

Peptide array: Histone H3 (phospho S28) (ab10543)

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

Cell Ref	AUC	Peptide
A1	21.98780938	ab10112 - H4 K91ac
A2	20.01255	ab10121 - H2B K85ac
A3	19.44241875	ab10127 - H2A K95me2
A4	21.85720417	ab10138 - H1.4 T17p
A5	31.44377708	ab10139 - H1.4 T146p
A6	19.78381563	ab113997 - H2A.X Y142p
A7	23.40675521	ab11477 - H3 S10p
A8	13.90663229	ab12946 - H2B K5me3
A9	21.38384792	ab12949 - H2B K23me2
A10	19.16521458	ab12952 - H2A K127me1
A11	14.68102813	ab13211 - H2B K5me1
A12	20.71901667	ab1340 - H3 K4me1
A13	23.54715521	ab13570 - H2B K43me1
A14	22.78864792	ab13835 - HA
A15	23.10469896	ab13837 - Myc
A16	21.35563333	ab140408 - ScH2B T129p
A17	20.65007917	ab14103 - H3 T6p
A18	25.27024583	ab14724 - H4 S1p
A19	57.59126563	ab14799 - H3 T32p
A20	21.55713958	ab14943 - His
A21	20.31218229	ab14952 - H2A R77me2s
A22	26.83189479	ab14964 - H4 K20me2
A23	20.5706	ab14997 - ScH2A K21ac
A24	23.34772813	ab154782 - H3 R26me2s
A25	19.69578646	ab15591 - H3 K14ac
B1	19.13041667	ab15644 - H3 K9me3 + S10p
B2	17.27567396	ab15645 - H2A.X S139p
B3	23.58945417	ab15824 - H4 K8ac
B4	20.39628958	ab15829 - V5
B5	16.06282396	ab16635 - H3 K9ac
B6	32.78959792	ab166685 - H2B S112GlcNAc
B7	19.04268021	ab166688 - H2B S36GlcNAc
B8	20.85854896	ab16935 - H3 R17me2a
B9	26.03669063	ab17043 - H4 K20me1
B10	23.43306042	ab17162 - H3 P16hyd + P30hyd
B11	23.28116354	ab17567 - H4 K20me3
B12	13.93233854	ab17587 - H1.4 K25me3
B13	16.43823646	ab175905 - H2B K11ac
B14	21.55857292	ab176211 - H2B K23ac
B15	33.11118958	ab17632 - H4 K5bio
B16	21.02287604	ab1771 - H3 K9me1
B17	22.08357917	ab1773 - H3 K9me3
B18	22.71040521	ab17770 - H4 R3me1
B19	20.93570208	ab1781 - H3 K27me2
B20	21.25596146	ab1783 - H3 K36me1
B21	22.80595938	ab1784 - H3 K36me2
B22	22.15533542	ab1785 - H3 K36me3
B23	28.90120938	ab178667 - ScH2A unmod(1-30)
B24	27.84495521	ab178668 - ScH2A unmod(102-132)
B25	26.32457813	ab178670 - H2A.Z unmod(1-30)
C1	20.54827396	ab178833 - Flag
C2	16.96144063	ab178835 - H2B R99mutO
C3	24.91168333	ab178836 - H4 R23mutO
C4	24.16536458	ab178837 - H4 G48mutD
C5	19.6427875	ab178987 - H1.4 unmod(1-30)
C6	20.39600833	ab178988 - H1.4 S1p
C7	21.46936354	ab178989 - H1.4 T3p
C8	22.5387125	ab178990 - H1.4 S1p + T3p
C9	17.85279375	ab178991 - H1.4 K25me1
C10	136.6858177	ab178992 - H1.4 K25me3 + S26p
C11	20.28439896	ab178993 - H1.4 unmod(130-160)
C12	27.24031563	ab178994 - H2A unmod(1-30)
C13	46.60791354	ab178995 - H2A S1p
C14	19.59036146	ab178996 - H2A S1p + K5ac
C15	18.79283438	ab178997 - H2A S1p + R3me2a + K5ac
C16	28.89573958	ab178998 - H2A S1p + R3me2s + K5ac
C17	18.01609583	ab178999 - H2A S1p + R3cit + K5ac
C18	26.64298854	ab179000 - H2A S1ac + S1p + K5ac + K9ac + K13ac + K15ac
C19	17.55917604	ab179001 - H2A S1p + K5ac + K9ac + K13ac + K15ac
C20	46.31592708	ab179002 - H2A R3cit

C21	11.90274167	ab179003 - H2A R3cit + K5ac
C22	16.80156146	ab179004 - H2A R3me2a + K5ac
C23	15.56580729	ab179005 - H2A R3me2s + K5ac
C24	29.71299271	ab179006 - H2A S1ac + K5ac + K9ac + K13ac + K15ac
C25	17.60376042	ab179007 - H2A K5ac + K9ac + K13ac + K15ac
D1	19.66618229	ab179008 - H2A K9me2
D2	18.19966146	ab179009 - H2A K9me3
D3	21.72224479	ab179010 - H2A R11me1
D4	21.71719375	ab179011 - H2A R11me2s
D5	19.21795104	ab179012 - H2A R11me2a
D6	20.09107396	ab179013 - H2A K13ac
D7	22.34004063	ab179014 - H2A K15ac
D8	19.70565938	ab179015 - H2A R17me1
D9	15.55229896	ab179016 - H2A R17me2s
D10	19.10111979	ab179017 - H2A R17me2a
D11	25.34291875	ab179018 - H2A unmod(20-50)
D12	19.62394583	ab179019 - H2A P26hyd
D13	20.48640104	ab179020 - H2A R29me1
D14	18.80212083	ab179021 - H2A R29me2s
D15	19.47937604	ab179022 - H2A R29me2a
D16	18.52108125	ab179023 - H2A K36ac
D17	23.4290375	ab179024 - H2A K36cr
D18	18.01924167	ab179025 - H2A Y39hyd
D19	26.83538229	ab179026 - H2A R42me1
D20	16.50044792	ab179027 - H2A R42me2s
D21	25.03117292	ab179028 - H2A R42me2a
D22	15.19938958	ab179029 - H2A unmod(71-103)
D23	22.59550208	ab179030 - H2A R77me1
D24	14.16012396	ab179031 - H2A R77me2a
D25	16.31835833	ab179032 - H2A R88me1
E1	17.43486354	ab179033 - H2A R88me2s
E2	17.15962604	ab179034 - H2A R88me2a
E3	17.08995625	ab179035 - H2A K95me1
E4	17.21814271	ab179036 - H2A K95me3
E5	21.11937813	ab179037 - H2A unmod(92-123)
E6	18.59124688	ab179038 - H2A T101GlcNAc
E7	21.20810417	ab179039 - H2A Q104me1
E8	20.39521354	ab179040 - H2A unmod(112-129)
E9	23.38167396	ab179041 - H2A K118me1
E10	21.46692813	ab179042 - H2A K118me2
E11	20.45973854	ab179043 - H2A K118me3
E12	21.04032813	ab179044 - H2A K118fo
E13	21.42719583	ab179045 - H2A K118cr
E14	18.94646979	ab179046 - H2A K119cr
E15	17.77266979	ab179047 - H2A K118ubiq
E16	20.3063375	ab179048 - H2A K119ubiq
E17	20.83854688	ab179049 - H2A T120p
E18	20.10595938	ab179050 - H2A K125me1
E19	18.71946563	ab179051 - H2A K125me2
E20	23.99580313	ab179052 - H2A K125cr
E21	19.03534375	ab179053 - H2A.X unmod(1-30)
E22	19.91333333	ab179054 - H2A.X K5ac
E23	16.22561042	ab179055 - H2A.X unmod(114-142)
E24	17.38225417	ab179056 - H2A.X K118ubiq
E25	15.00753229	ab179057 - H2A.X K119ubiq
F1	19.77655104	ab179058 - H2A.X T120p
F2	17.04665104	ab179059 - H2A.X K127ubiq
F3	18.89699896	ab179060 - H2A.X S139p + Y142p
F4	17.16353229	ab179061 - H2A.Z K4ac
F5	16.99448021	ab179062 - H2A.Z K7ac
F6	17.21324271	ab179063 - H2A.Z K11ac
F7	16.06864479	ab179064 - H2A.Z K4ac + K7ac + K11ac
F8	18.63631458	ab179065 - H2B unmod(1-30)
F9	15.50232604	ab179066 - H2B K5ac
F10	20.45225208	ab179067 - H2B K5cr
F11	15.88319792	ab179068 - H2B K5fo
F12	14.26420833	ab179069 - H2B P10hyd
F13	20.65713854	ab179070 - H2B K11me1
F14	20.15583958	ab179071 - H2B K11me2
F15	20.84753542	ab179072 - H2B K11me3
F16	17.36672813	ab179073 - H2B K11cr
F17	21.12228438	ab179074 - H2B K12ac
F18	19.40085833	ab179075 - H2B K12cr
F19	21.42103438	ab179076 - H2B K12me1

F20	16.42360417	ab179077 - H2B K12me2
F21	22.29971458	ab179078 - H2B K12me3
F22	59.75503333	ab179079 - H2B K12ac + K15ac
F23	21.33565	ab179080 - H2B S14p
F24	26.24074479	ab179081 - H2B K15ac
F25	20.58101563	ab179082 - H2B K15cr
G1	21.68808958	ab179083 - H2B K15me1
G2	20.20229375	ab179084 - H2B K15me2
G3	20.64831354	ab179085 - H2B K15me3
G4	16.69070938	ab179086 - H2B K16ac
G5	20.6552	ab179087 - H2B K16cr
G6	19.67457604	ab179088 - H2B K20ac
G7	20.75551146	ab179089 - H2B K20cr
G8	20.98696979	ab179090 - H2B K20me1
G9	21.02703854	ab179091 - H2B K20me2
G10	18.70938438	ab179092 - H2B K20me3
G11	17.35450208	ab179093 - H2B K23cr
G12	16.215175	ab179094 - H2B K23me1
G13	16.70071146	ab179095 - H2B K23me3
G14	19.16948021	ab179096 - H2B K24ac
G15	26.40193229	ab179097 - H2B unmod(26-56)
G16	118.7128521	ab179098 - H2B K34cr
G17	22.31888958	ab179099 - H2B K34fo
G18	24.51557813	ab179100 - H2B Y37hyd
G19	19.55804896	ab179101 - H2B Y37p
G20	32.69878125	ab179102 - H2B K43me2
G21	25.02768542	ab179103 - H2B K43me3
G22	25.62477917	ab179104 - H2B K46fo
G23	19.88154063	ab179105 - H2B K46ac
G24	20.5539	ab179106 - H2B K46me1
G25	18.40313542	ab179107 - H2B K46me2
H1	21.43147188	ab179108 - H2B K46me3
H2	22.83569479	ab179109 - H2B unmod(51-81)
H3	21.77469271	ab179110 - H2B K57me1
H4	22.54729167	ab179111 - H2B K57me2
H5	22.80762813	ab179112 - H2B K57me3
H6	18.62011875	ab179113 - H2B unmod(73-103)
H7	20.81501979	ab179114 - H2B R79me1
H8	19.26805625	ab179115 - H2B R79me2s
H9	21.16374167	ab179116 - H2B R79me2a
H10	20.4756375	ab179117 - H2B Y83hyd
H11	16.68151354	ab179118 - H2B K85me1
H12	16.08315521	ab179119 - H2B K85me2
H13	17.70133438	ab179120 - H2B K85me3
H14	16.60320104	ab179121 - H2B unmod(94-125)
H15	17.31776563	ab179122 - H2B R99me1
H16	19.17468333	ab179123 - H2B R99me2s
H17	16.75488229	ab179124 - H2B R99me2a
H18	17.9721375	ab179125 - H2B K108fo
H19	21.77263229	ab179126 - H2B K116fo
H20	18.68911458	ab179127 - H2B K116me1
H21	23.41142396	ab179128 - H2B K116me2
H22	16.62689792	ab179129 - H2B K116me3
H23	25.2615375	ab179130 - H2B K116ubiq
H24	20.76674792	ab179131 - H2B K120ac
H25	23.00568438	ab179132 - H2B K120fo
I1	17.63230313	ab179133 - H2B K120ubiq
I2	19.82907813	ab179134 - H2B K125ubiq
I3	22.17688958	ab179135 - Sch2B unmod(102-132)
I4	21.26298646	ab179136 - H3 unmod(1-33)
I5	20.48433542	ab179137 - H3 R2cit
I6	20.18131771	ab179138 - H3 R2cit + K4me2
I7	20.09959792	ab179139 - H3 R2cit + K4me3
I8	20.42383021	ab179140 - H3 R2cit + K4me3 + K9ac + K14ac + K18ac
I9	18.87935938	ab179141 - H3 R2cit + K4ac + K9ac + K14ac + K18ac
I10	21.07813958	ab179142 - H3 R2me1
I11	22.82445208	ab179143 - H3 R2me1 + K4me2
I12	20.65202188	ab179144 - H3 R2me1 + K4me3
I13	22.48950208	ab179145 - H3 R2me1 + K4me3 + K9ac + K14ac + K18ac
I14	20.98175833	ab179146 - H3 R2me2s
I15	21.89212188	ab179147 - H3 R2me2s + K4me2
I16	21.40396979	ab179148 - H3 R2me2s + K4me3
I17	24.29892813	ab179149 - H3 R2me2s + K4me3 + K9ac + K14ac + K18ac
I18	22.40459896	ab179150 - H3 R2me2a

I19	20.87054792	ab179151 - H3 R2me2a + K4me2
I20	22.54120833	ab179152 - H3 R2me2a + K4me3
I21	28.35396563	ab179153 - H3 R2me2a + K4me3 + S10p
I22	19.29664271	ab179154 - H3 R2me2a + T3p + K4me3
I23	22.78461146	ab179155 - H3 R2me2a + K4ac + K9ac + K14ac + K18ac
I24	19.91985938	ab179156 - H3 R2me2a + K4me3 + K9ac + K14ac + K18ac
I25	20.21772813	ab179157 - H3 R2me2a + T3p + K4me3 + K9ac + K14ac + K18ac
J1	21.37465625	ab179158 - H3 R2me2a + K4me3 + K9ac + S10p + K14ac + K18ac
J2	20.75872813	ab179159 - H3 T3p
J3	19.69333542	ab179160 - H3 T3p + K4me1
J4	21.09851458	ab179161 - H3 T3p + K4me2
J5	20.35961563	ab179162 - H3 T3p + K4me3 + K9ac + K14ac + K18ac
J6	21.09598229	ab179163 - H3 K4ac
J7	21.61910938	ab179164 - H3 K4ac + K9ac
J8	20.35306667	ab179165 - H3 K4ac + K14ac
J9	23.57124688	ab179166 - H3 K4ac + K18ac
J10	20.70735521	ab179167 - H3 K4ac + K9me3
J11	20.86686146	ab179168 - H3 K4ac + K9ac + K14ac
J12	21.47965417	ab179169 - H3 K4ac + K9ac + K18ac
J13	22.00529479	ab179170 - H3 K4ac + K14ac + K18ac
J14	19.51055729	ab179171 - H3 K4ac + K9ac + K14ac + K18ac
J15	23.54644583	ab179172 - H3 K4ac + K9me1 + K14ac + K18ac
J16	19.72086771	ab179173 - H3 K4ac + K9me2 + K14ac + K18ac
J17	23.75660938	ab179174 - H3 K4ac + K9me3 + K14ac + K18ac
J18	19.89841667	ab179175 - H3 K4ac + K9ac + S10p + K14ac + K18ac
J19	23.53263229	ab179176 - H3 K4cr
J20	25.24621042	ab179177 - H3 K4me1 + K9me2
J21	23.87635313	ab179178 - H3 K4me1 + K9ac + K18ac
J22	19.37847083	ab179179 - H3 K4me1 + K9ac + K14ac + K18ac
J23	21.97269375	ab179180 - H3 K4me2 + T6p
J24	24.34725729	ab179181 - H3 K4me2 + K9me2
J25	20.63288125	ab179182 - H3 K4me2 + K9ac + K18ac
K1	20.42199688	ab179183 - H3 K4me2 + K9ac + K14ac + K18ac
K2	20.15293854	ab179184 - H3 K4me3 + T6p
K3	21.42285313	ab179185 - H3 K4me3 + K9ac
K4	21.1472375	ab179186 - H3 K4me3 + K9me2
K5	23.60541771	ab179187 - H3 K4me3 + K9me3
K6	20.24229063	ab179188 - H3 K4me3 + S10p
K7	19.0950625	ab179189 - H3 K4me3 + K14ac
K8	22.56629063	ab179190 - H3 K4me3 + K18ac
K9	18.61504583	ab179191 - H3 K4me3 + R8me2s + K9me3
K10	17.72603646	ab179192 - H3 K4me3 + R8me2a + K9me3
K11	25.40674271	ab179193 - H3 K4me3 + K9ac + S10p
K12	23.47001979	ab179194 - H3 K4me3 + K9ac + K14ac
K13	25.73706042	ab179195 - H3 K4me3 + K9ac + K18ac
K14	27.29207917	ab179196 - H3 K4me3 + K14ac + K18ac
K15	25.15418958	ab179197 - H3 K4me3 + K9ac + K14ac + K18ac
K16	27.29030833	ab179198 - H3 K4me3 + K9ac + S10p + K14ac + K18ac
K17	23.79482396	ab179199 - H3 K4me3 + T6p + K9ac + K14ac + K18ac
K18	17.78941042	ab179200 - H3 K4me2 + T6p + K9ac + K14ac + K18ac
K19	24.35269063	ab179201 - H3 K4bio + K9bio + K18bio
K20	20.56813646	ab179202 - H3 T6p + K9me3
K21	18.81782083	ab179203 - H3 T6p + R8me2a + K9me3
K22	20.48561563	ab179204 - H3 T6p + K9ac + K14ac + K18ac
K23	19.70101667	ab179205 - H3 R8me1
K24	19.97871667	ab179206 - H3 R8me1 + K9me1
K25	17.28358542	ab179207 - H3 R8me1 + K9me2
L1	32.70177917	ab179208 - H3 R8me1 + K9me3
L2	22.38414479	ab179209 - H3 R8me2s
L3	27.42098229	ab179210 - H3 R8me2s + K9me1
L4	23.82277917	ab179211 - H3 R8me2s + K9me2
L5	17.45237708	ab179212 - H3 R8me2s + K9me3
L6	18.94567813	ab179213 - H3 R8me2a
L7	19.53741146	ab179214 - H3 R8me2a + K9me1
L8	18.29654792	ab179215 - H3 R8me2a + K9me2
L9	15.99471771	ab179216 - H3 R8me2a + K9me3
L10	21.51312604	ab179217 - H3 R8cit
L11	24.07034063	ab179218 - H3 K9ac + K14ac
L12	24.23136458	ab179219 - H3 K9ac + K18ac
L13	17.94338646	ab179220 - H3 K9ac + K14ac + K18ac
L14	19.04408125	ab179221 - H3 K9ac + K14ac + K18ac + K23ac + K27ac
L15	16.82391771	ab179222 - H3 K9ac + S10p
L16	18.73394792	ab179223 - H3 K9cr
L17	19.66932604	ab179224 - H3 K9me2

L18	20.4618	ab179225 - H3 K9me2 + S10p
L19	22.61801042	ab179226 - H3 K9me2 + K27me2
L20	5533.445564	ab179227 - H3 S10p + S28p
L21	20.41649688	ab179228 - H3 S10p K14ac
L22	21.84723229	ab179229 - H3 K14me1
L23	23.92826042	ab179230 - H3 K14me2
L24	22.26088021	ab179231 - H3 K14me3
L25	21.02960417	ab179232 - H3 K14ac + K18ac
M1	19.35950208	ab179233 - H3 R17cit
M2	21.05340729	ab179234 - H3 R17me1
M3	22.32156667	ab179235 - H3 R17me2s
M4	22.35276146	ab179236 - H3 R17me2a + K18ac
M5	23.71666458	ab179237 - H3 K18cr
M6	22.57137292	ab179238 - H3 K18me1
M7	24.00679896	ab179239 - H3 K18me2
M8	19.03813333	ab179240 - H3 K18me3 + K36me3
M9	27.10304688	ab179241 - H3 K23cr
M10	18.44468542	ab179242 - H3 K23me1
M11	24.31117708	ab179243 - H3 K23me2
M12	15.8434125	ab179244 - H3 K23me3
M13	21.61825208	ab179245 - H3 T22p + K23me3
M14	21.72770313	ab179246 - H3 K23pr
M15	28.73642083	ab179247 - H3 unmod(20-51)
M16	23.80802188	ab179248 - H3 R26me1
M17	20.848675	ab179249 - H3 R26me2a
M18	19.93535625	ab179250 - H3 R26me2a + K27me1
M19	22.15265625	ab179251 - H3 R26me2a + K27me3
M20	3740.924774	ab179252 - H3 K27ac + S28p
M21	22.86713021	ab179253 - H3 K27cr
M22	34.14912813	ab179254 - H3 K27me1
M23	3177.601832	ab179255 - H3 K27me2 + S28p
M24	18.90113646	ab179256 - H3 K27me3
M25	3490.655333	ab179257 - H3 K27me3 + S28p
N1	22.67907292	ab179258 - H3 P30hpro
N2	209.7644313	ab179259 - H3 P30(5,5me2)
N3	22.96114896	ab179260 - H3 P30
N4	27.62887083	ab179261 - H3 P30mutA
N5	20.51	ab179262 - H3 T32GlcNAc
N6	21.85911146	ab179263 - H3 K36ac
N7	21.23764375	ab179264 - H3 P38hpro
N8	23.64393542	ab179265 - H3 P38(5,5me2)
N9	19.21901458	ab179266 - H3 P38
N10	29.77854583	ab179267 - H3 P38mutA
N11	21.82875	ab179268 - H3 Y41p
N12	24.37415521	ab179269 - H3 unmod(50-80)
N13	21.75895313	ab179270 - H3 K56ac
N14	20.50188646	ab179271 - H3 K56cr
N15	21.47476458	ab179272 - H3 K56fo
N16	12.76338542	ab179273 - H3 K56me1
N17	22.10427396	ab179274 - H3 K56me2
N18	17.80733646	ab179275 - H3 K56me3
N19	22.57413646	ab179276 - H3 R63me1
N20	19.23715313	ab179277 - H3 R63me2s
N21	21.61303125	ab179278 - H3 R63me2a
N22	17.39743125	ab179279 - H3 K64ac
N23	19.71629479	ab179280 - H3 K64me2
N24	29.25991458	ab179281 - H3 K64me3
N25	22.77368021	ab179282 - H3 unmod(73-103)
O1	23.92089792	ab179283 - H3 K79ac
O2	23.16490625	ab179284 - H3 K79fo
O3	17.48901979	ab179285 - H3 T80p
O4	21.24050313	ab179286 - H3 K79me3 + T80p
O5	23.54186979	ab179287 - H3 unmod(99-135)
O6	26.98817083	ab179288 - H3 T118O-ac
O7	18.51319479	ab179289 - H3 K122fo
O8	21.59796042	ab179290 - H3 K122me1
O9	25.58158438	ab179291 - H3 R128me1
O10	21.41017292	ab179292 - H3 R128me2s
O11	29.21474688	ab179293 - H3.3 unmod(10-40)
O12	2108.415742	ab179294 - H3.3 S28p
O13	2186.317942	ab179295 - H3.3 S28p + S31p
O14	25.73178958	ab179296 - H3.3 S31p
O15	34.09879583	ab179297 - H4 unmod(1-30)
O16	26.07387083	ab179298 - H4 S1p + R3me1

O17	24.10008333	ab179299 - H4 S1p + R3me2s
O18	24.05701667	ab179300 - H4 S1p + R3me2a
O19	26.13350417	ab179301 - H4 S1p + K5ac + K8ac + K12ac + K16ac
O20	26.55054375	ab179302 - H4 R3cit
O21	32.83345625	ab179303 - H4 R3me1 + K5ac + K8ac + K12ac + K16ac + K20ac
O22	22.83659063	ab179304 - H4 R3me2s
O23	27.79285833	ab179305 - H4 R3me2s + K5ac + K8ac + K12ac + K16ac + K20ac
O24	24.09163229	ab179306 - H4 R3me2a
O25	24.97465417	ab179307 - H4 R3me2a + K5ac
P1	30.3596875	ab179308 - H4 R3me2a + K5ac + K8ac + K12ac + K16ac + K20ac
P2	22.77504375	ab179309 - H4 K5ac
P3	25.173125	ab179310 - H4 K5ac + K8ac
P4	21.27208958	ab179311 - H4 K5ac + K12ac
P5	31.36092813	ab179312 - H4 K5ac + K16ac
P6	23.90205104	ab179313 - H4 K5ac + K8ac + K12ac
P7	25.56938438	ab179314 - H4 K5ac + K8me1 + K12ac
P8	30.29821146	ab179315 - H4 K5ac + K8ac + K12ac + K16ac
P9	30.4794125	ab179316 - H4 K5ac + K8me1 + K12ac + K16ac
P10	36.38909271	ab179317 - H4 K5ac + K8ac + K12me1 + K16ac
P11	26.02781042	ab179318 - H4 K5me1
P12	24.95992188	ab179319 - H4 K5me1 + K8ac + K12me1
P13	24.63386563	ab179320 - H4 K5me1 + K8me1 + K12me1
P14	30.21349792	ab179321 - H4 K5me1 + K8ac + K12ac + K16ac
P15	25.60160625	ab179322 - H4 K5me2
P16	23.53342188	ab179323 - H4 K5me3
P17	23.68054167	ab179324 - H4 K5cr
P18	21.97370833	ab179325 - H4 K8ac + K12ac
P19	25.35570938	ab179326 - H4 K8ac + K16ac
P20	24.28151354	ab179327 - H4 K8cr
P21	23.78346563	ab179328 - H4 K8me1
P22	23.52045417	ab179329 - H4 K12ac
P23	22.52142917	ab179330 - H4 K12ac + K16ac
P24	28.59162604	ab179331 - H4 K12ac + K16ac + K20me2
P25	24.52380208	ab179332 - H4 K12ac + K16ac + K20me3
Q1	31.72806146	ab179333 - H4 K5mutQ + K8mutQ + K12ac + K16mutQ + K20mutQ
Q2	24.24881667	ab179334 - H4 K12cr
Q3	21.76254271	ab179335 - H4 K12fo
Q4	22.01054688	ab179336 - H4 K12me1
Q5	24.41890938	ab179337 - H4 K12me2
Q6	20.68743229	ab179338 - H4 K12me3
Q7	22.492775	ab179339 - H4 K16ac
Q8	24.48857396	ab179340 - H4 K16ac + K20ac
Q9	21.91607396	ab179341 - H4 K16cr
Q10	22.82902813	ab179342 - H4 K16me1
Q11	24.71694792	ab179343 - H4 K16me3
Q12	27.27765729	ab179344 - H4 R17me1
Q13	29.20293021	ab179345 - H4 R17me2s
Q14	26.35425417	ab179346 - H4 R17me2a
Q15	23.42527083	ab179347 - H4 R19me1
Q16	26.75905833	ab179348 - H4 R19me2s
Q17	26.33328021	ab179349 - H4 R19me2a
Q18	25.74350313	ab179350 - H4 K20ac
Q19	39.69552813	ab179351 - H4 R23me1
Q20	25.38408021	ab179352 - H4 R23me2s
Q21	23.74155	ab179353 - H4 R23me2a
Q22	27.35943854	ab179354 - H4 unmod(24-54)
Q23	23.10905833	ab179355 - H4 K31fo
Q24	22.70741354	ab179356 - H4 K31me1
Q25	21.1846125	ab179357 - H4 K31me2
R1	23.23443021	ab179358 - H4 K31me3
R2	28.16639375	ab179359 - H4 P32hyd
R3	23.38179479	ab179360 - H4 R35me1
R4	24.10717292	ab179361 - H4 R35me2s
R5	24.50459896	ab179362 - H4 R35me2a
R6	20.63640417	ab179363 - H4 S47O-ac
R7	24.74194063	ab179364 - H4 S47GlcNAc
R8	20.90608125	ab179365 - H4 S47p
R9	22.20648542	ab179366 - H4 unmod(45-75)
R10	24.06177604	ab179367 - H4 Y51hyd
R11	27.04468958	ab179368 - H4 R55me1
R12	26.12463021	ab179369 - H4 R55me2s
R13	22.98996042	ab179370 - H4 R55me2a
R14	27.16923854	ab179371 - H4 K59fo
R15	21.60334792	ab179372 - H4 K59me1

R16	18.73437708	ab179373 - H4 K59me2
R17	19.70447604	ab179374 - H4 R67me1
R18	19.79838542	ab179375 - H4 R67me2s
R19	17.81292292	ab179376 - H4 R67me2a
R20	20.13561042	ab179377 - H4 unmod(71-102)
R21	20.56829375	ab179378 - H4 K77me1
R22	20.62437604	ab179379 - H4 K77me2
R23	22.48018646	ab179380 - H4 K77me3
R24	23.56613229	ab179381 - H4 K79fo
R25	21.47417188	ab179382 - H4 K79me1
S1	23.19	ab179383 - H4 K79me3
S2	27.3487	ab179384 - H4 T82O-ac
S3	26.6946125	ab179385 - H4 Y88hyd
S4	22.42874583	ab179386 - H4 Y88nitro
S5	22.68819167	ab179387 - H4 K91fo
S6	17.48879479	ab179403 - H2A K5ac
S7	14.50287604	ab179404 - H2A K9me1
S8	18.33011771	ab18504 - H2B S32p
S9	17.52485833	ab19828 - ScH2A S129p
S10	22.67155313	ab20631 - H3 R26cit
S11	18.54580833	ab20635 - H3 T3p + K4me3
S12	22.59549167	ab21043 - H3 K37me1
S13	20.00598333	ab21998 - H1.4 K25me2
S14	18.38349479	ab22079 - H4 K59me3
S15	22.33706563	ab22080 - H3 K122me2
S16	20.96532188	ab22082 - H3 K122me3
S17	27.05999583	ab22214 - H3 K37me2
S18	27.69225417	ab22398 - H2A R3me2a
S19	16.636075	ab22399 - H2A R3me2s
S20	39.11041146	ab23539 - H3 T22p
S21	20.18530625	ab24003 - H3 K18ac
S22	18.19358542	ab24404 - H3 K27ac
S23	25.29616667	ab24417 - H3 K37me3
S24	706.6970302	ab24444 - H3 T11p
S25	17.30634583	ab24659 - H2A K127me3
T1	17.30283542	ab24660 - H2A K125me3
T2	20.73335417	ab26307 - H3 T45p
T3	18.57428958	ab26342 - H2B K5me2
T4	25.35954688	ab26739 - H4 T30p
T5	24.81301667	ab27896 - H4 T96p
T6	22.34569792	ab27898 - H4 T80p
T7	27.52609479	ab28756 - H2A R3me1
T8	22.80522188	ab32876 - H3 R2cit + R8cit + R17cit
T9	29.15060313	ab33047 - H2A K127me2
T10	23.79279271	ab33505 - H3 T118p
T11	25.12214792	ab33594 - H3 R128me2a
T12	24.05316458	ab34466 - H3 K122ac
T13	20.50476042	ab36927 - H2A S122p
T14	23.84060625	ab4555 - H3 K79me1
T15	23.78681146	ab4556 - H3 K79me2
T16	24.27518333	ab4557 - H3 K79me3
T17	22.96826146	ab4560 - H4 K79me2
T18	21.19427604	ab46854 - H3 K18me3
T19	18.85002813	ab48359 - H3 K23ac
T20	24.60954583	ab53244 - H3 K64me1
T21	39.95238854	ab54016 - H2A K9ac
T22	4786.227945	ab5499 - H3 S28p
T23	24.44459167	ab7768 - H3 K4me2
T24	1067.312396	ab89018 - H1.4 S26p
T25	21.40108854	ab92374 - H3 K4me3 (HEX1)
U1	27.09324792	ab95444 - H4 K16me2