
D13	29.69304688	ab179020 - H2A R29me1
D14	21.027625	ab179021 - H2A R29me2s
D15	25.15008958	ab179022 - H2A R29me2a
D16	24.21560313	ab179023 - H2A K36ac
D17	29.94605208	ab179024 - H2A K36cr
D18	21.96554271	ab179025 - H2A Y39hyd
D19	29.45583375	ab179026 - H2A R42me1
D20	21.25075	ab179027 - H2A R42me2s
D21	28.51830729	ab179028 - H2A R42me2a
D22	30.42382188	ab179029 - H2A unmod(71-103)
D23	28.37344375	ab179030 - H2A R77me1
D24	27.80694688	ab179031 - H2A R77me2a
D25	23.76982917	ab179032 - H2A R88me1
E1	25.09104896	ab179033 - H2A R88me2s
E2	25.79714375	ab179034 - H2A R88me2a
E3	29.94533021	ab179035 - H2A K95me1
E4	28.08759896	ab179036 - H2A K95me3
E5	22.64292292	ab179037 - H2A unmod(92-123)
E6	21.28470521	ab179038 - H2A T101GlcNAc
E7	20.23732083	ab179039 - H2A Q104me1
E8	21.55106042	ab179040 - H2A unmod(112-129)
E9	28.94952396	ab179041 - H2A K118me1
E10	21.95983438	ab179042 - H2A K118me2
E11	21.37289271	ab179043 - H2A K118me3
E12	23.13192604	ab179044 - H2A K118fo
E13	17.6173625	ab179045 - H2A K118cr
E14	28.78450417	ab179046 - H2A K119cr
E15	22.13557396	ab179047 - H2A K118ubiq
E16	25.0645125	ab179048 - H2A K119ubiq
E17	20.03633021	ab179049 - H2A T120p
E18	20.77351979	ab179050 - H2A K125me1
E19	26.83918958	ab179051 - H2A K125me2
E20	28.64177292	ab179052 - H2A K125cr
E21	32.84385729	ab179053 - H2A.X unmod(1-30)
E22	21.42846771	ab179054 - H2A.X K5ac
E23	44.94209792	ab179055 - H2A.X unmod(114-142)
E24	22.554	ab179056 - H2A.X K118ubiq
E25	25.92078958	ab179057 - H2A.X K119ubiq
F1	21.62043542	ab179058 - H2A.X T120p
F2	28.11683125	ab179059 - H2A.X K127ubiq
F3	28.85460208	ab179060 - H2A.X S139p + Y142p
F4	18.49929583	ab179061 - H2A.Z K4ac
F5	25.15107292	ab179062 - H2A.Z K7ac
F6	15.9344125	ab179063 - H2A.Z K11ac
F7	26.51662083	ab179064 - H2A.Z K4ac + K7ac + K11ac
F8	21.37414375	ab179065 - H2B unmod(1-30)
F9	24.575575	ab179066 - H2B K5ac
F10	22.82919896	ab179067 - H2B K5cr
F11	26.85807396	ab179068 - H2B K5fo
F12	21.98067708	ab179069 - H2B P10hyd
F13	16.81701979	ab179070 - H2B K11me1
F14	29.94336458	ab179071 - H2B K11me2
F15	19.55211771	ab179072 - H2B K11me3
F16	22.60729792	ab179073 - H2B K11cr
F17	19.4454875	ab179074 - H2B K12ac
F18	30.62425208	ab179075 - H2B K12cr
F19	17.62354896	ab179076 - H2B K12me1

F20	20.40752813	ab179077 - H2B K12me2
F21	19.64855313	ab179078 - H2B K12me3
F22	18.05706146	ab179079 - H2B K12ac + K15ac
F23	22.80521146	ab179080 - H2B S14p
F24	20.70623958	ab179081 - H2B K15ac
F25	25.02491146	ab179082 - H2B K15cr
G1	18.63254375	ab179083 - H2B K15me1
G2	23.03848333	ab179084 - H2B K15me2
G3	19.58236875	ab179085 - H2B K15me3
G4	31.61996979	ab179086 - H2B K16ac
G5	21.56411979	ab179087 - H2B K16cr
G6	25.59057813	ab179088 - H2B K20ac
G7	21.29276042	ab179089 - H2B K20cr
G8	31.16020208	ab179090 - H2B K20me1
G9	18.55382292	ab179091 - H2B K20me2
G10	30.8161375	ab179092 - H2B K20me3
G11	24.29329167	ab179093 - H2B K23cr
G12	20.35970104	ab179094 - H2B K23me1
G13	21.22806146	ab179095 - H2B K23me3
G14	23.82849375	ab179096 - H2B K24ac
G15	24.60159688	ab179097 - H2B unmod(26-56)
G16	20.09103125	ab179098 - H2B K34cr
G17	21.29385729	ab179099 - H2B K34fo
G18	18.14266979	ab179100 - H2B Y37hyd
G19	22.84531042	ab179101 - H2B Y37p
G20	23.61993125	ab179102 - H2B K43me2
G21	19.28024375	ab179103 - H2B K43me3
G22	21.51534167	ab179104 - H2B K46fo
G23	20.34216667	ab179105 - H2B K46ac
G24	17.80785104	ab179106 - H2B K46me1
G25	22.13904792	ab179107 - H2B K46me2
H1	17.68582396	ab179108 - H2B K46me3
H2	20.92940729	ab179109 - H2B unmod(51-81)
H3	20.04999896	ab179110 - H2B K57me1
H4	21.03649583	ab179111 - H2B K57me2
H5	19.46676563	ab179112 - H2B K57me3
H6	29.77579167	ab179113 - H2B unmod(73-103)
H7	22.53927188	ab179114 - H2B R79me1
H8	30.28938438	ab179115 - H2B R79me2s
H9	20.88614896	ab179116 - H2B R79me2a
H10	28.76848229	ab179117 - H2B Y83hyd
H11	28.47767292	ab179118 - H2B K85me1
H12	21.23365208	ab179119 - H2B K85me2
H13	23.43705729	ab179120 - H2B K85me3
H14	26.21392813	ab179121 - H2B unmod(94-125)
H15	25.33350729	ab179122 - H2B R99me1
H16	19.08612292	ab179123 - H2B R99me2s
H17	27.60941146	ab179124 - H2B R99me2a
H18	24.03639583	ab179125 - H2B K108fo
H19	17.20253438	ab179126 - H2B K116fo
H20	23.19443958	ab179127 - H2B K116me1
H21	19.9381625	ab179128 - H2B K116me2
H22	19.2425125	ab179129 - H2B K116me3
H23	20.60062604	ab179130 - H2B K116ubiq
H24	26.08532813	ab179131 - H2B K120ac
H25	18.09541563	ab179132 - H2B K120fo
I1	27.21955938	ab179133 - H2B K120ubiq
I2	23.35026563	ab179134 - H2B K125ubiq
I3	22.50593229	ab179135 - Sch2B unmod(102-132)
I4	18.54859896	ab179136 - H3 unmod(1-33)
I5	22.78807188	ab179137 - H3 R2cit
I6	17.63820208	ab179138 - H3 R2cit + K4me2
I7	25.02508646	ab179139 - H3 R2cit + K4me3
I8	19.10753958	ab179140 - H3 R2cit + K4me3 + K9ac + K14ac + K18ac
I9	20.53875938	ab179141 - H3 R2cit + K4ac + K9ac + K14ac + K18ac
I10	24.80931042	ab179142 - H3 R2me1
I11	21.12379063	ab179143 - H3 R2me1 + K4me2
I12	23.68772083	ab179144 - H3 R2me1 + K4me3
I13	19.92042604	ab179145 - H3 R2me1 + K4me3 + K9ac + K14ac + K18ac
I14	22.10538333	ab179146 - H3 R2me2s
I15	21.02160625	ab179147 - H3 R2me2s + K4me2
I16	29.15571875	ab179148 - H3 R2me2s + K4me3
I17	21.03712083	ab179149 - H3 R2me2s + K4me3 + K9ac + K14ac + K18ac
I18	27.13250938	ab179150 - H3 R2me2a

I19	21.78451042	ab179151 - H3 R2me2a + K4me2
I20	28.29882917	ab179152 - H3 R2me2a + K4me3
I21	21.47895625	ab179153 - H3 R2me2a + K4me3 + S10p
I22	24.84709167	ab179154 - H3 R2me2a + T3p + K4me3
I23	22.0301125	ab179155 - H3 R2me2a + K4ac + K9ac + K14ac + K18ac
I24	20.02995	ab179156 - H3 R2me2a + K4me3 + K9ac + K14ac + K18ac
I25	19.83889583	ab179157 - H3 R2me2a + T3p + K4me3 + K9ac + K14ac + K18ac
J1	19.93056042	ab179158 - H3 R2me2a + K4me3 + K9ac + S10p + K14ac + K18ac
J2	20.21403021	ab179159 - H3 T3p
J3	20.2919125	ab179160 - H3 T3p + K4me1
J4	25.2191375	ab179161 - H3 T3p + K4me2
J5	22.23490417	ab179162 - H3 T3p + K4me3 + K9ac + K14ac + K18ac
J6	23.30848229	ab179163 - H3 K4ac
J7	21.11202604	ab179164 - H3 K4ac + K9ac
J8	22.80060938	ab179165 - H3 K4ac + K14ac
J9	23.79719688	ab179166 - H3 K4ac + K18ac
J10	21.08772396	ab179167 - H3 K4ac + K9me3
J11	21.17189583	ab179168 - H3 K4ac + K9ac + K14ac
J12	23.76649167	ab179169 - H3 K4ac + K9ac + K18ac
J13	18.53393854	ab179170 - H3 K4ac + K14ac + K18ac
J14	28.69117604	ab179171 - H3 K4ac + K9ac + K14ac + K18ac
J15	20.57155208	ab179172 - H3 K4ac + K9me1 + K14ac + K18ac
J16	28.30051146	ab179173 - H3 K4ac + K9me2 + K14ac + K18ac
J17	21.87980313	ab179174 - H3 K4ac + K9me3 + K14ac + K18ac
J18	30.32125729	ab179175 - H3 K4ac + K9ac + S10p + K14ac + K18ac
J19	21.98540417	ab179176 - H3 K4cr
J20	21.58283438	ab179177 - H3 K4me1 + K9me2
J21	21.35271458	ab179178 - H3 K4me1 + K9ac + K18ac
J22	21.54546771	ab179179 - H3 K4me1 + K9ac + K14ac + K18ac
J23	23.61628125	ab179180 - H3 K4me2 + T6p
J24	21.02601771	ab179181 - H3 K4me2 + K9me2
J25	21.62718333	ab179182 - H3 K4me2 + K9ac + K18ac
K1	20.82461042	ab179183 - H3 K4me2 + K9ac + K14ac + K18ac
K2	23.01817292	ab179184 - H3 K4me3 + T6p
K3	20.70977917	ab179185 - H3 K4me3 + K9ac
K4	23.29195938	ab179186 - H3 K4me3 + K9me2
K5	22.6339	ab179187 - H3 K4me3 + K9me3
K6	21.37067188	ab179188 - H3 K4me3 + S10p
K7	23.21611146	ab179189 - H3 K4me3 + K14ac
K8	25.727525	ab179190 - H3 K4me3 + K18ac
K9	20.89543958	ab179191 - H3 K4me3 + R8me2s + K9me3
K10	27.08402813	ab179192 - H3 K4me3 + R8me2a + K9me3
K11	21.01624167	ab179193 - H3 K4me3 + K9ac + S10p
K12	28.28730521	ab179194 - H3 K4me3 + K9ac + K14ac
K13	26.98844896	ab179195 - H3 K4me3 + K9ac + K18ac
K14	28.27766146	ab179196 - H3 K4me3 + K14ac + K18ac
K15	22.77858333	ab179197 - H3 K4me3 + K9ac + K14ac + K18ac
K16	25.14620417	ab179198 - H3 K4me3 + K9ac + S10p + K14ac + K18ac
K17	22.5677125	ab179199 - H3 K4me3 + T6p + K9ac + K14ac + K18ac
K18	21.89885833	ab179200 - H3 K4me2 + T6p + K9ac + K14ac + K18ac
K19	20.98181458	ab179201 - H3 K4bio + K9bio + K18bio
K20	19.68787292	ab179202 - H3 T6p + K9me3
K21	19.06365417	ab179203 - H3 T6p + R8me2a + K9me3
K22	17.92478333	ab179204 - H3 T6p + K9ac + K14ac + K18ac
K23	24.32064063	ab179205 - H3 R8me1
K24	21.32565313	ab179206 - H3 R8me1 + K9me1
K25	18.29707188	ab179207 - H3 R8me1 + K9me2
L1	20.29017083	ab179208 - H3 R8me1 + K9me3
L2	20.21196146	ab179209 - H3 R8me2s
L3	19.79919375	ab179210 - H3 R8me2s + K9me1
L4	24.45046875	ab179211 - H3 R8me2s + K9me2
L5	22.85658125	ab179212 - H3 R8me2s + K9me3
L6	23.73816771	ab179213 - H3 R8me2a
L7	22.58610521	ab179214 - H3 R8me2a + K9me1
L8	25.25609479	ab179215 - H3 R8me2a + K9me2
L9	19.14749583	ab179216 - H3 R8me2a + K9me3
L10	24.95626146	ab179217 - H3 R8cit
L11	34.30202917	ab179218 - H3 K9ac + K14ac
L12	24.87060521	ab179219 - H3 K9ac + K18ac
L13	26.42460729	ab179220 - H3 K9ac + K14ac + K18ac
L14	22.88046667	ab179221 - H3 K9ac + K14ac + K18ac + K23ac + K27ac
L15	21.12057396	ab179222 - H3 K9ac + S10p
L16	20.01620729	ab179223 - H3 K9cr
L17	33.61066042	ab179224 - H3 K9me2

L18	26.34525104	ab179225 - H3 K9me2 + S10p
L19	18.96572708	ab179226 - H3 K9me2 + K27me2
L20	24.55749479	ab179227 - H3 S10p + S28p
L21	29.0099625	ab179228 - H3 S10p K14ac
L22	29.77212292	ab179229 - H3 K14me1
L23	21.84650417	ab179230 - H3 K14me2
L24	26.47332604	ab179231 - H3 K14me3
L25	27.91812604	ab179232 - H3 K14ac + K18ac
M1	25.92397083	ab179233 - H3 R17cit
M2	21.65769167	ab179234 - H3 R17me1
M3	24.41118438	ab179235 - H3 R17me2s
M4	23.22850104	ab179236 - H3 R17me2a + K18ac
M5	25.32016458	ab179237 - H3 K18cr
M6	20.50351354	ab179238 - H3 K18me1
M7	25.08049271	ab179239 - H3 K18me2
M8	25.69542708	ab179240 - H3 K18me3 + K36me3
M9	26.1734875	ab179241 - H3 K23cr
M10	24.83345729	ab179242 - H3 K23me1
M11	20.80305938	ab179243 - H3 K23me2
M12	26.02563229	ab179244 - H3 K23me3
M13	22.29225521	ab179245 - H3 T22p + K23me3
M14	27.37343542	ab179246 - H3 K23pr
M15	8027.083985	ab179247 - H3 unmod(20-51)
M16	3972.925809	ab179248 - H3 R26me1
M17	98.21777917	ab179249 - H3 R26me2a
M18	22.93216458	ab179250 - H3 R26me2a + K27me1
M19	32.1519125	ab179251 - H3 R26me2a + K27me3
M20	20.92869583	ab179252 - H3 K27ac + S28p
M21	22.94941979	ab179253 - H3 K27cr
M22	20.28028333	ab179254 - H3 K27me1
M23	19.5866	ab179255 - H3 K27me2 + S28p
M24	20.53749688	ab179256 - H3 K27me3
M25	20.687775	ab179257 - H3 K27me3 + S28p
N1	17.97009792	ab179258 - H3 P30hpro
N2	23.79278333	ab179259 - H3 P30(5,5me2)
N3	23.50453438	ab179260 - H3 P30
N4	28.05238229	ab179261 - H3 P30mutA
N5	18.49477604	ab179262 - H3 T32GlcNAc
N6	24.14704063	ab179263 - H3 K36ac
N7	27.24277083	ab179264 - H3 P38hpro
N8	30.55848958	ab179265 - H3 P38(5,5me2)
N9	22.14692083	ab179266 - H3 P38
N10	20.72336042	ab179267 - H3 P38mutA
N11	8481.789281	ab179268 - H3 Y41p
N12	1474.264555	ab179269 - H3 unmod(50-80)
N13	23.20439479	ab179270 - H3 K56ac
N14	34.48767292	ab179271 - H3 K56cr
N15	22.60652708	ab179272 - H3 K56fo
N16	24.74694063	ab179273 - H3 K56me1
N17	23.98	ab179274 - H3 K56me2
N18	24.07885313	ab179275 - H3 K56me3
N19	22.64075313	ab179276 - H3 R63me1
N20	25.03136667	ab179277 - H3 R63me2s
N21	21.96552083	ab179278 - H3 R63me2a
N22	25.32442813	ab179279 - H3 K64ac
N23	25.49155104	ab179280 - H3 K64me2
N24	31.56857083	ab179281 - H3 K64me3
N25	29.32381875	ab179282 - H3 unmod(73-103)
O1	21.02818125	ab179283 - H3 K79ac
O2	27.53301667	ab179284 - H3 K79fo
O3	22.18107708	ab179285 - H3 T80p
O4	22.64775521	ab179286 - H3 K79me3 + T80p
O5	2379.017616	ab179287 - H3 unmod(99-135)
O6	19.5827	ab179288 - H3 T118O-ac
O7	24.52834896	ab179289 - H3 K122fo
O8	22.64586042	ab179290 - H3 K122me1
O9	18.97676667	ab179291 - H3 R128me1
O10	21.96495729	ab179292 - H3 R128me2s
O11	24.39971979	ab179293 - H3.3 unmod(10-40)
O12	19.16110208	ab179294 - H3.3 S28p
O13	24.84499271	ab179295 - H3.3 S28p + S31p
O14	24.90570417	ab179296 - H3.3 S31p
O15	35.15711563	ab179297 - H4 unmod(1-30)
O16	26.73931458	ab179298 - H4 S1p + R3me1

O17	24.44729583	ab179299 - H4 S1p + R3me2s
O18	26.36045521	ab179300 - H4 S1p + R3me2a
O19	28.59991667	ab179301 - H4 S1p + K5ac + K8ac + K12ac + K16ac
O20	28.24543854	ab179302 - H4 R3cit
O21	19.93648854	ab179303 - H4 R3me1 + K5ac + K8ac + K12ac + K16ac + K20ac
O22	23.68443958	ab179304 - H4 R3me2s
O23	21.08767188	ab179305 - H4 R3me2s + K5ac + K8ac + K12ac + K16ac + K20ac
O24	26.13278333	ab179306 - H4 R3me2a
O25	18.22662292	ab179307 - H4 R3me2a + K5ac
P1	25.48469063	ab179308 - H4 R3me2a + K5ac + K8ac + K12ac + K16ac + K20ac
P2	21.07189688	ab179309 - H4 K5ac
P3	25.04807604	ab179310 - H4 K5ac + K8ac
P4	26.70764896	ab179311 - H4 K5ac + K12ac
P5	24.97289896	ab179312 - H4 K5ac + K16ac
P6	26.22320729	ab179313 - H4 K5ac + K8ac + K12ac
P7	24.77554792	ab179314 - H4 K5ac + K8me1 + K12ac
P8	26.91664063	ab179315 - H4 K5ac + K8ac + K12ac + K16ac
P9	28.22732917	ab179316 - H4 K5ac + K8me1 + K12ac + K16ac
P10	26.18408854	ab179317 - H4 K5ac + K8ac + K12me1 + K16ac
P11	23.34759479	ab179318 - H4 K5me1
P12	21.39831354	ab179319 - H4 K5me1 + K8ac + K12me1
P13	23.27101563	ab179320 - H4 K5me1 + K8me1 + K12me1
P14	22.67784688	ab179321 - H4 K5me1 + K8ac + K12ac + K16ac
P15	22.95847396	ab179322 - H4 K5me2
P16	21.93722396	ab179323 - H4 K5me3
P17	23.52658438	ab179324 - H4 K5cr
P18	25.53342813	ab179325 - H4 K8ac + K12ac
P19	27.36685313	ab179326 - H4 K8ac + K16ac
P20	24.1420625	ab179327 - H4 K8cr
P21	27.66034896	ab179328 - H4 K8me1
P22	25.56112188	ab179329 - H4 K12ac
P23	27.14742604	ab179330 - H4 K12ac + K16ac
P24	27.70226667	ab179331 - H4 K12ac + K16ac + K20me2
P25	19.60447813	ab179332 - H4 K12ac + K16ac + K20me3
Q1	27.14480938	ab179333 - H4 K5mutQ + K8mutQ + K12ac + K16mutQ + K20mutQ
Q2	22.33927396	ab179334 - H4 K12cr
Q3	26.60723125	ab179335 - H4 K12fo
Q4	22.22731979	ab179336 - H4 K12me1
Q5	23.15142813	ab179337 - H4 K12me2
Q6	19.61710521	ab179338 - H4 K12me3
Q7	23.0675375	ab179339 - H4 K16ac
Q8	24.82580938	ab179340 - H4 K16ac + K20ac
Q9	24.45053229	ab179341 - H4 K16cr
Q10	25.55500104	ab179342 - H4 K16me1
Q11	23.39557708	ab179343 - H4 K16me3
Q12	27.43279896	ab179344 - H4 R17me1
Q13	24.94617708	ab179345 - H4 R17me2s
Q14	26.56150313	ab179346 - H4 R17me2a
Q15	23.86848646	ab179347 - H4 R19me1
Q16	19.06640729	ab179348 - H4 R19me2s
Q17	24.03040104	ab179349 - H4 R19me2a
Q18	20.44259688	ab179350 - H4 K20ac
Q19	40.02120938	ab179351 - H4 R23me1
Q20	24.00086563	ab179352 - H4 R23me2s
Q21	23.16208854	ab179353 - H4 R23me2a
Q22	28.63788125	ab179354 - H4 unmod(24-54)
Q23	27.18739271	ab179355 - H4 K31fo
Q24	25.06395313	ab179356 - H4 K31me1
Q25	25.96334479	ab179357 - H4 K31me2
R1	20.19070521	ab179358 - H4 K31me3
R2	29.0510875	ab179359 - H4 P32hyd
R3	20.72714479	ab179360 - H4 R35me1
R4	24.60975313	ab179361 - H4 R35me2s
R5	22.94391771	ab179362 - H4 R35me2a
R6	24.77672708	ab179363 - H4 S47O-ac
R7	22.37905208	ab179364 - H4 S47GlcNAc
R8	26.61961771	ab179365 - H4 S47p
R9	26.03693958	ab179366 - H4 unmod(45-75)
R10	25.80074063	ab179367 - H4 Y51hyd
R11	23.77859375	ab179368 - H4 R55me1
R12	26.98185313	ab179369 - H4 R55me2s
R13	25.83044063	ab179370 - H4 R55me2a
R14	26.79844479	ab179371 - H4 K59fo
R15	23.60818854	ab179372 - H4 K59me1

R16	25.60875625	ab179373 - H4 K59me2
R17	24.77957396	ab179374 - H4 R67me1
R18	27.52656979	ab179375 - H4 R67me2s
R19	25.74652292	ab179376 - H4 R67me2a
R20	28.45185208	ab179377 - H4 unmod(71-102)
R21	26.66643125	ab179378 - H4 K77me1
R22	26.59236667	ab179379 - H4 K77me2
R23	25.10762604	ab179380 - H4 K77me3
R24	27.50699479	ab179381 - H4 K79fo
R25	25.65753021	ab179382 - H4 K79me1
S1	26.58709583	ab179383 - H4 K79me3
S2	25.66983542	ab179384 - H4 T82O-ac
S3	26.66683229	ab179385 - H4 Y88hyd
S4	26.06234063	ab179386 - H4 Y88nitro
S5	24.229225	ab179387 - H4 K91fo
S6	24.3133	ab179403 - H2A K5ac
S7	20.38586563	ab179404 - H2A K9me1
S8	23.08885625	ab18504 - H2B S32p
S9	21.83036979	ab19828 - ScH2A S129p
S10	25.20643333	ab20631 - H3 R26cit
S11	18.03217292	ab20635 - H3 T3p + K4me3
S12	170.1163708	ab21043 - H3 K37me1
S13	21.4899625	ab21998 - H1.4 K25me2
S14	27.06451979	ab22079 - H4 K59me3
S15	20.73873958	ab22080 - H3 K122me2
S16	24.76760104	ab22082 - H3 K122me3
S17	26.87657917	ab22214 - H3 K37me2
S18	27.28237604	ab22398 - H2A R3me2a
S19	31.7688375	ab22399 - H2A R3me2s
S20	24.28643229	ab23539 - H3 T22p
S21	19.25868958	ab24003 - H3 K18ac
S22	19.52719375	ab24404 - H3 K27ac
S23	77.52236667	ab24417 - H3 K37me3
S24	18.69352188	ab24444 - H3 T11p
S25	34.61214583	ab24659 - H2A K127me3
T1	24.4361125	ab24660 - H2A K125me3
T2	370.2037063	ab26307 - H3 T45p
T3	21.48380521	ab26342 - H2B K5me2
T4	25.83334479	ab26739 - H4 T30p
T5	23.51254479	ab27896 - H4 T96p
T6	26.1370625	ab27898 - H4 T80p
T7	31.3986125	ab28756 - H2A R3me1
T8	21.88102083	ab32876 - H3 R2cit + R8cit + R17cit
T9	28.79618854	ab33047 - H2A K127me2
T10	41.98420417	ab33505 - H3 T118p
T11	19.940225	ab33594 - H3 R128me2a
T12	21.4225875	ab34466 - H3 K122ac
T13	26.4496375	ab36927 - H2A S122p
T14	23.12032708	ab4555 - H3 K79me1
T15	25.29565938	ab4556 - H3 K79me2
T16	21.56341667	ab4557 - H3 K79me3
T17	25.42798021	ab4560 - H4 K79me2
T18	19.04641667	ab46854 - H3 K18me3
T19	27.78582813	ab48359 - H3 K23ac
T20	21.82667708	ab53244 - H3 K64me1
T21	24.74278229	ab54016 - H2A K9ac
T22	25.59889167	ab5499 - H3 S28p
T23	21.94964479	ab7768 - H3 K4me2
T24	20.6315125	ab89018 - H1.4 S26p
T25	23.08368229	ab92374 - H3 K4me3 (HEX1)
U1	25.5333125	ab95444 - H4 K16me2