

Peptide array: Histone H3 (ab309551)

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

Cell Ref	AUC	Peptide
A1	16.82941771	ab10112 - H4 K91ac
A2	17.45375417	ab10121 - H2B K85ac
A3	18.35754688	ab10127 - H2A K95me2
A4	24.14633229	ab10138 - H1.4 T17p
A5	15.78250729	ab10139 - H1.4 T146p
A6	20.41273438	ab113997 - H2A.X Y142p
A7	17.31871042	ab11477 - H3 S10p
A8	16.68683646	ab12946 - H2B K5me3
A9	14.96067188	ab12949 - H2B K23me2
A10	19.10258125	ab12952 - H2A K127me1
A11	18.17192292	ab13211 - H2B K5me1
A12	23.80687917	ab1340 - H3 K4me1
A13	16.62283229	ab13570 - H2B K43me1
A14	14.07104167	ab13835 - HA
A15	16.52939271	ab13837 - Myc
A16	18.10541042	ab140408 - ScH2B T129p
A17	17.14933229	ab14103 - H3 T6p
A18	23.60830729	ab14724 - H4 S1p
A19	14.11496771	ab14799 - H3 T32p
A20	18.03208333	ab14943 - His
A21	23.5499	ab14952 - H2A R77me2s
A22	22.04369167	ab14964 - H4 K20me2
A23	18.68584167	ab14997 - ScH2A K21ac
A24	19.23958333	ab154782 - H3 R26me2s
A25	15.91529271	ab15591 - H3 K14ac
B1	16.57565833	ab15644 - H3 K9me3 + S10p
B2	20.67181771	ab15645 - H2A.X S139p
B3	23.39936458	ab15824 - H4 K8ac
B4	18.07785417	ab15829 - V5
B5	18.90080938	ab16635 - H3 K9ac
B6	16.69465938	ab166685 - H2B S112GlcNAc
B7	18.92333438	ab166688 - H2B S36GlcNAc
B8	22.38851563	ab16935 - H3 R17me2a
B9	37.84709688	ab17043 - H4 K20me1
B10	21.7786625	ab17162 - H3 P16hyd + P30hyd
B11	26.62209167	ab17567 - H4 K20me3
B12	16.11063854	ab17587 - H1.4 K25me3
B13	31.277025	ab175905 - H2B K11ac
B14	20.00269583	ab176211 - H2B K23ac
B15	22.1554625	ab17632 - H4 K5bio
B16	17.03117188	ab1771 - H3 K9me1
B17	19.21239375	ab1773 - H3 K9me3
B18	26.95069167	ab17770 - H4 R3me1
B19	17.19802292	ab1781 - H3 K27me2
B20	17.30157708	ab1783 - H3 K36me1
B21	21.53693125	ab1784 - H3 K36me2
B22	21.93781458	ab1785 - H3 K36me3
B23	21.17986354	ab178667 - ScH2A unmod(1-30)
B24	17.65728229	ab178668 - ScH2A unmod(102-132)
B25	26.07744479	ab178670 - H2A.Z unmod(1-30)
C1	18.54798125	ab178833 - Flag
C2	19.2901375	ab178835 - H2B R99mutO
C3	18.82724271	ab178836 - H4 R23mutO
C4	21.05977708	ab178837 - H4 G48mutD
C5	17.75254688	ab178987 - H1.4 unmod(1-30)
C6	14.87448333	ab178988 - H1.4 S1p
C7	16.88641979	ab178989 - H1.4 T3p
C8	16.09169167	ab178990 - H1.4 S1p + T3p
C9	17.01242813	ab178991 - H1.4 K25me1
C10	14.67116875	ab178992 - H1.4 K25me3 + S26p
C11	16.17662396	ab178993 - H1.4 unmod(130-160)
C12	24.85962188	ab178994 - H2A unmod(1-30)
C13	18.10588646	ab178995 - H2A S1p
C14	18.39273958	ab178996 - H2A S1p + K5ac
C15	13.10128958	ab178997 - H2A S1p + R3me2a + K5ac
C16	20.60088542	ab178998 - H2A S1p + R3me2s + K5ac
C17	16.90640208	ab178999 - H2A S1p + R3cit + K5ac
C18	16.43539167	ab179000 - H2A S1ac + S1p + K5ac + K9ac + K13ac + K15ac
C19	20.34262396	ab179001 - H2A S1p + K5ac + K9ac + K13ac + K15ac
C20	19.2508	ab179002 - H2A R3cit

C21	15.44335625	ab179003 - H2A R3cit + K5ac
C22	19.50013542	ab179004 - H2A R3me2a + K5ac
C23	18.86319583	ab179005 - H2A R3me2s + K5ac
C24	23.97605833	ab179006 - H2A S1ac + K5ac + K9ac + K13ac + K15ac
C25	17.44912604	ab179007 - H2A K5ac + K9ac + K13ac + K15ac
D1	16.8162875	ab179008 - H2A K9me2
D2	16.07414583	ab179009 - H2A K9me3
D3	17.71110313	ab179010 - H2A R11me1
D4	13.96710833	ab179011 - H2A R11me2s
D5	20.36205313	ab179012 - H2A R11me2a
D6	13.2341375	ab179013 - H2A K13ac
D7	19.39414688	ab179014 - H2A K15ac
D8	12.53507917	ab179015 - H2A R17me1
D9	16.31000417	ab179016 - H2A R17me2s
D10	15.46199375	ab179017 - H2A R17me2a
D11	18.06784896	ab179018 - H2A unmod(20-50)
D12	15.91301458	ab179019 - H2A P26hyd
D13	15.11452917	ab179020 - H2A R29me1
D14	13.06825833	ab179021 - H2A R29me2s
D15	19.54547292	ab179022 - H2A R29me2a
D16	14.71621875	ab179023 - H2A K36ac
D17	15.40229375	ab179024 - H2A K36cr
D18	17.38043854	ab179025 - H2A Y39hyd
D19	17.687775	ab179026 - H2A R42me1
D20	18.47908854	ab179027 - H2A R42me2s
D21	15.26136354	ab179028 - H2A R42me2a
D22	19.12574063	ab179029 - H2A unmod(71-103)
D23	17.00937813	ab179030 - H2A R77me1
D24	22.01763333	ab179031 - H2A R77me2a
D25	20.75935417	ab179032 - H2A R88me1
E1	18.44559792	ab179033 - H2A R88me2s
E2	19.20026875	ab179034 - H2A R88me2a
E3	17.26078021	ab179035 - H2A K95me1
E4	19.69418646	ab179036 - H2A K95me3
E5	17.81502188	ab179037 - H2A unmod(92-123)
E6	17.6944125	ab179038 - H2A T101GlcNAc
E7	13.76663854	ab179039 - H2A Q104me1
E8	28.48489167	ab179040 - H2A unmod(112-129)
E9	26.49947708	ab179041 - H2A K118me1
E10	19.22619479	ab179042 - H2A K118me2
E11	16.09793542	ab179043 - H2A K118me3
E12	20.09319479	ab179044 - H2A K118fo
E13	14.57369271	ab179045 - H2A K118cr
E14	17.75220417	ab179046 - H2A K119cr
E15	18.24590104	ab179047 - H2A K118ubiq
E16	17.3783375	ab179048 - H2A K119ubiq
E17	12.97353333	ab179049 - H2A T120p
E18	13.99988646	ab179050 - H2A K125me1
E19	13.28172083	ab179051 - H2A K125me2
E20	14.62981875	ab179052 - H2A K125cr
E21	35.96505313	ab179053 - H2A.X unmod(1-30)
E22	20.39024479	ab179054 - H2A.X K5ac
E23	32.43432396	ab179055 - H2A.X unmod(114-142)
E24	24.25117396	ab179056 - H2A.X K118ubiq
E25	21.76323542	ab179057 - H2A.X K119ubiq
F1	18.88846979	ab179058 - H2A.X T120p
F2	18.83543333	ab179059 - H2A.X K127ubiq
F3	23.04306875	ab179060 - H2A.X S139p + Y142p
F4	20.59856979	ab179061 - H2A.Z K4ac
F5	18.31784792	ab179062 - H2A.Z K7ac
F6	18.25434792	ab179063 - H2A.Z K11ac
F7	17.36057292	ab179064 - H2A.Z K4ac + K7ac + K11ac
F8	45.00379896	ab179065 - H2B unmod(1-30)
F9	14.84991458	ab179066 - H2B K5ac
F10	21.20533021	ab179067 - H2B K5cr
F11	19.75026875	ab179068 - H2B K5fo
F12	12.22251771	ab179069 - H2B P10hyd
F13	11.50325729	ab179070 - H2B K11me1
F14	20.44985833	ab179071 - H2B K11me2
F15	15.43933125	ab179072 - H2B K11me3
F16	18.23893125	ab179073 - H2B K11cr
F17	14.94923542	ab179074 - H2B K12ac
F18	15.96326875	ab179075 - H2B K12cr
F19	16.40172917	ab179076 - H2B K12me1

F20	19.54145625	ab179077 - H2B K12me2
F21	15.77092604	ab179078 - H2B K12me3
F22	5.333185417	ab179079 - H2B K12ac + K15ac
F23	12.49973958	ab179080 - H2B S14p
F24	10.25398438	ab179081 - H2B K15ac
F25	14.90199479	ab179082 - H2B K15cr
G1	3.688532292	ab179083 - H2B K15me1
G2	16.425675	ab179084 - H2B K15me2
G3	5.932645833	ab179085 - H2B K15me3
G4	15.46773229	ab179086 - H2B K16ac
G5	19.31280833	ab179087 - H2B K16cr
G6	17.89507708	ab179088 - H2B K20ac
G7	22.643425	ab179089 - H2B K20cr
G8	20.87832917	ab179090 - H2B K20me1
G9	21.95018542	ab179091 - H2B K20me2
G10	15.19070729	ab179092 - H2B K20me3
G11	17.17035521	ab179093 - H2B K23cr
G12	25.14335625	ab179094 - H2B K23me1
G13	19.8198625	ab179095 - H2B K23me3
G14	13.72389896	ab179096 - H2B K24ac
G15	17.0279125	ab179097 - H2B unmod(26-56)
G16	21.84835729	ab179098 - H2B K34cr
G17	15.65873646	ab179099 - H2B K34fo
G18	17.44170208	ab179100 - H2B Y37hyd
G19	16.52605313	ab179101 - H2B Y37p
G20	14.95868438	ab179102 - H2B K43me2
G21	15.34951146	ab179103 - H2B K43me3
G22	20.14119375	ab179104 - H2B K46fo
G23	15.94866771	ab179105 - H2B K46ac
G24	5.162848958	ab179106 - H2B K46me1
G25	13.64141563	ab179107 - H2B K46me2
H1	8.667015625	ab179108 - H2B K46me3
H2	14.41174063	ab179109 - H2B unmod(51-81)
H3	6.270648958	ab179110 - H2B K57me1
H4	16.86971146	ab179111 - H2B K57me2
H5	6.49823125	ab179112 - H2B K57me3
H6	12.97948542	ab179113 - H2B unmod(73-103)
H7	21.68232083	ab179114 - H2B R79me1
H8	17.8039625	ab179115 - H2B R79me2s
H9	15.10880938	ab179116 - H2B R79me2a
H10	15.36454688	ab179117 - H2B Y83hyd
H11	21.12587917	ab179118 - H2B K85me1
H12	21.16333646	ab179119 - H2B K85me2
H13	14.10923125	ab179120 - H2B K85me3
H14	20.73721875	ab179121 - H2B unmod(94-125)
H15	14.54737292	ab179122 - H2B R99me1
H16	17.07434271	ab179123 - H2B R99me2s
H17	17.32101979	ab179124 - H2B R99me2a
H18	12.21509896	ab179125 - H2B K108fo
H19	17.11183438	ab179126 - H2B K116fo
H20	18.3474625	ab179127 - H2B K116me1
H21	17.17076146	ab179128 - H2B K116me2
H22	16.76328542	ab179129 - H2B K116me3
H23	16.20693854	ab179130 - H2B K116ubiq
H24	14.58656875	ab179131 - H2B K120ac
H25	15.19741667	ab179132 - H2B K120fo
I1	20.24706771	ab179133 - H2B K120ubiq
I2	20.27470729	ab179134 - H2B K125ubiq
I3	5.568092708	ab179135 - Sch2B unmod(102-132)
I4	11.72697083	ab179136 - H3 unmod(1-33)
I5	16.43182604	ab179137 - H3 R2cit
I6	12.70269063	ab179138 - H3 R2cit + K4me2
I7	14.60106354	ab179139 - H3 R2cit + K4me3
I8	16.33393333	ab179140 - H3 R2cit + K4me3 + K9ac + K14ac + K18ac
I9	9.3185375	ab179141 - H3 R2cit + K4ac + K9ac + K14ac + K18ac
I10	10.35000625	ab179142 - H3 R2me1
I11	16.54300208	ab179143 - H3 R2me1 + K4me2
I12	11.98507708	ab179144 - H3 R2me1 + K4me3
I13	13.69334688	ab179145 - H3 R2me1 + K4me3 + K9ac + K14ac + K18ac
I14	10.88629063	ab179146 - H3 R2me2s
I15	12.88456979	ab179147 - H3 R2me2s + K4me2
I16	23.37075313	ab179148 - H3 R2me2s + K4me3
I17	14.96414792	ab179149 - H3 R2me2s + K4me3 + K9ac + K14ac + K18ac
I18	23.87962708	ab179150 - H3 R2me2a

I19	17.87041875	ab179151 - H3 R2me2a + K4me2
I20	21.95258125	ab179152 - H3 R2me2a + K4me3
I21	15.26214167	ab179153 - H3 R2me2a + K4me3 + S10p
I22	23.47180833	ab179154 - H3 R2me2a + T3p + K4me3
I23	16.30215313	ab179155 - H3 R2me2a + K4ac + K9ac + K14ac + K18ac
I24	20.47424479	ab179156 - H3 R2me2a + K4me3 + K9ac + K14ac + K18ac
I25	19.20754271	ab179157 - H3 R2me2a + T3p + K4me3 + K9ac + K14ac + K18ac
J1	17.77146979	ab179158 - H3 R2me2a + K4me3 + K9ac + S10p + K14ac + K18ac
J2	19.36751875	ab179159 - H3 T3p
J3	18.49604583	ab179160 - H3 T3p + K4me1
J4	33.07911875	ab179161 - H3 T3p + K4me2
J5	22.38529583	ab179162 - H3 T3p + K4me3 + K9ac + K14ac + K18ac
J6	11.79900313	ab179163 - H3 K4ac
J7	15.94366354	ab179164 - H3 K4ac + K9ac
J8	9.353380208	ab179165 - H3 K4ac + K14ac
J9	15.81525625	ab179166 - H3 K4ac + K18ac
J10	12.85810625	ab179167 - H3 K4ac + K9me3
J11	17.07275313	ab179168 - H3 K4ac + K9ac + K14ac
J12	14.60344896	ab179169 - H3 K4ac + K9ac + K18ac
J13	14.51418021	ab179170 - H3 K4ac + K14ac + K18ac
J14	24.06856354	ab179171 - H3 K4ac + K9ac + K14ac + K18ac
J15	20.48894167	ab179172 - H3 K4ac + K9me1 + K14ac + K18ac
J16	24.40581146	ab179173 - H3 K4ac + K9me2 + K14ac + K18ac
J17	16.06371875	ab179174 - H3 K4ac + K9me3 + K14ac + K18ac
J18	23.3877125	ab179175 - H3 K4ac + K9ac + S10p + K14ac + K18ac
J19	14.78472917	ab179176 - H3 K4cr
J20	15.37986667	ab179177 - H3 K4me1 + K9me2
J21	20.54942604	ab179178 - H3 K4me1 + K9ac + K18ac
J22	18.64648438	ab179179 - H3 K4me1 + K9ac + K14ac + K18ac
J23	17.28972917	ab179180 - H3 K4me2 + T6p
J24	22.353175	ab179181 - H3 K4me2 + K9me2
J25	21.27358229	ab179182 - H3 K4me2 + K9ac + K18ac
K1	20.70985	ab179183 - H3 K4me2 + K9ac + K14ac + K18ac
K2	8.923509375	ab179184 - H3 K4me3 + T6p
K3	13.60499583	ab179185 - H3 K4me3 + K9ac
K4	10.74910521	ab179186 - H3 K4me3 + K9me2
K5	12.93231771	ab179187 - H3 K4me3 + K9me3
K6	11.04658854	ab179188 - H3 K4me3 + S10p
K7	14.37413125	ab179189 - H3 K4me3 + K14ac
K8	12.76849167	ab179190 - H3 K4me3 + K18ac
K9	14.38451042	ab179191 - H3 K4me3 + R8me2s + K9me3
K10	21.7502375	ab179192 - H3 K4me3 + R8me2a + K9me3
K11	16.51424688	ab179193 - H3 K4me3 + K9ac + S10p
K12	24.46198021	ab179194 - H3 K4me3 + K9ac + K14ac
K13	18.07553021	ab179195 - H3 K4me3 + K9ac + K18ac
K14	21.43744479	ab179196 - H3 K4me3 + K14ac + K18ac
K15	16.99269896	ab179197 - H3 K4me3 + K9ac + K14ac + K18ac
K16	23.45019479	ab179198 - H3 K4me3 + K9ac + S10p + K14ac + K18ac
K17	15.94725729	ab179199 - H3 K4me3 + T6p + K9ac + K14ac + K18ac
K18	21.99515	ab179200 - H3 K4me2 + T6p + K9ac + K14ac + K18ac
K19	21.42329896	ab179201 - H3 K4bio + K9bio + K18bio
K20	15.6286	ab179202 - H3 T6p + K9me3
K21	14.74323646	ab179203 - H3 T6p + R8me2a + K9me3
K22	18.58858438	ab179204 - H3 T6p + K9ac + K14ac + K18ac
K23	20.59127188	ab179205 - H3 R8me1
K24	19.14697604	ab179206 - H3 R8me1 + K9me1
K25	17.22208021	ab179207 - H3 R8me1 + K9me2
L1	15.92953333	ab179208 - H3 R8me1 + K9me3
L2	17.72684479	ab179209 - H3 R8me2s
L3	15.9458875	ab179210 - H3 R8me2s + K9me1
L4	21.84871146	ab179211 - H3 R8me2s + K9me2
L5	13.77223646	ab179212 - H3 R8me2s + K9me3
L6	54.40526667	ab179213 - H3 R8me2a
L7	18.21760625	ab179214 - H3 R8me2a + K9me1
L8	32.66934167	ab179215 - H3 R8me2a + K9me2
L9	15.06487604	ab179216 - H3 R8me2a + K9me3
L10	34.00900833	ab179217 - H3 R8cit
L11	17.56400313	ab179218 - H3 K9ac + K14ac
L12	17.51638333	ab179219 - H3 K9ac + K18ac
L13	12.2594125	ab179220 - H3 K9ac + K14ac + K18ac
L14	16.54641771	ab179221 - H3 K9ac + K14ac + K18ac + K23ac + K27ac
L15	14.68312604	ab179222 - H3 K9ac + S10p
L16	22.06242396	ab179223 - H3 K9cr
L17	20.95073021	ab179224 - H3 K9me2

L18	15.4711625	ab179225 - H3 K9me2 + S10p
L19	14.22292396	ab179226 - H3 K9me2 + K27me2
L20	18.95750417	ab179227 - H3 S10p + S28p
L21	38.35166354	ab179228 - H3 S10p K14ac
L22	16.38342708	ab179229 - H3 K14me1
L23	20.00929896	ab179230 - H3 K14me2
L24	17.08181979	ab179231 - H3 K14me3
L25	18.69765	ab179232 - H3 K14ac + K18ac
M1	15.06443854	ab179233 - H3 R17cit
M2	18.31082604	ab179234 - H3 R17me1
M3	36.869225	ab179235 - H3 R17me2s
M4	20.72456771	ab179236 - H3 R17me2a + K18ac
M5	17.10256875	ab179237 - H3 K18cr
M6	18.19965833	ab179238 - H3 K18me1
M7	18.61050625	ab179239 - H3 K18me2
M8	18.37860208	ab179240 - H3 K18me3 + K36me3
M9	17.95746458	ab179241 - H3 K23cr
M10	19.66751354	ab179242 - H3 K23me1
M11	17.12608646	ab179243 - H3 K23me2
M12	18.63039271	ab179244 - H3 K23me3
M13	14.0541125	ab179245 - H3 T22p + K23me3
M14	13.65487083	ab179246 - H3 K23pr
M15	23.2968875	ab179247 - H3 unmod(20-51)
M16	19.53414688	ab179248 - H3 R26me1
M17	20.65803854	ab179249 - H3 R26me2a
M18	16.24143021	ab179250 - H3 R26me2a + K27me1
M19	22.0217875	ab179251 - H3 R26me2a + K27me3
M20	13.86917396	ab179252 - H3 K27ac + S28p
M21	19.09510417	ab179253 - H3 K27cr
M22	15.57469167	ab179254 - H3 K27me1
M23	11.067625	ab179255 - H3 K27me2 + S28p
M24	14.50596563	ab179256 - H3 K27me3
M25	12.20736563	ab179257 - H3 K27me3 + S28p
N1	16.61217813	ab179258 - H3 P30hpro
N2	15.6826875	ab179259 - H3 P30(5,5me2)
N3	15.80022292	ab179260 - H3 P30
N4	14.20159375	ab179261 - H3 P30mutA
N5	19.30293021	ab179262 - H3 T32GlcNAc
N6	16.10644167	ab179263 - H3 K36ac
N7	20.53055	ab179264 - H3 P38hpro
N8	20.97543438	ab179265 - H3 P38(5,5me2)
N9	15.11759896	ab179266 - H3 P38
N10	12.77975208	ab179267 - H3 P38mutA
N11	12.07825	ab179268 - H3 Y41p
N12	31.14108438	ab179269 - H3 unmod(50-80)
N13	18.52395	ab179270 - H3 K56ac
N14	15.48255313	ab179271 - H3 K56cr
N15	17.83547188	ab179272 - H3 K56fo
N16	20.79774271	ab179273 - H3 K56me1
N17	20.43971042	ab179274 - H3 K56me2
N18	19.47892292	ab179275 - H3 K56me3
N19	21.65484688	ab179276 - H3 R63me1
N20	18.2524	ab179277 - H3 R63me2s
N21	17.90189479	ab179278 - H3 R63me2a
N22	19.34146563	ab179279 - H3 K64ac
N23	18.91529375	ab179280 - H3 K64me2
N24	27.53280833	ab179281 - H3 K64me3
N25	19.49412396	ab179282 - H3 unmod(73-103)
O1	26.08532396	ab179283 - H3 K79ac
O2	19.39969063	ab179284 - H3 K79fo
O3	11.69658125	ab179285 - H3 T80p
O4	15.51666042	ab179286 - H3 K79me3 + T80p
O5	7294.456318	ab179287 - H3 unmod(99-135)
O6	19.81487188	ab179288 - H3 T118O-ac
O7	99.99289583	ab179289 - H3 K122fo
O8	17.19782813	ab179290 - H3 K122me1
O9	31.46023854	ab179291 - H3 R128me1
O10	62.94159792	ab179292 - H3 R128me2s
O11	25.86470313	ab179293 - H3.3 unmod(10-40)
O12	19.22044479	ab179294 - H3.3 S28p
O13	18.84905729	ab179295 - H3.3 S28p + S31p
O14	24.39017083	ab179296 - H3.3 S31p
O15	25.25104167	ab179297 - H4 unmod(1-30)
O16	33.54370104	ab179298 - H4 S1p + R3me1

O17	25.12614271	ab179299 - H4 S1p + R3me2s
O18	25.34244688	ab179300 - H4 S1p + R3me2a
O19	26.46793958	ab179301 - H4 S1p + K5ac + K8ac + K12ac + K16ac
O20	22.41920729	ab179302 - H4 R3cit
O21	19.83466458	ab179303 - H4 R3me1 + K5ac + K8ac + K12ac + K16ac + K20ac
O22	21.33342708	ab179304 - H4 R3me2s
O23	20.07791667	ab179305 - H4 R3me2s + K5ac + K8ac + K12ac + K16ac + K20ac
O24	24.74021875	ab179306 - H4 R3me2a
O25	17.67426771	ab179307 - H4 R3me2a + K5ac
P1	24.45079479	ab179308 - H4 R3me2a + K5ac + K8ac + K12ac + K16ac + K20ac
P2	19.2735375	ab179309 - H4 K5ac
P3	23.42766354	ab179310 - H4 K5ac + K8ac
P4	21.83098333	ab179311 - H4 K5ac + K12ac
P5	26.00874583	ab179312 - H4 K5ac + K16ac
P6	25.26993021	ab179313 - H4 K5ac + K8ac + K12ac
P7	31.21863958	ab179314 - H4 K5ac + K8me1 + K12ac
P8	21.81464896	ab179315 - H4 K5ac + K8ac + K12ac + K16ac
P9	24.41350417	ab179316 - H4 K5ac + K8me1 + K12ac + K16ac
P10	25.31771771	ab179317 - H4 K5ac + K8ac + K12me1 + K16ac
P11	23.56113646	ab179318 - H4 K5me1
P12	19.7825375	ab179319 - H4 K5me1 + K8ac + K12me1
P13	20.33031042	ab179320 - H4 K5me1 + K8me1 + K12me1
P14	22.4835875	ab179321 - H4 K5me1 + K8ac + K12ac + K16ac
P15	19.49476667	ab179322 - H4 K5me2
P16	21.72378646	ab179323 - H4 K5me3
P17	25.704575	ab179324 - H4 K5cr
P18	26.99735313	ab179325 - H4 K8ac + K12ac
P19	26.05740313	ab179326 - H4 K8ac + K16ac
P20	27.01656979	ab179327 - H4 K8cr
P21	22.91388021	ab179328 - H4 K8me1
P22	23.82244792	ab179329 - H4 K12ac
P23	24.66348333	ab179330 - H4 K12ac + K16ac
P24	25.69055833	ab179331 - H4 K12ac + K16ac + K20me2
P25	23.00081667	ab179332 - H4 K12ac + K16ac + K20me3
Q1	28.52007188	ab179333 - H4 K5mutQ + K8mutQ + K12ac + K16mutQ + K20mutQ
Q2	20.80489479	ab179334 - H4 K12cr
Q3	28.13640417	ab179335 - H4 K12fo
Q4	18.73851875	ab179336 - H4 K12me1
Q5	29.14561458	ab179337 - H4 K12me2
Q6	22.78453438	ab179338 - H4 K12me3
Q7	25.71602292	ab179339 - H4 K16ac
Q8	19.71225521	ab179340 - H4 K16ac + K20ac
Q9	20.82901354	ab179341 - H4 K16cr
Q10	17.24039271	ab179342 - H4 K16me1
Q11	21.14898333	ab179343 - H4 K16me3
Q12	21.58886146	ab179344 - H4 R17me1
Q13	24.67556042	ab179345 - H4 R17me2s
Q14	23.36001146	ab179346 - H4 R17me2a
Q15	34.42535208	ab179347 - H4 R19me1
Q16	20.70290729	ab179348 - H4 R19me2s
Q17	39.05425104	ab179349 - H4 R19me2a
Q18	21.01059896	ab179350 - H4 K20ac
Q19	22.13990417	ab179351 - H4 R23me1
Q20	23.52199688	ab179352 - H4 R23me2s
Q21	20.91947708	ab179353 - H4 R23me2a
Q22	20.97012604	ab179354 - H4 unmod(24-54)
Q23	22.16688958	ab179355 - H4 K31fo
Q24	24.75390521	ab179356 - H4 K31me1
Q25	19.98817292	ab179357 - H4 K31me2
R1	20.39401667	ab179358 - H4 K31me3
R2	25.53273125	ab179359 - H4 P32hyd
R3	20.73607396	ab179360 - H4 R35me1
R4	23.14279792	ab179361 - H4 R35me2s
R5	20.74459792	ab179362 - H4 R35me2a
R6	18.59754167	ab179363 - H4 S47O-ac
R7	19.58524167	ab179364 - H4 S47GlcNAc
R8	18.25934271	ab179365 - H4 S47p
R9	23.62802396	ab179366 - H4 unmod(45-75)
R10	21.63505	ab179367 - H4 Y51hyd
R11	22.51769688	ab179368 - H4 R55me1
R12	23.57286875	ab179369 - H4 R55me2s
R13	23.17443958	ab179370 - H4 R55me2a
R14	22.18090313	ab179371 - H4 K59fo
R15	24.4559	ab179372 - H4 K59me1

R16	15.07296771	ab179373 - H4 K59me2
R17	15.19520417	ab179374 - H4 R67me1
R18	14.24291563	ab179375 - H4 R67me2s
R19	17.19156354	ab179376 - H4 R67me2a
R20	14.23343958	ab179377 - H4 unmod(71-102)
R21	20.72345521	ab179378 - H4 K77me1
R22	13.81194896	ab179379 - H4 K77me2
R23	17.18807917	ab179380 - H4 K77me3
R24	18.69165938	ab179381 - H4 K79fo
R25	16.72359896	ab179382 - H4 K79me1
S1	16.66578854	ab179383 - H4 K79me3
S2	17.42959375	ab179384 - H4 T82O-ac
S3	17.93265	ab179385 - H4 Y88hyd
S4	17.10720313	ab179386 - H4 Y88nitro
S5	19.3425125	ab179387 - H4 K91fo
S6	19.24451563	ab179403 - H2A K5ac
S7	19.09841667	ab179404 - H2A K9me1
S8	22.10391875	ab18504 - H2B S32p
S9	21.74740521	ab19828 - ScH2A S129p
S10	14.14322083	ab20631 - H3 R26cit
S11	17.6352125	ab20635 - H3 T3p + K4me3
S12	18.04735521	ab21043 - H3 K37me1
S13	18.32795625	ab21998 - H1.4 K25me2
S14	14.98853958	ab22079 - H4 K59me3
S15	22.16850104	ab22080 - H3 K122me2
S16	21.20396771	ab22082 - H3 K122me3
S17	19.75944896	ab22214 - H3 K37me2
S18	16.18790208	ab22398 - H2A R3me2a
S19	22.4463125	ab22399 - H2A R3me2s
S20	18.53724583	ab23539 - H3 T22p
S21	16.45839063	ab24003 - H3 K18ac
S22	18.90644896	ab24404 - H3 K27ac
S23	17.08771667	ab24417 - H3 K37me3
S24	20.59277396	ab24444 - H3 T11p
S25	26.83653646	ab24659 - H2A K127me3
T1	15.96651146	ab24660 - H2A K125me3
T2	17.54871563	ab26307 - H3 T45p
T3	29.27548958	ab26342 - H2B K5me2
T4	21.02194375	ab26739 - H4 T30p
T5	17.66770313	ab27896 - H4 T96p
T6	16.95420521	ab27898 - H4 T80p
T7	20.71764167	ab28756 - H2A R3me1
T8	14.83300521	ab32876 - H3 R2cit + R8cit + R17cit
T9	25.93871979	ab33047 - H2A K127me2
T10	1532.668655	ab33505 - H3 T118p
T11	24.41726042	ab33594 - H3 R128me2a
T12	19.23460625	ab34466 - H3 K122ac
T13	18.52596667	ab36927 - H2A S122p
T14	20.54078958	ab4555 - H3 K79me1
T15	18.48294688	ab4556 - H3 K79me2
T16	29.89621354	ab4557 - H3 K79me3
T17	17.39063542	ab4560 - H4 K79me2
T18	20.72477708	ab46854 - H3 K18me3
T19	17.00065521	ab48359 - H3 K23ac
T20	20.05685104	ab53244 - H3 K64me1
T21	16.51535313	ab54016 - H2A K9ac
T22	16.40884688	ab5499 - H3 S28p
T23	20.54250208	ab7768 - H3 K4me2
T24	15.61700208	ab89018 - H1.4 S26p
T25	14.64879583	ab92374 - H3 K4me3 (HEX1)
U1	23.70422292	ab95444 - H4 K16me2