

Peptide array: Histone H3.1 (propionylated K23)(ab317819)

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

Cell Ref	AUC	Peptide
A1	24.74731354	ab10112 - H4 K91ac
A2	18.10313333	ab10121 - H2B K85ac
A3	20.02137708	ab10127 - H2A K95me2
A4	19.62795208	ab10138 - H1.4 T17p
A5	22.54920729	ab10139 - H1.4 T146p
A6	22.49718333	ab113997 - H2A.X Y142p
A7	18.56032083	ab11477 - H3 S10p
A8	20.44414167	ab12946 - H2B K5me3
A9	18.60741771	ab12949 - H2B K23me2
A10	25.79175521	ab12952 - H2A K127me1
A11	19.98147083	ab13211 - H2B K5me1
A12	23.76399375	ab1340 - H3 K4me1
A13	24.64175625	ab13570 - H2B K43me1
A14	24.15951771	ab13835 - HA
A15	24.93963646	ab13837 - Myc
A16	16.10700625	ab140408 - ScH2B T129p
A17	22.83591354	ab14103 - H3 T6p
A18	22.65382917	ab14724 - H4 S1p
A19	21.13558333	ab14799 - H3 T32p
A20	24.25286354	ab14943 - His
A21	27.56504896	ab14952 - H2A R77me2s
A22	24.06490938	ab14964 - H4 K20me2
A23	32.07716979	ab14997 - ScH2A K21ac
A24	25.66637813	ab154782 - H3 R26me2s
A25	19.88375729	ab15591 - H3 K14ac
B1	21.90684063	ab15644 - H3 K9me3 + S10p
B2	16.14805313	ab15645 - H2A.X S139p
B3	25.82691563	ab15824 - H4 K8ac
B4	18.73189375	ab15829 - V5
B5	20.85559896	ab16635 - H3 K9ac
B6	20.72395521	ab166685 - H2B S112GlcNAc
B7	25.62664792	ab166688 - H2B S36GlcNAc
B8	26.02304167	ab16935 - H3 R17me2a
B9	26.31859063	ab17043 - H4 K20me1
B10	25.61375521	ab17162 - H3 P16hyd + P30hyd
B11	27.86468958	ab17567 - H4 K20me3
B12	12.69246667	ab17587 - H1.4 K25me3
B13	23.94436458	ab175905 - H2B K11ac
B14	33.04541875	ab176211 - H2B K23ac
B15	26.37567188	ab17632 - H4 K5bio
B16	20.53338854	ab1771 - H3 K9me1
B17	20.52651875	ab1773 - H3 K9me3
B18	25.83009271	ab17770 - H4 R3me1
B19	16.81338021	ab1781 - H3 K27me2
B20	23.74642083	ab1783 - H3 K36me1
B21	22.78486667	ab1784 - H3 K36me2
B22	24.55368854	ab1785 - H3 K36me3
B23	31.62012083	ab178667 - ScH2A unmod(1-30)
B24	29.08001458	ab178668 - ScH2A unmod(102-132)
B25	31.81396771	ab178670 - H2A.Z unmod(1-30)
C1	23.57341354	ab178833 - Flag
C2	37.8244125	ab178835 - H2B R99mutO
C3	22.69976667	ab178836 - H4 R23mutO
C4	22.08061042	ab178837 - H4 G48mutD
C5	21.75774479	ab178987 - H1.4 unmod(1-30)
C6	17.90485833	ab178988 - H1.4 S1p
C7	21.948	ab178989 - H1.4 T3p
C8	23.68218646	ab178990 - H1.4 S1p + T3p
C9	16.79098021	ab178991 - H1.4 K25me1
C10	19.94795833	ab178992 - H1.4 K25me3 + S26p
C11	19.29329479	ab178993 - H1.4 unmod(130-160)
C12	24.28075521	ab178994 - H2A unmod(1-30)
C13	21.200425	ab178995 - H2A S1p
C14	20.00095208	ab178996 - H2A S1p + K5ac
C15	22.13940938	ab178997 - H2A S1p + R3me2a + K5ac
C16	25.24446354	ab178998 - H2A S1p + R3me2s + K5ac
C17	21.43088854	ab178999 - H2A S1p + R3cit + K5ac
C18	26.66969896	ab179000 - H2A S1ac + S1p + K5ac + K9ac + K13ac + K15ac
C19	22.88369896	ab179001 - H2A S1p + K5ac + K9ac + K13ac + K15ac
C20	23.653925	ab179002 - H2A R3cit

C21	20.31564792	ab179003 - H2A R3cit + K5ac
C22	18.77445938	ab179004 - H2A R3me2a + K5ac
C23	23.4029625	ab179005 - H2A R3me2s + K5ac
C24	42.35041667	ab179006 - H2A S1ac + K5ac + K9ac + K13ac + K15ac
C25	22.25352396	ab179007 - H2A K5ac + K9ac + K13ac + K15ac
D1	19.4463625	ab179008 - H2A K9me2
D2	17.79152604	ab179009 - H2A K9me3
D3	21.56604583	ab179010 - H2A R11me1
D4	18.97408854	ab179011 - H2A R11me2s
D5	20.54887708	ab179012 - H2A R11me2a
D6	18.32688646	ab179013 - H2A K13ac
D7	24.52601042	ab179014 - H2A K15ac
D8	16.58044583	ab179015 - H2A R17me1
D9	19.26145417	ab179016 - H2A R17me2s
D10	23.85593854	ab179017 - H2A R17me2a
D11	27.93339688	ab179018 - H2A unmod(20-50)
D12	22.02971771	ab179019 - H2A P26hyd
D13	21.83124063	ab179020 - H2A R29me1
D14	22.18865208	ab179021 - H2A R29me2s
D15	21.31516146	ab179022 - H2A R29me2a
D16	37.24837813	ab179023 - H2A K36ac
D17	25.75786563	ab179024 - H2A K36cr
D18	25.04604792	ab179025 - H2A Y39hyd
D19	23.49409167	ab179026 - H2A R42me1
D20	22.29272188	ab179027 - H2A R42me2s
D21	25.38332813	ab179028 - H2A R42me2a
D22	22.11425104	ab179029 - H2A unmod(71-103)
D23	26.79633438	ab179030 - H2A R77me1
D24	23.51940104	ab179031 - H2A R77me2a
D25	20.49379479	ab179032 - H2A R88me1
E1	24.52322813	ab179033 - H2A R88me2s
E2	19.85310938	ab179034 - H2A R88me2a
E3	21.09610313	ab179035 - H2A K95me1
E4	23.50112083	ab179036 - H2A K95me3
E5	20.91111458	ab179037 - H2A unmod(92-123)
E6	18.766275	ab179038 - H2A T101GlcNAc
E7	19.04421563	ab179039 - H2A Q104me1
E8	22.15065521	ab179040 - H2A unmod(112-129)
E9	18.80947188	ab179041 - H2A K118me1
E10	20.92283438	ab179042 - H2A K118me2
E11	21.19515417	ab179043 - H2A K118me3
E12	23.65867292	ab179044 - H2A K118fo
E13	30.29274896	ab179045 - H2A K118cr
E14	60.28169271	ab179046 - H2A K119cr
E15	28.93139479	ab179047 - H2A K118ubiq
E16	25.91303542	ab179048 - H2A K119ubiq
E17	24.7890875	ab179049 - H2A T120p
E18	20.68241354	ab179050 - H2A K125me1
E19	20.74842604	ab179051 - H2A K125me2
E20	26.34129896	ab179052 - H2A K125cr
E21	26.78521042	ab179053 - H2A.X unmod(1-30)
E22	23.95798333	ab179054 - H2A.X K5ac
E23	60.15373958	ab179055 - H2A.X unmod(114-142)
E24	22.18960625	ab179056 - H2A.X K118ubiq
E25	23.90244583	ab179057 - H2A.X K119ubiq
F1	19.35651875	ab179058 - H2A.X T120p
F2	23.25291979	ab179059 - H2A.X K127ubiq
F3	18.09454375	ab179060 - H2A.X S139p + Y142p
F4	19.10645104	ab179061 - H2A.Z K4ac
F5	21.84841667	ab179062 - H2A.Z K7ac
F6	17.89927396	ab179063 - H2A.Z K11ac
F7	22.60210938	ab179064 - H2A.Z K4ac + K7ac + K11ac
F8	19.60356146	ab179065 - H2B unmod(1-30)
F9	18.65538333	ab179066 - H2B K5ac
F10	21.4774625	ab179067 - H2B K5cr
F11	23.11430417	ab179068 - H2B K5fo
F12	16.74134792	ab179069 - H2B P10hyd
F13	21.93329375	ab179070 - H2B K11me1
F14	25.60233854	ab179071 - H2B K11me2
F15	22.506125	ab179072 - H2B K11me3
F16	25.40626979	ab179073 - H2B K11cr
F17	29.87307917	ab179074 - H2B K12ac
F18	23.92773229	ab179075 - H2B K12cr
F19	24.11571458	ab179076 - H2B K12me1

F20	24.29875833	ab179077 - H2B K12me2
F21	15.01433646	ab179078 - H2B K12me3
F22	18.55505938	ab179079 - H2B K12ac + K15ac
F23	15.52920521	ab179080 - H2B S14p
F24	16.79474167	ab179081 - H2B K15ac
F25	19.11358646	ab179082 - H2B K15cr
G1	13.56846354	ab179083 - H2B K15me1
G2	17.02549063	ab179084 - H2B K15me2
G3	12.50498542	ab179085 - H2B K15me3
G4	20.29836667	ab179086 - H2B K16ac
G5	17.72370313	ab179087 - H2B K16cr
G6	32.63554375	ab179088 - H2B K20ac
G7	21.64155729	ab179089 - H2B K20cr
G8	23.2565375	ab179090 - H2B K20me1
G9	19.52674063	ab179091 - H2B K20me2
G10	22.49229479	ab179092 - H2B K20me3
G11	16.15301563	ab179093 - H2B K23cr
G12	17.72941979	ab179094 - H2B K23me1
G13	18.93037396	ab179095 - H2B K23me3
G14	129.1896156	ab179096 - H2B K24ac
G15	27.22757083	ab179097 - H2B unmod(26-56)
G16	22.82840833	ab179098 - H2B K34cr
G17	23.64818958	ab179099 - H2B K34fo
G18	25.56326146	ab179100 - H2B Y37hyd
G19	26.91746563	ab179101 - H2B Y37p
G20	21.82132396	ab179102 - H2B K43me2
G21	22.88804896	ab179103 - H2B K43me3
G22	22.47236979	ab179104 - H2B K46fo
G23	18.76775417	ab179105 - H2B K46ac
G24	14.68322917	ab179106 - H2B K46me1
G25	18.13805	ab179107 - H2B K46me2
H1	19.08370104	ab179108 - H2B K46me3
H2	19.06969792	ab179109 - H2B unmod(51-81)
H3	14.80126146	ab179110 - H2B K57me1
H4	14.23681458	ab179111 - H2B K57me2
H5	15.11094479	ab179112 - H2B K57me3
H6	22.34394688	ab179113 - H2B unmod(73-103)
H7	18.22788438	ab179114 - H2B R79me1
H8	22.61712604	ab179115 - H2B R79me2s
H9	19.75403438	ab179116 - H2B R79me2a
H10	21.31984688	ab179117 - H2B Y83hyd
H11	18.8655875	ab179118 - H2B K85me1
H12	20.37977813	ab179119 - H2B K85me2
H13	20.710025	ab179120 - H2B K85me3
H14	23.01325104	ab179121 - H2B unmod(94-125)
H15	21.63536042	ab179122 - H2B R99me1
H16	22.23507396	ab179123 - H2B R99me2s
H17	18.0337	ab179124 - H2B R99me2a
H18	17.20156146	ab179125 - H2B K108fo
H19	21.38081875	ab179126 - H2B K116fo
H20	24.76167292	ab179127 - H2B K116me1
H21	23.86616667	ab179128 - H2B K116me2
H22	19.18543958	ab179129 - H2B K116me3
H23	21.15273333	ab179130 - H2B K116ubiq
H24	23.63458542	ab179131 - H2B K120ac
H25	23.20960104	ab179132 - H2B K120fo
I1	20.68964167	ab179133 - H2B K120ubiq
I2	16.369425	ab179134 - H2B K125ubiq
I3	13.51858854	ab179135 - Sch2B unmod(102-132)
I4	12.46314271	ab179136 - H3 unmod(1-33)
I5	22.87006354	ab179137 - H3 R2cit
I6	13.79164167	ab179138 - H3 R2cit + K4me2
I7	23.06951667	ab179139 - H3 R2cit + K4me3
I8	15.57678854	ab179140 - H3 R2cit + K4me3 + K9ac + K14ac + K18ac
I9	28.35521146	ab179141 - H3 R2cit + K4ac + K9ac + K14ac + K18ac
I10	20.22833333	ab179142 - H3 R2me1
I11	26.23811354	ab179143 - H3 R2me1 + K4me2
I12	19.54528542	ab179144 - H3 R2me1 + K4me3
I13	22.28904479	ab179145 - H3 R2me1 + K4me3 + K9ac + K14ac + K18ac
I14	20.88748646	ab179146 - H3 R2me2s
I15	23.69982813	ab179147 - H3 R2me2s + K4me2
I16	20.25316667	ab179148 - H3 R2me2s + K4me3
I17	25.38051979	ab179149 - H3 R2me2s + K4me3 + K9ac + K14ac + K18ac
I18	20.65625521	ab179150 - H3 R2me2a

I19	26.28817604	ab179151 - H3 R2me2a + K4me2
I20	21.86163125	ab179152 - H3 R2me2a + K4me3
I21	22.07755104	ab179153 - H3 R2me2a + K4me3 + S10p
I22	20.88250938	ab179154 - H3 R2me2a + T3p + K4me3
I23	23.93912708	ab179155 - H3 R2me2a + K4ac + K9ac + K14ac + K18ac
I24	24.92980417	ab179156 - H3 R2me2a + K4me3 + K9ac + K14ac + K18ac
I25	20.22288229	ab179157 - H3 R2me2a + T3p + K4me3 + K9ac + K14ac + K18ac
J1	23.39501771	ab179158 - H3 R2me2a + K4me3 + K9ac + S10p + K14ac + K18ac
J2	20.573825	ab179159 - H3 T3p
J3	21.75867708	ab179160 - H3 T3p + K4me1
J4	20.01408021	ab179161 - H3 T3p + K4me2
J5	20.39215938	ab179162 - H3 T3p + K4me3 + K9ac + K14ac + K18ac
J6	17.65583021	ab179163 - H3 K4ac
J7	26.18844479	ab179164 - H3 K4ac + K9ac
J8	21.34575313	ab179165 - H3 K4ac + K14ac
J9	37.39010625	ab179166 - H3 K4ac + K18ac
J10	17.02824063	ab179167 - H3 K4ac + K9me3
J11	24.647075	ab179168 - H3 K4ac + K9ac + K14ac
J12	30.10731146	ab179169 - H3 K4ac + K9ac + K18ac
J13	32.12801667	ab179170 - H3 K4ac + K14ac + K18ac
J14	30.14229479	ab179171 - H3 K4ac + K9ac + K14ac + K18ac
J15	26.99219792	ab179172 - H3 K4ac + K9me1 + K14ac + K18ac
J16	29.34163854	ab179173 - H3 K4ac + K9me2 + K14ac + K18ac
J17	26.7299375	ab179174 - H3 K4ac + K9me3 + K14ac + K18ac
J18	34.82347396	ab179175 - H3 K4ac + K9ac + S10p + K14ac + K18ac
J19	25.01373333	ab179176 - H3 K4cr
J20	25.16974688	ab179177 - H3 K4me1 + K9me2
J21	35.93717917	ab179178 - H3 K4me1 + K9ac + K18ac
J22	24.50626042	ab179179 - H3 K4me1 + K9ac + K14ac + K18ac
J23	25.18460833	ab179180 - H3 K4me2 + T6p
J24	21.48086146	ab179181 - H3 K4me2 + K9me2
J25	24.44561146	ab179182 - H3 K4me2 + K9ac + K18ac
K1	25.2032125	ab179183 - H3 K4me2 + K9ac + K14ac + K18ac
K2	16.73085208	ab179184 - H3 K4me3 + T6p
K3	24.80095208	ab179185 - H3 K4me3 + K9ac
K4	17.45185313	ab179186 - H3 K4me3 + K9me2
K5	24.29274583	ab179187 - H3 K4me3 + K9me3
K6	21.52381979	ab179188 - H3 K4me3 + S10p
K7	24.42841979	ab179189 - H3 K4me3 + K14ac
K8	19.41012188	ab179190 - H3 K4me3 + K18ac
K9	21.36391979	ab179191 - H3 K4me3 + R8me2s + K9me3
K10	20.37905208	ab179192 - H3 K4me3 + R8me2a + K9me3
K11	25.05493542	ab179193 - H3 K4me3 + K9ac + S10p
K12	22.49230729	ab179194 - H3 K4me3 + K9ac + K14ac
K13	27.8788125	ab179195 - H3 K4me3 + K9ac + K18ac
K14	28.72349792	ab179196 - H3 K4me3 + K14ac + K18ac
K15	26.40499792	ab179197 - H3 K4me3 + K9ac + K14ac + K18ac
K16	24.59343542	ab179198 - H3 K4me3 + K9ac + S10p + K14ac + K18ac
K17	27.33198854	ab179199 - H3 K4me3 + T6p + K9ac + K14ac + K18ac
K18	21.03819583	ab179200 - H3 K4me2 + T6p + K9ac + K14ac + K18ac
K19	29.57341146	ab179201 - H3 K4bio + K9bio + K18bio
K20	20.18613958	ab179202 - H3 T6p + K9me3
K21	41.13002708	ab179203 - H3 T6p + R8me2a + K9me3
K22	20.29603229	ab179204 - H3 T6p + K9ac + K14ac + K18ac
K23	24.51197188	ab179205 - H3 R8me1
K24	20.67517188	ab179206 - H3 R8me1 + K9me1
K25	20.80808229	ab179207 - H3 R8me1 + K9me2
L1	21.63709375	ab179208 - H3 R8me1 + K9me3
L2	20.17994896	ab179209 - H3 R8me2s
L3	19.15287708	ab179210 - H3 R8me2s + K9me1
L4	20.30392708	ab179211 - H3 R8me2s + K9me2
L5	16.78351667	ab179212 - H3 R8me2s + K9me3
L6	18.30681667	ab179213 - H3 R8me2a
L7	19.45956771	ab179214 - H3 R8me2a + K9me1
L8	20.38757083	ab179215 - H3 R8me2a + K9me2
L9	22.04650938	ab179216 - H3 R8me2a + K9me3
L10	19.58141771	ab179217 - H3 R8cit
L11	33.6829375	ab179218 - H3 K9ac + K14ac
L12	35.57276771	ab179219 - H3 K9ac + K18ac
L13	18.50948542	ab179220 - H3 K9ac + K14ac + K18ac
L14	96.64608229	ab179221 - H3 K9ac + K14ac + K18ac + K23ac + K27ac
L15	17.95525104	ab179222 - H3 K9ac + S10p
L16	30.14671042	ab179223 - H3 K9cr
L17	21.97256667	ab179224 - H3 K9me2

L18	20.43595521	ab179225 - H3 K9me2 + S10p
L19	20.21680729	ab179226 - H3 K9me2 + K27me2
L20	22.35723542	ab179227 - H3 S10p + S28p
L21	22.58946667	ab179228 - H3 S10p K14ac
L22	22.48065208	ab179229 - H3 K14me1
L23	21.86745417	ab179230 - H3 K14me2
L24	23.80500313	ab179231 - H3 K14me3
L25	22.99654479	ab179232 - H3 K14ac + K18ac
M1	21.42672083	ab179233 - H3 R17cit
M2	21.8735625	ab179234 - H3 R17me1
M3	18.22696042	ab179235 - H3 R17me2s
M4	20.04176875	ab179236 - H3 R17me2a + K18ac
M5	19.85411146	ab179237 - H3 K18cr
M6	25.30824167	ab179238 - H3 K18me1
M7	20.65979375	ab179239 - H3 K18me2
M8	23.62560625	ab179240 - H3 K18me3 + K36me3
M9	23.9863125	ab179241 - H3 K23cr
M10	19.69450729	ab179242 - H3 K23me1
M11	22.42336979	ab179243 - H3 K23me2
M12	16.64469375	ab179244 - H3 K23me3
M13	20.90783854	ab179245 - H3 T22p + K23me3
M14	1654.146181	ab179246 - H3 K23pr
M15	27.91206771	ab179247 - H3 unmod(20-51)
M16	25.34131771	ab179248 - H3 R26me1
M17	25.66723229	ab179249 - H3 R26me2a
M18	20.97756563	ab179250 - H3 R26me2a + K27me1
M19	24.9118375	ab179251 - H3 R26me2a + K27me3
M20	24.89839167	ab179252 - H3 K27ac + S28p
M21	16.04445	ab179253 - H3 K27cr
M22	21.12555521	ab179254 - H3 K27me1
M23	20.09185729	ab179255 - H3 K27me2 + S28p
M24	16.29084167	ab179256 - H3 K27me3
M25	19.98970417	ab179257 - H3 K27me3 + S28p
N1	22.4793625	ab179258 - H3 P30hpro
N2	17.90549271	ab179259 - H3 P30(5,5me2)
N3	21.78569792	ab179260 - H3 P30
N4	22.96535	ab179261 - H3 P30mutA
N5	22.03450417	ab179262 - H3 T32GlcNAc
N6	25.25376458	ab179263 - H3 K36ac
N7	24.40239583	ab179264 - H3 P38hpro
N8	23.44426771	ab179265 - H3 P38(5,5me2)
N9	16.37562188	ab179266 - H3 P38
N10	21.64556875	ab179267 - H3 P38mutA
N11	20.59168021	ab179268 - H3 Y41p
N12	21.57175313	ab179269 - H3 unmod(50-80)
N13	22.47392083	ab179270 - H3 K56ac
N14	16.80858125	ab179271 - H3 K56cr
N15	19.75541979	ab179272 - H3 K56fo
N16	15.8959625	ab179273 - H3 K56me1
N17	25.90084375	ab179274 - H3 K56me2
N18	19.81780833	ab179275 - H3 K56me3
N19	23.61202083	ab179276 - H3 R63me1
N20	16.69798542	ab179277 - H3 R63me2s
N21	20.18052604	ab179278 - H3 R63me2a
N22	21.62939167	ab179279 - H3 K64ac
N23	20.10846042	ab179280 - H3 K64me2
N24	32.20384688	ab179281 - H3 K64me3
N25	26.63457396	ab179282 - H3 unmod(73-103)
O1	29.07227292	ab179283 - H3 K79ac
O2	23.35889688	ab179284 - H3 K79fo
O3	15.622275	ab179285 - H3 T80p
O4	27.2039625	ab179286 - H3 K79me3 + T80p
O5	19.69332604	ab179287 - H3 unmod(99-135)
O6	24.63340729	ab179288 - H3 T118O-ac
O7	20.48126354	ab179289 - H3 K122fo
O8	23.07819167	ab179290 - H3 K122me1
O9	22.15268958	ab179291 - H3 R128me1
O10	21.1300375	ab179292 - H3 R128me2s
O11	27.53774479	ab179293 - H3.3 unmod(10-40)
O12	20.79608021	ab179294 - H3.3 S28p
O13	22.75184479	ab179295 - H3.3 S28p + S31p
O14	24.7744875	ab179296 - H3.3 S31p
O15	33.76255833	ab179297 - H4 unmod(1-30)
O16	23.98446563	ab179298 - H4 S1p + R3me1

O17	22.08699167	ab179299 - H4 S1p + R3me2s
O18	25.54536979	ab179300 - H4 S1p + R3me2a
O19	31.66227917	ab179301 - H4 S1p + K5ac + K8ac + K12ac + K16ac
O20	23.80580833	ab179302 - H4 R3cit
O21	134.9210646	ab179303 - H4 R3me1 + K5ac + K8ac + K12ac + K16ac + K20ac
O22	25.60916354	ab179304 - H4 R3me2s
O23	120.4048438	ab179305 - H4 R3me2s + K5ac + K8ac + K12ac + K16ac + K20ac
O24	24.69128542	ab179306 - H4 R3me2a
O25	24.89998646	ab179307 - H4 R3me2a + K5ac
P1	84.82021979	ab179308 - H4 R3me2a + K5ac + K8ac + K12ac + K16ac + K20ac
P2	21.08565313	ab179309 - H4 K5ac
P3	27.53018021	ab179310 - H4 K5ac + K8ac
P4	25.4034	ab179311 - H4 K5ac + K12ac
P5	34.918925	ab179312 - H4 K5ac + K16ac
P6	25.22213646	ab179313 - H4 K5ac + K8ac + K12ac
P7	35.30065625	ab179314 - H4 K5ac + K8me1 + K12ac
P8	39.95635938	ab179315 - H4 K5ac + K8ac + K12ac + K16ac
P9	35.97284271	ab179316 - H4 K5ac + K8me1 + K12ac + K16ac
P10	54.63962292	ab179317 - H4 K5ac + K8ac + K12me1 + K16ac
P11	30.31893021	ab179318 - H4 K5me1
P12	26.162275	ab179319 - H4 K5me1 + K8ac + K12me1
P13	33.20836875	ab179320 - H4 K5me1 + K8me1 + K12me1
P14	26.37093229	ab179321 - H4 K5me1 + K8ac + K12ac + K16ac
P15	28.68815938	ab179322 - H4 K5me2
P16	23.21178542	ab179323 - H4 K5me3
P17	26.40106458	ab179324 - H4 K5cr
P18	19.47386042	ab179325 - H4 K8ac + K12ac
P19	24.52429792	ab179326 - H4 K8ac + K16ac
P20	25.6351125	ab179327 - H4 K8cr
P21	23.10808021	ab179328 - H4 K8me1
P22	23.6352125	ab179329 - H4 K12ac
P23	22.40815521	ab179330 - H4 K12ac + K16ac
P24	27.35755521	ab179331 - H4 K12ac + K16ac + K20me2
P25	22.82205313	ab179332 - H4 K12ac + K16ac + K20me3
Q1	34.27928125	ab179333 - H4 K5mutQ + K8mutQ + K12ac + K16mutQ + K20mutQ
Q2	21.78429896	ab179334 - H4 K12cr
Q3	24.02235729	ab179335 - H4 K12fo
Q4	20.49315208	ab179336 - H4 K12me1
Q5	21.79337083	ab179337 - H4 K12me2
Q6	22.04568021	ab179338 - H4 K12me3
Q7	22.69977813	ab179339 - H4 K16ac
Q8	42.255025	ab179340 - H4 K16ac + K20ac
Q9	23.60690313	ab179341 - H4 K16cr
Q10	21.9934875	ab179342 - H4 K16me1
Q11	23.60735313	ab179343 - H4 K16me3
Q12	33.92941042	ab179344 - H4 R17me1
Q13	27.48960521	ab179345 - H4 R17me2s
Q14	28.90644167	ab179346 - H4 R17me2a
Q15	26.45593229	ab179347 - H4 R19me1
Q16	22.35351875	ab179348 - H4 R19me2s
Q17	27.960125	ab179349 - H4 R19me2a
Q18	26.30765417	ab179350 - H4 K20ac
Q19	24.958975	ab179351 - H4 R23me1
Q20	20.01033958	ab179352 - H4 R23me2s
Q21	24.93805833	ab179353 - H4 R23me2a
Q22	25.26985729	ab179354 - H4 unmod(24-54)
Q23	21.98507813	ab179355 - H4 K31fo
Q24	22.57457083	ab179356 - H4 K31me1
Q25	19.45768021	ab179357 - H4 K31me2
R1	22.03079792	ab179358 - H4 K31me3
R2	32.68879063	ab179359 - H4 P32hyd
R3	21.76745208	ab179360 - H4 R35me1
R4	26.33315521	ab179361 - H4 R35me2s
R5	19.53829167	ab179362 - H4 R35me2a
R6	21.89105208	ab179363 - H4 S47O-ac
R7	19.61377083	ab179364 - H4 S47GlcNAc
R8	21.95824271	ab179365 - H4 S47p
R9	22.29063021	ab179366 - H4 unmod(45-75)
R10	21.74551458	ab179367 - H4 Y51hyd
R11	21.29415625	ab179368 - H4 R55me1
R12	24.68701458	ab179369 - H4 R55me2s
R13	20.54691458	ab179370 - H4 R55me2a
R14	25.22054583	ab179371 - H4 K59fo
R15	22.64440208	ab179372 - H4 K59me1

R16	20.83290104	ab179373 - H4 K59me2
R17	20.44726771	ab179374 - H4 R67me1
R18	22.88673958	ab179375 - H4 R67me2s
R19	24.09325104	ab179376 - H4 R67me2a
R20	22.76139375	ab179377 - H4 unmod(71-102)
R21	25.40272188	ab179378 - H4 K77me1
R22	19.62821458	ab179379 - H4 K77me2
R23	21.51851875	ab179380 - H4 K77me3
R24	24.92148646	ab179381 - H4 K79fo
R25	23.40234792	ab179382 - H4 K79me1
S1	20.1271625	ab179383 - H4 K79me3
S2	25.71199688	ab179384 - H4 T82O-ac
S3	24.41076458	ab179385 - H4 Y88hyd
S4	25.80185	ab179386 - H4 Y88nitro
S5	25.81788229	ab179387 - H4 K91fo
S6	20.18751042	ab179403 - H2A K5ac
S7	21.41411667	ab179404 - H2A K9me1
S8	21.39514479	ab18504 - H2B S32p
S9	23.97382813	ab19828 - ScH2A S129p
S10	47.795825	ab20631 - H3 R26cit
S11	22.57948021	ab20635 - H3 T3p + K4me3
S12	23.17534375	ab21043 - H3 K37me1
S13	21.88150625	ab21998 - H1.4 K25me2
S14	20.98265625	ab22079 - H4 K59me3
S15	19.93329375	ab22080 - H3 K122me2
S16	21.56389583	ab22082 - H3 K122me3
S17	26.08975625	ab22214 - H3 K37me2
S18	28.40521563	ab22398 - H2A R3me2a
S19	28.29642396	ab22399 - H2A R3me2s
S20	21.06679896	ab23539 - H3 T22p
S21	25.33674688	ab24003 - H3 K18ac
S22	16.64760625	ab24404 - H3 K27ac
S23	23.49524688	ab24417 - H3 K37me3
S24	19.49283438	ab24444 - H3 T11p
S25	63.11779271	ab24659 - H2A K127me3
T1	22.06217292	ab24660 - H2A K125me3
T2	21.03440521	ab26307 - H3 T45p
T3	20.11262396	ab26342 - H2B K5me2
T4	24.32491458	ab26739 - H4 T30p
T5	22.98786146	ab27896 - H4 T96p
T6	20.54361875	ab27898 - H4 T80p
T7	34.76565625	ab28756 - H2A R3me1
T8	24.16146458	ab32876 - H3 R2cit + R8cit + R17cit
T9	58.98641979	ab33047 - H2A K127me2
T10	17.68629271	ab33505 - H3 T118p
T11	20.615225	ab33594 - H3 R128me2a
T12	26.81399583	ab34466 - H3 K122ac
T13	20.05767188	ab36927 - H2A S122p
T14	23.9949375	ab4555 - H3 K79me1
T15	23.19839375	ab4556 - H3 K79me2
T16	23.69047083	ab4557 - H3 K79me3
T17	21.20232604	ab4560 - H4 K79me2
T18	22.73326458	ab46854 - H3 K18me3
T19	171.310326	ab48359 - H3 K23ac
T20	22.84677604	ab53244 - H3 K64me1
T21	40.89833333	ab54016 - H2A K9ac
T22	16.32476458	ab5499 - H3 S28p
T23	25.35831042	ab7768 - H3 K4me2
T24	25.71350104	ab89018 - H1.4 S26p
T25	22.64245	ab92374 - H3 K4me3 (HEX1)
U1	31.06208542	ab95444 - H4 K16me2