

Peptide array: Histone H3 (ab201456)

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

Cell Ref	AUC	Peptide
A1	5.493486458	ab10112 - H4 K91ac
A2	3.09116875	ab10121 - H2B K85ac
A3	4.411754167	ab10127 - H2A K95me2
A4	7.870710417	ab10138 - H1.4 T17p
A5	4.28805	ab10139 - H1.4 T146p
A6	5.134846875	ab113997 - H2A.X Y142p
A7	4.979896875	ab11477 - H3 S10p
A8	4.73565625	ab12946 - H2B K5me3
A9	4.92995	ab12949 - H2B K23me2
A10	4.822030208	ab12952 - H2A K127me1
A11	4.195184375	ab13211 - H2B K5me1
A12	39.04384271	ab1340 - H3 K4me1
A13	4.403760417	ab13570 - H2B K43me1
A14	5.8369	ab13835 - HA
A15	3.48214375	ab13837 - Myc
A16	4.772207292	ab140408 - ScH2B T129p
A17	5.355244792	ab14103 - H3 T6p
A18	5.470977083	ab14724 - H4 S1p
A19	5.637186458	ab14799 - H3 T32p
A20	5.277932292	ab14943 - His
A21	4.365358333	ab14952 - H2A R77me2s
A22	6.078375	ab14964 - H4 K20me2
A23	3.746480208	ab14997 - ScH2A K21ac
A24	4.285154167	ab154782 - H3 R26me2s
A25	4.606986458	ab15591 - H3 K14ac
B1	5.568301042	ab15644 - H3 K9me3 + S10p
B2	4.700397917	ab15645 - H2A.X S139p
B3	4.447158333	ab15824 - H4 K8ac
B4	4.722698958	ab15829 - V5
B5	3.961992708	ab16635 - H3 K9ac
B6	3.872335417	ab166685 - H2B S112GlcNAc
B7	11.331975	ab166688 - H2B S36GlcNAc
B8	4.202727083	ab16935 - H3 R17me2a
B9	5.795814583	ab17043 - H4 K20me1
B10	3.880564583	ab17162 - H3 P16hyd + P30hyd
B11	4.902823958	ab17567 - H4 K20me3
B12	4.940057292	ab17587 - H1.4 K25me3
B13	3.28248125	ab175905 - H2B K11ac
B14	4.337509375	ab176211 - H2B K23ac
B15	5.435314583	ab17632 - H4 K5bio
B16	3.946815625	ab1771 - H3 K9me1
B17	5.060359375	ab1773 - H3 K9me3
B18	5.227141667	ab17770 - H4 R3me1
B19	3.646457292	ab1781 - H3 K27me2
B20	4.074513542	ab1783 - H3 K36me1
B21	7.123291146	ab1784 - H3 K36me2
B22	4.701653125	ab1785 - H3 K36me3
B23	5.487545833	ab178667 - ScH2A unmod(1-30)
B24	5.118621875	ab178668 - ScH2A unmod(102-132)
B25	5.316447917	ab178670 - H2A.Z unmod(1-30)
C1	4.702711458	ab178833 - Flag
C2	3.4261875	ab178835 - H2B R99mutO
C3	6.5484625	ab178836 - H4 R23mutO
C4	4.590569792	ab178837 - H4 G48mutD
C5	4.94393125	ab178987 - H1.4 unmod(1-30)
C6	4.829029167	ab178988 - H1.4 S1p
C7	5.920653125	ab178989 - H1.4 T3p
C8	166.7449771	ab178990 - H1.4 S1p + T3p
C9	4.377714583	ab178991 - H1.4 K25me1
C10	5.214676042	ab178992 - H1.4 K25me3 + S26p
C11	3.603203125	ab178993 - H1.4 unmod(130-160)
C12	4.893855208	ab178994 - H2A unmod(1-30)
C13	4.796560417	ab178995 - H2A S1p
C14	4.518294792	ab178996 - H2A S1p + K5ac
C15	3.856551042	ab178997 - H2A S1p + R3me2a + K5ac
C16	4.852672917	ab178998 - H2A S1p + R3me2s + K5ac
C17	4.938405208	ab178999 - H2A S1p + R3cit + K5ac
C18	5.796486458	ab179000 - H2A S1ac + S1p + K5ac + K9ac + K13ac + K15ac
C19	5.055725	ab179001 - H2A S1p + K5ac + K9ac + K13ac + K15ac
C20	5.204889583	ab179002 - H2A R3cit

C21	4.424342708	ab179003 - H2A R3cit + K5ac
C22	8.406107292	ab179004 - H2A R3me2a + K5ac
C23	4.846873958	ab179005 - H2A R3me2s + K5ac
C24	7.332715625	ab179006 - H2A S1ac + K5ac + K9ac + K13ac + K15ac
C25	4.858508333	ab179007 - H2A K5ac + K9ac + K13ac + K15ac
D1	5.737051042	ab179008 - H2A K9me2
D2	4.505355208	ab179009 - H2A K9me3
D3	5.86451875	ab179010 - H2A R11me1
D4	15.61956979	ab179011 - H2A R11me2s
D5	3.997680208	ab179012 - H2A R11me2a
D6	6.033622917	ab179013 - H2A K13ac
D7	4.334461458	ab179014 - H2A K15ac
D8	5.301836458	ab179015 - H2A R17me1
D9	3.749432292	ab179016 - H2A R17me2s
D10	5.740519792	ab179017 - H2A R17me2a
D11	5.806616667	ab179018 - H2A unmod(20-50)
D12	3.870073958	ab179019 - H2A P26hyd
D13	4.612040625	ab179020 - H2A R29me1
D14	5.138959375	ab179021 - H2A R29me2s
D15	7.283972917	ab179022 - H2A R29me2a
D16	4.771646875	ab179023 - H2A K36ac
D17	4.800997917	ab179024 - H2A K36cr
D18	5.142791667	ab179025 - H2A Y39hyd
D19	4.437134375	ab179026 - H2A R42me1
D20	3.579957292	ab179027 - H2A R42me2s
D21	5.232490625	ab179028 - H2A R42me2a
D22	4.788619792	ab179029 - H2A unmod(71-103)
D23	4.272788542	ab179030 - H2A R77me1
D24	15.66313646	ab179031 - H2A R77me2a
D25	4.778519792	ab179032 - H2A R88me1
E1	3.707355208	ab179033 - H2A R88me2s
E2	5.097710417	ab179034 - H2A R88me2a
E3	4.846604167	ab179035 - H2A K95me1
E4	3.959675	ab179036 - H2A K95me3
E5	5.682376042	ab179037 - H2A unmod(92-123)
E6	6.064588542	ab179038 - H2A T101GlcNAc
E7	4.375952083	ab179039 - H2A Q104me1
E8	4.171482292	ab179040 - H2A unmod(112-129)
E9	4.636302083	ab179041 - H2A K118me1
E10	31.33834167	ab179042 - H2A K118me2
E11	4.907548958	ab179043 - H2A K118me3
E12	5.234615625	ab179044 - H2A K118fo
E13	5.744090625	ab179045 - H2A K118cr
E14	5.436297917	ab179046 - H2A K119cr
E15	3.800954167	ab179047 - H2A K118ubiq
E16	4.51835	ab179048 - H2A K119ubiq
E17	4.025846875	ab179049 - H2A T120p
E18	3.945161458	ab179050 - H2A K125me1
E19	4.022011458	ab179051 - H2A K125me2
E20	4.004666667	ab179052 - H2A K125cr
E21	17.79359583	ab179053 - H2A.X unmod(1-30)
E22	4.130816667	ab179054 - H2A.X K5ac
E23	6.41768125	ab179055 - H2A.X unmod(114-142)
E24	5.687764583	ab179056 - H2A.X K118ubiq
E25	11.23938854	ab179057 - H2A.X K119ubiq
F1	4.545922917	ab179058 - H2A.X T120p
F2	6.254165625	ab179059 - H2A.X K127ubiq
F3	5.406901042	ab179060 - H2A.X S139p + Y142p
F4	4.866903125	ab179061 - H2A.Z K4ac
F5	5.305879167	ab179062 - H2A.Z K7ac
F6	4.585678125	ab179063 - H2A.Z K11ac
F7	4.805004167	ab179064 - H2A.Z K4ac + K7ac + K11ac
F8	3.318498958	ab179065 - H2B unmod(1-30)
F9	5.543916667	ab179066 - H2B K5ac
F10	4.144410417	ab179067 - H2B K5cr
F11	5.147267708	ab179068 - H2B K5fo
F12	5.051740625	ab179069 - H2B P10hyd
F13	5.178373958	ab179070 - H2B K11me1
F14	8.268948958	ab179071 - H2B K11me2
F15	5.217955208	ab179072 - H2B K11me3
F16	6.199154167	ab179073 - H2B K11cr
F17	4.66843125	ab179074 - H2B K12ac
F18	18.64166771	ab179075 - H2B K12cr
F19	5.296659375	ab179076 - H2B K12me1

F20	10.08408229	ab179077 - H2B K12me2
F21	4.50988125	ab179078 - H2B K12me3
F22	3.686234375	ab179079 - H2B K12ac + K15ac
F23	4.754817708	ab179080 - H2B S14p
F24	4.087047917	ab179081 - H2B K15ac
F25	5.638011458	ab179082 - H2B K15cr
G1	4.394242708	ab179083 - H2B K15me1
G2	4.78945	ab179084 - H2B K15me2
G3	4.149807292	ab179085 - H2B K15me3
G4	5.987142708	ab179086 - H2B K16ac
G5	4.826245833	ab179087 - H2B K16cr
G6	5.625813542	ab179088 - H2B K20ac
G7	3.462447917	ab179089 - H2B K20cr
G8	4.596214583	ab179090 - H2B K20me1
G9	4.459789583	ab179091 - H2B K20me2
G10	4.290294792	ab179092 - H2B K20me3
G11	4.204953125	ab179093 - H2B K23cr
G12	4.203411458	ab179094 - H2B K23me1
G13	3.438314583	ab179095 - H2B K23me3
G14	4.458751042	ab179096 - H2B K24ac
G15	3.757111458	ab179097 - H2B unmod(26-56)
G16	3.562044792	ab179098 - H2B K34cr
G17	4.263085417	ab179099 - H2B K34fo
G18	4.224602083	ab179100 - H2B Y37hyd
G19	3.9946	ab179101 - H2B Y37p
G20	5.190363542	ab179102 - H2B K43me2
G21	4.759902083	ab179103 - H2B K43me3
G22	9.029605208	ab179104 - H2B K46fo
G23	5.090625	ab179105 - H2B K46ac
G24	4.505645833	ab179106 - H2B K46me1
G25	5.226629167	ab179107 - H2B K46me2
H1	4.430051042	ab179108 - H2B K46me3
H2	4.473942708	ab179109 - H2B unmod(51-81)
H3	2.74698125	ab179110 - H2B K57me1
H4	4.000857292	ab179111 - H2B K57me2
H5	3.66320625	ab179112 - H2B K57me3
H6	4.55645625	ab179113 - H2B unmod(73-103)
H7	4.008366667	ab179114 - H2B R79me1
H8	3.920065625	ab179115 - H2B R79me2s
H9	4.333511458	ab179116 - H2B R79me2a
H10	3.774955208	ab179117 - H2B Y83hyd
H11	5.692401042	ab179118 - H2B K85me1
H12	3.219955208	ab179119 - H2B K85me2
H13	4.474258333	ab179120 - H2B K85me3
H14	4.005198958	ab179121 - H2B unmod(94-125)
H15	4.999096875	ab179122 - H2B R99me1
H16	3.631609375	ab179123 - H2B R99me2s
H17	5.056128125	ab179124 - H2B R99me2a
H18	11.26286354	ab179125 - H2B K108fo
H19	3.792020833	ab179126 - H2B K116fo
H20	5.901814583	ab179127 - H2B K116me1
H21	3.669009375	ab179128 - H2B K116me2
H22	4.405757292	ab179129 - H2B K116me3
H23	4.31338125	ab179130 - H2B K116ubiq
H24	3.735	ab179131 - H2B K120ac
H25	4.421561458	ab179132 - H2B K120fo
I1	13.39465938	ab179133 - H2B K120ubiq
I2	5.852783333	ab179134 - H2B K125ubiq
I3	4.840858333	ab179135 - Sch2B unmod(102-132)
I4	5.836311458	ab179136 - H3 unmod(1-33)
I5	4.913897917	ab179137 - H3 R2cit
I6	3.530864583	ab179138 - H3 R2cit + K4me2
I7	4.256080208	ab179139 - H3 R2cit + K4me3
I8	5.623202083	ab179140 - H3 R2cit + K4me3 + K9ac + K14ac + K18ac
I9	4.714647917	ab179141 - H3 R2cit + K4ac + K9ac + K14ac + K18ac
I10	5.684685417	ab179142 - H3 R2me1
I11	5.270848958	ab179143 - H3 R2me1 + K4me2
I12	3.947002083	ab179144 - H3 R2me1 + K4me3
I13	4.453963542	ab179145 - H3 R2me1 + K4me3 + K9ac + K14ac + K18ac
I14	4.30896875	ab179146 - H3 R2me2s
I15	4.665107292	ab179147 - H3 R2me2s + K4me2
I16	4.95145	ab179148 - H3 R2me2s + K4me3
I17	3.810782292	ab179149 - H3 R2me2s + K4me3 + K9ac + K14ac + K18ac
I18	5.355119792	ab179150 - H3 R2me2a

I19	5.094579167	ab179151 - H3 R2me2a + K4me2
I20	4.124552083	ab179152 - H3 R2me2a + K4me3
I21	4.169988542	ab179153 - H3 R2me2a + K4me3 + S10p
I22	3.741142708	ab179154 - H3 R2me2a + T3p + K4me3
I23	4.432451042	ab179155 - H3 R2me2a + K4ac + K9ac + K14ac + K18ac
I24	4.344678125	ab179156 - H3 R2me2a + K4me3 + K9ac + K14ac + K18ac
I25	3.957452083	ab179157 - H3 R2me2a + T3p + K4me3 + K9ac + K14ac + K18ac
J1	3.748898958	ab179158 - H3 R2me2a + K4me3 + K9ac + S10p + K14ac + K18ac
J2	4.782326042	ab179159 - H3 T3p
J3	4.806684375	ab179160 - H3 T3p + K4me1
J4	3.625136458	ab179161 - H3 T3p + K4me2
J5	4.21404375	ab179162 - H3 T3p + K4me3 + K9ac + K14ac + K18ac
J6	4.19260625	ab179163 - H3 K4ac
J7	4.842439583	ab179164 - H3 K4ac + K9ac
J8	4.556796875	ab179165 - H3 K4ac + K14ac
J9	3.603660417	ab179166 - H3 K4ac + K18ac
J10	3.595873958	ab179167 - H3 K4ac + K9me3
J11	4.507461458	ab179168 - H3 K4ac + K9ac + K14ac
J12	6.104426042	ab179169 - H3 K4ac + K9ac + K18ac
J13	4.46469375	ab179170 - H3 K4ac + K14ac + K18ac
J14	7.7951875	ab179171 - H3 K4ac + K9ac + K14ac + K18ac
J15	4.642476042	ab179172 - H3 K4ac + K9me1 + K14ac + K18ac
J16	7.70271875	ab179173 - H3 K4ac + K9me2 + K14ac + K18ac
J17	5.320053125	ab179174 - H3 K4ac + K9me3 + K14ac + K18ac
J18	4.767076042	ab179175 - H3 K4ac + K9ac + S10p + K14ac + K18ac
J19	4.319457292	ab179176 - H3 K4cr
J20	4.551664583	ab179177 - H3 K4me1 + K9me2
J21	4.6762875	ab179178 - H3 K4me1 + K9ac + K18ac
J22	4.3859625	ab179179 - H3 K4me1 + K9ac + K14ac + K18ac
J23	3.690841667	ab179180 - H3 K4me2 + T6p
J24	5.614601042	ab179181 - H3 K4me2 + K9me2
J25	4.32913125	ab179182 - H3 K4me2 + K9ac + K18ac
K1	4.352205208	ab179183 - H3 K4me2 + K9ac + K14ac + K18ac
K2	4.56004375	ab179184 - H3 K4me3 + T6p
K3	5.829042708	ab179185 - H3 K4me3 + K9ac
K4	5.033071875	ab179186 - H3 K4me3 + K9me2
K5	4.675028125	ab179187 - H3 K4me3 + K9me3
K6	3.45361875	ab179188 - H3 K4me3 + S10p
K7	6.823427083	ab179189 - H3 K4me3 + K14ac
K8	5.038086458	ab179190 - H3 K4me3 + K18ac
K9	5.540686458	ab179191 - H3 K4me3 + R8me2s + K9me3
K10	4.505684375	ab179192 - H3 K4me3 + R8me2a + K9me3
K11	5.057108333	ab179193 - H3 K4me3 + K9ac + S10p
K12	4.586077083	ab179194 - H3 K4me3 + K9ac + K14ac
K13	4.790567708	ab179195 - H3 K4me3 + K9ac + K18ac
K14	9.599588542	ab179196 - H3 K4me3 + K14ac + K18ac
K15	4.282758333	ab179197 - H3 K4me3 + K9ac + K14ac + K18ac
K16	3.758469792	ab179198 - H3 K4me3 + K9ac + S10p + K14ac + K18ac
K17	3.910610417	ab179199 - H3 K4me3 + T6p + K9ac + K14ac + K18ac
K18	4.596011458	ab179200 - H3 K4me2 + T6p + K9ac + K14ac + K18ac
K19	8.546851042	ab179201 - H3 K4bio + K9bio + K18bio
K20	4.319271875	ab179202 - H3 T6p + K9me3
K21	4.648663542	ab179203 - H3 T6p + R8me2a + K9me3
K22	3.66279375	ab179204 - H3 T6p + K9ac + K14ac + K18ac
K23	4.699578125	ab179205 - H3 R8me1
K24	4.650379167	ab179206 - H3 R8me1 + K9me1
K25	4.565714583	ab179207 - H3 R8me1 + K9me2
L1	6.892892708	ab179208 - H3 R8me1 + K9me3
L2	6.130442708	ab179209 - H3 R8me2s
L3	3.979003125	ab179210 - H3 R8me2s + K9me1
L4	5.796265625	ab179211 - H3 R8me2s + K9me2
L5	4.076985417	ab179212 - H3 R8me2s + K9me3
L6	4.519873958	ab179213 - H3 R8me2a
L7	4.452696875	ab179214 - H3 R8me2a + K9me1
L8	4.195445833	ab179215 - H3 R8me2a + K9me2
L9	4.446775	ab179216 - H3 R8me2a + K9me3
L10	4.389975	ab179217 - H3 R8cit
L11	7.125527083	ab179218 - H3 K9ac + K14ac
L12	7.444939583	ab179219 - H3 K9ac + K18ac
L13	4.7867125	ab179220 - H3 K9ac + K14ac + K18ac
L14	3.919621875	ab179221 - H3 K9ac + K14ac + K18ac + K23ac + K27ac
L15	4.516604167	ab179222 - H3 K9ac + S10p
L16	4.627867708	ab179223 - H3 K9cr
L17	3.882326042	ab179224 - H3 K9me2

L18	5.267458333	ab179225 - H3 K9me2 + S10p
L19	4.833641667	ab179226 - H3 K9me2 + K27me2
L20	4.2953875	ab179227 - H3 S10p + S28p
L21	4.820297917	ab179228 - H3 S10p K14ac
L22	4.600664583	ab179229 - H3 K14me1
L23	3.848441667	ab179230 - H3 K14me2
L24	3.822977083	ab179231 - H3 K14me3
L25	4.211473958	ab179232 - H3 K14ac + K18ac
M1	3.641727083	ab179233 - H3 R17cit
M2	3.446220833	ab179234 - H3 R17me1
M3	3.4201625	ab179235 - H3 R17me2s
M4	3.14210625	ab179236 - H3 R17me2a + K18ac
M5	3.980383333	ab179237 - H3 K18cr
M6	3.276890625	ab179238 - H3 K18me1
M7	4.128286458	ab179239 - H3 K18me2
M8	9.339410417	ab179240 - H3 K18me3 + K36me3
M9	4.1474	ab179241 - H3 K23cr
M10	5.402588542	ab179242 - H3 K23me1
M11	4.739972917	ab179243 - H3 K23me2
M12	5.132514583	ab179244 - H3 K23me3
M13	4.914517708	ab179245 - H3 T22p + K23me3
M14	4.503276042	ab179246 - H3 K23pr
M15	9699.831522	ab179247 - H3 unmod(20-51)
M16	2085.448985	ab179248 - H3 R26me1
M17	152.578349	ab179249 - H3 R26me2a
M18	4.844109375	ab179250 - H3 R26me2a + K27me1
M19	50.28819948	ab179251 - H3 R26me2a + K27me3
M20	4.3691375	ab179252 - H3 K27ac + S28p
M21	4.682383333	ab179253 - H3 K27cr
M22	4.522401042	ab179254 - H3 K27me1
M23	4.099494792	ab179255 - H3 K27me2 + S28p
M24	4.85624375	ab179256 - H3 K27me3
M25	3.965332292	ab179257 - H3 K27me3 + S28p
N1	4.236142708	ab179258 - H3 P30hpro
N2	4.949017708	ab179259 - H3 P30(5,5me2)
N3	3.221579167	ab179260 - H3 P30
N4	5.564767708	ab179261 - H3 P30mutA
N5	4.116419792	ab179262 - H3 T32GlcNAc
N6	4.1898125	ab179263 - H3 K36ac
N7	4.158334375	ab179264 - H3 P38hpro
N8	19.21399167	ab179265 - H3 P38(5,5me2)
N9	2.9658	ab179266 - H3 P38
N10	4.16008125	ab179267 - H3 P38mutA
N11	3828.40522	ab179268 - H3 Y41p
N12	206.6614885	ab179269 - H3 unmod(50-80)
N13	4.362891667	ab179270 - H3 K56ac
N14	5.834317708	ab179271 - H3 K56cr
N15	4.402996875	ab179272 - H3 K56fo
N16	4.872592708	ab179273 - H3 K56me1
N17	6.0083625	ab179274 - H3 K56me2
N18	4.402691667	ab179275 - H3 K56me3
N19	5.19895625	ab179276 - H3 R63me1
N20	5.449277083	ab179277 - H3 R63me2s
N21	4.368733333	ab179278 - H3 R63me2a
N22	5.109339583	ab179279 - H3 K64ac
N23	4.80026875	ab179280 - H3 K64me2
N24	22.74338958	ab179281 - H3 K64me3
N25	6.009494792	ab179282 - H3 unmod(73-103)
O1	8.220475	ab179283 - H3 K79ac
O2	4.827672917	ab179284 - H3 K79fo
O3	5.081889583	ab179285 - H3 T80p
O4	7.769366667	ab179286 - H3 K79me3 + T80p
O5	5.596294792	ab179287 - H3 unmod(99-135)
O6	5.488261458	ab179288 - H3 T118O-ac
O7	5.514159375	ab179289 - H3 K122fo
O8	5.839827083	ab179290 - H3 K122me1
O9	7.307475	ab179291 - H3 R128me1
O10	5.825530208	ab179292 - H3 R128me2s
O11	5.741666667	ab179293 - H3.3 unmod(10-40)
O12	6.466211458	ab179294 - H3.3 S28p
O13	4.685883333	ab179295 - H3.3 S28p + S31p
O14	4.95510625	ab179296 - H3.3 S31p
O15	154.1007781	ab179297 - H4 unmod(1-30)
O16	8.145046875	ab179298 - H4 S1p + R3me1

O17	3.897148958	ab179299 - H4 S1p + R3me2s
O18	6.110461458	ab179300 - H4 S1p + R3me2a
O19	6.068317708	ab179301 - H4 S1p + K5ac + K8ac + K12ac + K16ac
O20	5.4872875	ab179302 - H4 R3cit
O21	16.295225	ab179303 - H4 R3me1 + K5ac + K8ac + K12ac + K16ac + K20ac
O22	5.203275	ab179304 - H4 R3me2s
O23	13.20512188	ab179305 - H4 R3me2s + K5ac + K8ac + K12ac + K16ac + K20ac
O24	6.047257292	ab179306 - H4 R3me2a
O25	7.570733333	ab179307 - H4 R3me2a + K5ac
P1	28.88804792	ab179308 - H4 R3me2a + K5ac + K8ac + K12ac + K16ac + K20ac
P2	5.870608333	ab179309 - H4 K5ac
P3	6.738963542	ab179310 - H4 K5ac + K8ac
P4	5.042479167	ab179311 - H4 K5ac + K12ac
P5	6.2656125	ab179312 - H4 K5ac + K16ac
P6	5.077871875	ab179313 - H4 K5ac + K8ac + K12ac
P7	15.64317396	ab179314 - H4 K5ac + K8me1 + K12ac
P8	5.361719792	ab179315 - H4 K5ac + K8ac + K12ac + K16ac
P9	20.85999167	ab179316 - H4 K5ac + K8me1 + K12ac + K16ac
P10	7.005240625	ab179317 - H4 K5ac + K8ac + K12me1 + K16ac
P11	5.135297917	ab179318 - H4 K5me1
P12	6.125588542	ab179319 - H4 K5me1 + K8ac + K12me1
P13	5.86010625	ab179320 - H4 K5me1 + K8me1 + K12me1
P14	5.653380208	ab179321 - H4 K5me1 + K8ac + K12ac + K16ac
P15	4.266404167	ab179322 - H4 K5me2
P16	5.356645833	ab179323 - H4 K5me3
P17	5.423564583	ab179324 - H4 K5cr
P18	5.045926042	ab179325 - H4 K8ac + K12ac
P19	5.625991667	ab179326 - H4 K8ac + K16ac
P20	21.39654792	ab179327 - H4 K8cr
P21	5.225913542	ab179328 - H4 K8me1
P22	6.134482292	ab179329 - H4 K12ac
P23	3.95236875	ab179330 - H4 K12ac + K16ac
P24	7.158673958	ab179331 - H4 K12ac + K16ac + K20me2
P25	5.906417708	ab179332 - H4 K12ac + K16ac + K20me3
Q1	5.903084375	ab179333 - H4 K5mutQ + K8mutQ + K12ac + K16mutQ + K20mutQ
Q2	5.9464375	ab179334 - H4 K12cr
Q3	4.808254167	ab179335 - H4 K12fo
Q4	5.033355208	ab179336 - H4 K12me1
Q5	3.7494375	ab179337 - H4 K12me2
Q6	4.445442708	ab179338 - H4 K12me3
Q7	4.766166667	ab179339 - H4 K16ac
Q8	5.213784375	ab179340 - H4 K16ac + K20ac
Q9	5.588742708	ab179341 - H4 K16cr
Q10	4.477770833	ab179342 - H4 K16me1
Q11	4.843238542	ab179343 - H4 K16me3
Q12	19.60636979	ab179344 - H4 R17me1
Q13	5.138527083	ab179345 - H4 R17me2s
Q14	14.64092604	ab179346 - H4 R17me2a
Q15	4.959780208	ab179347 - H4 R19me1
Q16	4.63030625	ab179348 - H4 R19me2s
Q17	5.059202083	ab179349 - H4 R19me2a
Q18	7.749875	ab179350 - H4 K20ac
Q19	6.066053125	ab179351 - H4 R23me1
Q20	4.63585625	ab179352 - H4 R23me2s
Q21	5.265192708	ab179353 - H4 R23me2a
Q22	6.016554167	ab179354 - H4 unmod(24-54)
Q23	6.041690625	ab179355 - H4 K31fo
Q24	5.319020833	ab179356 - H4 K31me1
Q25	5.911982292	ab179357 - H4 K31me2
R1	5.426577083	ab179358 - H4 K31me3
R2	6.228321875	ab179359 - H4 P32hyd
R3	6.974530208	ab179360 - H4 R35me1
R4	6.482509375	ab179361 - H4 R35me2s
R5	7.341641667	ab179362 - H4 R35me2a
R6	5.171497917	ab179363 - H4 S47O-ac
R7	5.967733333	ab179364 - H4 S47GlcNAc
R8	6.77458125	ab179365 - H4 S47p
R9	4.629059375	ab179366 - H4 unmod(45-75)
R10	5.467055208	ab179367 - H4 Y51hyd
R11	4.920701042	ab179368 - H4 R55me1
R12	4.740610417	ab179369 - H4 R55me2s
R13	5.908698958	ab179370 - H4 R55me2a
R14	4.874019792	ab179371 - H4 K59fo
R15	5.109825	ab179372 - H4 K59me1

R16	5.739171875	ab179373 - H4 K59me2
R17	4.875905208	ab179374 - H4 R67me1
R18	4.37353125	ab179375 - H4 R67me2s
R19	4.55230625	ab179376 - H4 R67me2a
R20	5.130166667	ab179377 - H4 unmod(71-102)
R21	4.783795833	ab179378 - H4 K77me1
R22	4.190134375	ab179379 - H4 K77me2
R23	5.213102083	ab179380 - H4 K77me3
R24	4.68178125	ab179381 - H4 K79fo
R25	6.206533333	ab179382 - H4 K79me1
S1	5.270588542	ab179383 - H4 K79me3
S2	6.111347917	ab179384 - H4 T82O-ac
S3	5.637326042	ab179385 - H4 Y88hyd
S4	5.188917708	ab179386 - H4 Y88nitro
S5	6.2909875	ab179387 - H4 K91fo
S6	4.988196875	ab179403 - H2A K5ac
S7	3.937992708	ab179404 - H2A K9me1
S8	4.29665625	ab18504 - H2B S32p
S9	5.490646875	ab19828 - ScH2A S129p
S10	3.429958333	ab20631 - H3 R26cit
S11	4.346139583	ab20635 - H3 T3p + K4me3
S12	12.30140313	ab21043 - H3 K37me1
S13	4.07839375	ab21998 - H1.4 K25me2
S14	4.766745833	ab22079 - H4 K59me3
S15	6.45634375	ab22080 - H3 K122me2
S16	6.0965	ab22082 - H3 K122me3
S17	4.830903125	ab22214 - H3 K37me2
S18	5.439841667	ab22398 - H2A R3me2a
S19	4.167034375	ab22399 - H2A R3me2s
S20	3.671496875	ab23539 - H3 T22p
S21	5.102877083	ab24003 - H3 K18ac
S22	5.451732292	ab24404 - H3 K27ac
S23	4.501673958	ab24417 - H3 K37me3
S24	4.617534375	ab24444 - H3 T11p
S25	8.986279167	ab24659 - H2A K127me3
T1	4.426740625	ab24660 - H2A K125me3
T2	32.34146146	ab26307 - H3 T45p
T3	5.064776042	ab26342 - H2B K5me2
T4	5.735022917	ab26739 - H4 T30p
T5	4.409245833	ab27896 - H4 T96p
T6	5.371096875	ab27898 - H4 T80p
T7	4.9097	ab28756 - H2A R3me1
T8	8.30980625	ab32876 - H3 R2cit + R8cit + R17cit
T9	3.384840625	ab33047 - H2A K127me2
T10	5.4124	ab33505 - H3 T118p
T11	6.370182292	ab33594 - H3 R128me2a
T12	4.832741667	ab34466 - H3 K122ac
T13	5.060279167	ab36927 - H2A S122p
T14	5.634764583	ab4555 - H3 K79me1
T15	4.403147917	ab4556 - H3 K79me2
T16	5.530223958	ab4557 - H3 K79me3
T17	5.968382292	ab4560 - H4 K79me2
T18	4.276897917	ab46854 - H3 K18me3
T19	6.293359375	ab48359 - H3 K23ac
T20	5.625820833	ab53244 - H3 K64me1
T21	5.521219792	ab54016 - H2A K9ac
T22	4.332626042	ab5499 - H3 S28p
T23	6.292228125	ab7768 - H3 K4me2
T24	5.561688542	ab89018 - H1.4 S26p
T25	5.288984375	ab92374 - H3 K4me3 (HEX1)
U1	5.137195833	ab95444 - H4 K16me2