

Peptide array: Histone H2B (acetyl K15) (ab62335)

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

Cell Ref	AUC	Peptide
ab10112	25.38115521	ab10112 - H4 K91ac
ab10121	25.88488021	ab10121 - H2B K85ac
ab10127	20.68347292	ab10127 - H2A K95me2
ab10138	15.78201354	ab10138 - H1.4 T17p
ab10139	27.16058438	ab10139 - H1.4 T146p
ab113997	21.55846458	ab113997 - H2A.X Y142p
ab11477	22.79747604	ab11477 - H3 S10p
ab12946	19.54957396	ab12946 - H2B K5me3
ab12949	21.29349167	ab12949 - H2B K23me2
ab12952	22.93077708	ab12952 - H2A K127me1
ab13211	23.32628646	ab13211 - H2B K5me1
ab1340	20.38025313	ab1340 - H3 K4me1
ab13570	22.33392604	ab13570 - H2B K43me1
ab13835	22.94088021	ab13835 - HA
ab13837	24.15411042	ab13837 - Myc
ab140408	23.44878125	ab140408 - ScH2B T129p
ab14103	22.54877396	ab14103 - H3 T6p
ab14724	21.19238854	ab14724 - H4 S1p
ab14799	22.67915521	ab14799 - H3 T32p
ab14943	29.99963542	ab14943 - His
ab14952	22.09029167	ab14952 - H2A R77me2s
ab14964	22.62451042	ab14964 - H4 K20me2
ab14997	22.9787875	ab14997 - ScH2A K21ac
ab154782	20.51108021	ab154782 - H3 R26me2s
ab15591	20.28778229	ab15591 - H3 K14ac
ab15644	21.62947813	ab15644 - H3 K9me3 + S10p
ab15645	20.25050833	ab15645 - H2A.X S139p
ab15824	20.87095417	ab15824 - H4 K8ac
ab15829	19.93479688	ab15829 - V5
ab16635	21.89145313	ab16635 - H3 K9ac
ab166685	19.7227125	ab166685 - H2B S112GlcNAc
ab166688	20.31616667	ab166688 - H2B S36GlcNAc
ab16935	20.63849271	ab16935 - H3 R17me2a
ab17043	24.74653333	ab17043 - H4 K20me1
ab17162	20.67473438	ab17162 - H3 P16hyd + P30hyd
ab17567	24.03007396	ab17567 - H4 K20me3
ab17587	24.29042083	ab17587 - H1.4 K25me3
ab175905	17.41700313	ab175905 - H2B K11ac
ab176211	23.42174688	ab176211 - H2B K23ac
ab17632	23.61919479	ab17632 - H4 K5bio
ab1771	20.38885521	ab1771 - H3 K9me1
ab1773	22.92644375	ab1773 - H3 K9me3
ab17770	24.28246042	ab17770 - H4 R3me1
ab1781	20.83988125	ab1781 - H3 K27me2
ab1783	20.89380833	ab1783 - H3 K36me1
ab1784	19.93196771	ab1784 - H3 K36me2
ab1785	20.39179896	ab1785 - H3 K36me3
ab178667	27.58523021	ab178667 - ScH2A unmod(1-30)
ab178668	25.89416875	ab178668 - ScH2A unmod(102-132)
ab178670	24.03723542	ab178670 - H2A.Z unmod(1-30)
ab178833	24.79284583	ab178833 - Flag
ab178835	21.26838229	ab178835 - H2B R99mutO
ab178836	23.69834583	ab178836 - H4 R23mutO
ab178837	21.42093229	ab178837 - H4 G48mutD
ab178987	21.08764792	ab178987 - H1.4 unmod(1-30)
ab178988	26.54435729	ab178988 - H1.4 S1p
ab178989	17.18075729	ab178989 - H1.4 T3p
ab178990	26.47654479	ab178990 - H1.4 S1p + T3p
ab178991	24.31183125	ab178991 - H1.4 K25me1
ab178992	19.64623542	ab178992 - H1.4 K25me3 + S26p
ab178993	17.42369167	ab178993 - H1.4 unmod(130-160)
ab178994	20.37529375	ab178994 - H2A unmod(1-30)
ab178995	23.06095521	ab178995 - H2A S1p
ab178996	19.74250729	ab178996 - H2A S1p + K5ac
ab178997	23.48600417	ab178997 - H2A S1p + R3me2a + K5ac
ab178998	25.64394167	ab178998 - H2A S1p + R3me2s + K5ac
ab178999	21.25930104	ab178999 - H2A S1p + R3cit + K5ac
ab179000	25.50266042	ab179000 - H2A S1ac + S1p + K5ac + K9ac + K13ac + K15ac
ab179001	20.77143333	ab179001 - H2A S1p + K5ac + K9ac + K13ac + K15ac
ab179002	24.18806875	ab179002 - H2A R3cit

ab179003	18.17176875	ab179003 - H2A R3cit + K5ac
ab179004	19.75553438	ab179004 - H2A R3me2a + K5ac
ab179005	19.08486563	ab179005 - H2A R3me2s + K5ac
ab179006	31.05120417	ab179006 - H2A S1ac + K5ac + K9ac + K13ac + K15ac
ab179007	17.76020729	ab179007 - H2A K5ac + K9ac + K13ac + K15ac
ab179008	17.09665625	ab179008 - H2A K9me2
ab179009	24.37989375	ab179009 - H2A K9me3
ab179010	19.22910104	ab179010 - H2A R11me1
ab179011	23.91562604	ab179011 - H2A R11me2s
ab179012	22.96279688	ab179012 - H2A R11me2a
ab179013	24.3584375	ab179013 - H2A K13ac
ab179014	17.76466042	ab179014 - H2A K15ac
ab179015	23.72459167	ab179015 - H2A R17me1
ab179016	21.46197917	ab179016 - H2A R17me2s
ab179017	22.44044479	ab179017 - H2A R17me2a
ab179018	22.4831125	ab179018 - H2A unmod(20-50)
ab179019	22.76567813	ab179019 - H2A P26hyd
ab179020	20.66601146	ab179020 - H2A R29me1
ab179021	23.71853542	ab179021 - H2A R29me2s
ab179022	21.97010417	ab179022 - H2A R29me2a
ab179023	22.6820875	ab179023 - H2A K36ac
ab179024	22.926325	ab179024 - H2A K36cr
ab179025	20.75396875	ab179025 - H2A Y39hyd
ab179026	23.86817292	ab179026 - H2A R42me1
ab179027	20.69840729	ab179027 - H2A R42me2s
ab179028	24.64973229	ab179028 - H2A R42me2a
ab179029	20.73168542	ab179029 - H2A unmod(71-103)
ab179030	27.18454375	ab179030 - H2A R77me1
ab179031	21.35217604	ab179031 - H2A R77me2a
ab179032	21.74597917	ab179032 - H2A R88me1
ab179033	23.12364896	ab179033 - H2A R88me2s
ab179034	20.45784063	ab179034 - H2A R88me2a
ab179035	24.80826042	ab179035 - H2A K95me1
ab179036	25.8697125	ab179036 - H2A K95me3
ab179037	17.51808958	ab179037 - H2A unmod(92-123)
ab179038	22.06538958	ab179038 - H2A T101GlcNAc
ab179039	26.56115313	ab179039 - H2A Q104me1
ab179040	16.11807396	ab179040 - H2A unmod(112-129)
ab179041	24.23553854	ab179041 - H2A K118me1
ab179042	18.69639167	ab179042 - H2A K118me2
ab179043	23.23518646	ab179043 - H2A K118me3
ab179044	21.361425	ab179044 - H2A K118fo
ab179045	23.64221875	ab179045 - H2A K118cr
ab179046	21.30720729	ab179046 - H2A K119cr
ab179047	25.8783	ab179047 - H2A K118ubiq
ab179048	21.78715625	ab179048 - H2A K119ubiq
ab179049	26.10269688	ab179049 - H2A T120p
ab179050	22.06211563	ab179050 - H2A K125me1
ab179051	23.74306979	ab179051 - H2A K125me2
ab179052	23.6430875	ab179052 - H2A K125cr
ab179053	20.93880417	ab179053 - H2A.X unmod(1-30)
ab179054	24.64032083	ab179054 - H2A.X K5ac
ab179055	20.86814688	ab179055 - H2A.X unmod(114-142)
ab179056	26.03289375	ab179056 - H2A.X K118ubiq
ab179057	23.21925938	ab179057 - H2A.X K119ubiq
ab179058	20.34767813	ab179058 - H2A.X T120p
ab179059	22.29744167	ab179059 - H2A.X K127ubiq
ab179060	16.47558438	ab179060 - H2A.X S139p + Y142p
ab179061	15.23525729	ab179061 - H2A.Z K4ac
ab179062	18.13445313	ab179062 - H2A.Z K7ac
ab179063	19.37001875	ab179063 - H2A.Z K11ac
ab179064	19.54726458	ab179064 - H2A.Z K4ac + K7ac + K11ac
ab179065	26.93768542	ab179065 - H2B unmod(1-30)
ab179066	21.22598333	ab179066 - H2B K5ac
ab179067	19.86391875	ab179067 - H2B K5cr
ab179068	17.99537083	ab179068 - H2B K5fo
ab179069	21.92748542	ab179069 - H2B P10hyd
ab179070	20.50212396	ab179070 - H2B K11me1
ab179071	17.93124792	ab179071 - H2B K11me2
ab179072	20.43878125	ab179072 - H2B K11me3
ab179073	18.66377708	ab179073 - H2B K11cr
ab179074	21.75524792	ab179074 - H2B K12ac
ab179075	18.70895625	ab179075 - H2B K12cr
ab179076	18.56280938	ab179076 - H2B K12me1

ab179077	21.33092083	ab179077 - H2B K12me2
ab179078	21.34241875	ab179078 - H2B K12me3
ab179079	689.9474469	ab179079 - H2B K12ac + K15ac
ab179080	20.84193229	ab179080 - H2B S14p
ab179081	1875.51694	ab179081 - H2B K15ac
ab179082	28.91256979	ab179082 - H2B K15cr
ab179083	169.5628708	ab179083 - H2B K15me1
ab179084	20.34360521	ab179084 - H2B K15me2
ab179085	20.14860833	ab179085 - H2B K15me3
ab179086	19.3069625	ab179086 - H2B K16ac
ab179087	20.13345625	ab179087 - H2B K16cr
ab179088	20.92595	ab179088 - H2B K20ac
ab179089	16.65567708	ab179089 - H2B K20cr
ab179090	19.21057813	ab179090 - H2B K20me1
ab179091	21.66506771	ab179091 - H2B K20me2
ab179092	20.38216458	ab179092 - H2B K20me3
ab179093	21.99809479	ab179093 - H2B K23cr
ab179094	20.43463438	ab179094 - H2B K23me1
ab179095	18.58756563	ab179095 - H2B K23me3
ab179096	24.29121458	ab179096 - H2B K24ac
ab179097	25.38347396	ab179097 - H2B unmod(26-56)
ab179098	22.80439271	ab179098 - H2B K34cr
ab179099	22.82015208	ab179099 - H2B K34fo
ab179100	26.76904583	ab179100 - H2B Y37hyd
ab179101	21.68314583	ab179101 - H2B Y37p
ab179102	22.52947083	ab179102 - H2B K43me2
ab179103	19.57113958	ab179103 - H2B K43me3
ab179104	17.931425	ab179104 - H2B K46fo
ab179105	21.98791667	ab179105 - H2B K46ac
ab179106	19.90949896	ab179106 - H2B K46me1
ab179107	17.31759792	ab179107 - H2B K46me2
ab179108	19.19006563	ab179108 - H2B K46me3
ab179109	20.546325	ab179109 - H2B unmod(51-81)
ab179110	17.58089063	ab179110 - H2B K57me1
ab179111	20.86551875	ab179111 - H2B K57me2
ab179112	21.06357813	ab179112 - H2B K57me3
ab179113	19.16215833	ab179113 - H2B unmod(73-103)
ab179114	19.67586146	ab179114 - H2B R79me1
ab179115	20.54181563	ab179115 - H2B R79me2s
ab179116	21.95095313	ab179116 - H2B R79me2a
ab179117	20.94073333	ab179117 - H2B Y83hyd
ab179118	18.05501979	ab179118 - H2B K85me1
ab179119	21.22078646	ab179119 - H2B K85me2
ab179120	21.47284583	ab179120 - H2B K85me3
ab179121	20.13955208	ab179121 - H2B unmod(94-125)
ab179122	23.27971979	ab179122 - H2B R99me1
ab179123	23.54648854	ab179123 - H2B R99me2s
ab179124	24.20983333	ab179124 - H2B R99me2a
ab179125	19.68225729	ab179125 - H2B K108fo
ab179126	21.55309167	ab179126 - H2B K116fo
ab179127	20.74578333	ab179127 - H2B K116me1
ab179128	25.48992292	ab179128 - H2B K116me2
ab179129	17.05276771	ab179129 - H2B K116me3
ab179130	19.91503646	ab179130 - H2B K116ubiq
ab179131	19.73006563	ab179131 - H2B K120ac
ab179132	23.00532604	ab179132 - H2B K120fo
ab179133	17.64978958	ab179133 - H2B K120ubiq
ab179134	21.03758021	ab179134 - H2B K125ubiq
ab179135	22.1539625	ab179135 - Sch2B unmod(102-132)
ab179136	20.21637396	ab179136 - H3 unmod(1-33)
ab179137	21.52805938	ab179137 - H3 R2cit
ab179138	20.13128125	ab179138 - H3 R2cit + K4me2
ab179139	20.51014167	ab179139 - H3 R2cit + K4me3
ab179140	20.91986563	ab179140 - H3 R2cit + K4me3 + K9ac + K14ac + K18ac
ab179141	20.37450313	ab179141 - H3 R2cit + K4ac + K9ac + K14ac + K18ac
ab179142	20.58801354	ab179142 - H3 R2me1
ab179143	22.87705417	ab179143 - H3 R2me1 + K4me2
ab179144	20.41178438	ab179144 - H3 R2me1 + K4me3
ab179145	21.26610729	ab179145 - H3 R2me1 + K4me3 + K9ac + K14ac + K18ac
ab179146	23.06512604	ab179146 - H3 R2me2s
ab179147	24.02591458	ab179147 - H3 R2me2s + K4me2
ab179148	18.31966458	ab179148 - H3 R2me2s + K4me3
ab179149	23.081725	ab179149 - H3 R2me2s + K4me3 + K9ac + K14ac + K18ac
ab179150	24.61460833	ab179150 - H3 R2me2a

ab179151	22.15307813	ab179151 - H3 R2me2a + K4me2
ab179152	22.76467917	ab179152 - H3 R2me2a + K4me3
ab179153	20.65927292	ab179153 - H3 R2me2a + K4me3 + S10p
ab179154	23.11282708	ab179154 - H3 R2me2a + T3p + K4me3
ab179155	21.61299896	ab179155 - H3 R2me2a + K4ac + K9ac + K14ac + K18ac
ab179156	19.31457917	ab179156 - H3 R2me2a + K4me3 + K9ac + K14ac + K18ac
ab179157	21.76105625	ab179157 - H3 R2me2a + T3p + K4me3 + K9ac + K14ac + K18ac
ab179158	23.32466563	ab179158 - H3 R2me2a + K4me3 + K9ac + S10p + K14ac + K18ac
ab179159	19.945425	ab179159 - H3 T3p
ab179160	18.62500833	ab179160 - H3 T3p + K4me1
ab179161	21.17671875	ab179161 - H3 T3p + K4me2
ab179162	22.26504063	ab179162 - H3 T3p + K4me3 + K9ac + K14ac + K18ac
ab179163	20.70747604	ab179163 - H3 K4ac
ab179164	24.96387813	ab179164 - H3 K4ac + K9ac
ab179165	25.17752396	ab179165 - H3 K4ac + K14ac
ab179166	25.54426771	ab179166 - H3 K4ac + K18ac
ab179167	21.25102708	ab179167 - H3 K4ac + K9me3
ab179168	21.37613438	ab179168 - H3 K4ac + K9ac + K14ac
ab179169	17.99125521	ab179169 - H3 K4ac + K9ac + K18ac
ab179170	23.54396563	ab179170 - H3 K4ac + K14ac + K18ac
ab179171	20.46065	ab179171 - H3 K4ac + K9ac + K14ac + K18ac
ab179172	21.94223333	ab179172 - H3 K4ac + K9me1 + K14ac + K18ac
ab179173	23.75580938	ab179173 - H3 K4ac + K9me2 + K14ac + K18ac
ab179174	24.93668542	ab179174 - H3 K4ac + K9me3 + K14ac + K18ac
ab179175	23.46909063	ab179175 - H3 K4ac + K9ac + S10p + K14ac + K18ac
ab179176	24.81497083	ab179176 - H3 K4cr
ab179177	26.12139792	ab179177 - H3 K4me1 + K9me2
ab179178	20.72552708	ab179178 - H3 K4me1 + K9ac + K18ac
ab179179	22.90560625	ab179179 - H3 K4me1 + K9ac + K14ac + K18ac
ab179180	24.02096042	ab179180 - H3 K4me2 + T6p
ab179181	23.94564688	ab179181 - H3 K4me2 + K9me2
ab179182	22.96647292	ab179182 - H3 K4me2 + K9ac + K18ac
ab179183	22.27978333	ab179183 - H3 K4me2 + K9ac + K14ac + K18ac
ab179184	22.57272083	ab179184 - H3 K4me3 + T6p
ab179185	23.52481875	ab179185 - H3 K4me3 + K9ac
ab179186	22.09434792	ab179186 - H3 K4me3 + K9me2
ab179187	24.33600313	ab179187 - H3 K4me3 + K9me3
ab179188	20.97344479	ab179188 - H3 K4me3 + S10p
ab179189	22.35588438	ab179189 - H3 K4me3 + K14ac
ab179190	22.60758333	ab179190 - H3 K4me3 + K18ac
ab179191	23.45399479	ab179191 - H3 K4me3 + R8me2s + K9me3
ab179192	24.03427396	ab179192 - H3 K4me3 + R8me2a + K9me3
ab179193	24.05155833	ab179193 - H3 K4me3 + K9ac + S10p
ab179194	23.38698854	ab179194 - H3 K4me3 + K9ac + K14ac
ab179195	24.7470375	ab179195 - H3 K4me3 + K9ac + K18ac
ab179196	22.92100938	ab179196 - H3 K4me3 + K14ac + K18ac
ab179197	23.49920938	ab179197 - H3 K4me3 + K9ac + K14ac + K18ac
ab179198	24.68136979	ab179198 - H3 K4me3 + K9ac + S10p + K14ac + K18ac
ab179199	22.92321771	ab179199 - H3 K4me3 + T6p + K9ac + K14ac + K18ac
ab179200	19.89497813	ab179200 - H3 K4me2 + T6p + K9ac + K14ac + K18ac
ab179201	23.89976979	ab179201 - H3 K4bio + K9bio + K18bio
ab179202	21.144125	ab179202 - H3 T6p + K9me3
ab179203	20.58652083	ab179203 - H3 T6p + R8me2a + K9me3
ab179204	22.54463125	ab179204 - H3 T6p + K9ac + K14ac + K18ac
ab179205	23.82314792	ab179205 - H3 R8me1
ab179206	22.38223021	ab179206 - H3 R8me1 + K9me1
ab179207	21.64713125	ab179207 - H3 R8me1 + K9me2
ab179208	19.60968958	ab179208 - H3 R8me1 + K9me3
ab179209	19.09376771	ab179209 - H3 R8me2s
ab179210	23.29035313	ab179210 - H3 R8me2s + K9me1
ab179211	19.67713542	ab179211 - H3 R8me2s + K9me2
ab179212	20.73960208	ab179212 - H3 R8me2s + K9me3
ab179213	18.95121563	ab179213 - H3 R8me2a
ab179214	23.26210521	ab179214 - H3 R8me2a + K9me1
ab179215	21.88138958	ab179215 - H3 R8me2a + K9me2
ab179216	20.17116875	ab179216 - H3 R8me2a + K9me3
ab179217	22.28674167	ab179217 - H3 R8cit
ab179218	27.32319479	ab179218 - H3 K9ac + K14ac
ab179219	30.19305729	ab179219 - H3 K9ac + K18ac
ab179220	19.89207813	ab179220 - H3 K9ac + K14ac + K18ac
ab179221	22.41160938	ab179221 - H3 K9ac + K14ac + K18ac + K23ac + K27ac
ab179222	22.04364271	ab179222 - H3 K9ac + S10p
ab179223	18.13553125	ab179223 - H3 K9cr
ab179224	18.64339688	ab179224 - H3 K9me2

ab179225	21.97888125	ab179225 - H3 K9me2 + S10p
ab179226	21.77197292	ab179226 - H3 K9me2 + K27me2
ab179227	27.71717708	ab179227 - H3 S10p + S28p
ab179228	24.83602083	ab179228 - H3 S10p K14ac
ab179229	19.319925	ab179229 - H3 K14me1
ab179230	19.61490833	ab179230 - H3 K14me2
ab179231	22.00561458	ab179231 - H3 K14me3
ab179232	19.95371354	ab179232 - H3 K14ac + K18ac
ab179233	20.70866563	ab179233 - H3 R17cit
ab179234	18.66418958	ab179234 - H3 R17me1
ab179235	18.64188646	ab179235 - H3 R17me2s
ab179236	21.20490521	ab179236 - H3 R17me2a + K18ac
ab179237	19.99206458	ab179237 - H3 K18cr
ab179238	19.33657917	ab179238 - H3 K18me1
ab179239	23.33970938	ab179239 - H3 K18me2
ab179240	19.819525	ab179240 - H3 K18me3 + K36me3
ab179241	21.55663646	ab179241 - H3 K23cr
ab179242	21.88048229	ab179242 - H3 K23me1
ab179243	20.32328438	ab179243 - H3 K23me2
ab179244	22.71705521	ab179244 - H3 K23me3
ab179245	21.00739375	ab179245 - H3 T22p + K23me3
ab179246	26.14427188	ab179246 - H3 K23pr
ab179247	25.63074375	ab179247 - H3 unmod(20-51)
ab179248	18.31414688	ab179248 - H3 R26me1
ab179249	18.06729167	ab179249 - H3 R26me2a
ab179250	17.69120208	ab179250 - H3 R26me2a + K27me1
ab179251	20.34058646	ab179251 - H3 R26me2a + K27me3
ab179252	21.50250208	ab179252 - H3 K27ac + S28p
ab179253	21.11369583	ab179253 - H3 K27cr
ab179254	20.17457708	ab179254 - H3 K27me1
ab179255	18.75397708	ab179255 - H3 K27me2 + S28p
ab179256	18.92077708	ab179256 - H3 K27me3
ab179257	29.26699688	ab179257 - H3 K27me3 + S28p
ab179258	21.83779479	ab179258 - H3 P30hpro
ab179259	19.06377813	ab179259 - H3 P30(5,5me2)
ab179260	21.08699271	ab179260 - H3 P30
ab179261	22.32933542	ab179261 - H3 P30mutA
ab179262	20.94788958	ab179262 - H3 T32GlcNAc
ab179263	46.56200938	ab179263 - H3 K36ac
ab179264	21.06884271	ab179264 - H3 P38hpro
ab179265	22.37496875	ab179265 - H3 P38(5,5me2)
ab179266	16.71303229	ab179266 - H3 P38
ab179267	20.359	ab179267 - H3 P38mutA
ab179268	21.093975	ab179268 - H3 Y41p
ab179269	21.65731042	ab179269 - H3 unmod(50-80)
ab179270	21.34162917	ab179270 - H3 K56ac
ab179271	21.11562917	ab179271 - H3 K56cr
ab179272	20.34459792	ab179272 - H3 K56fo
ab179273	18.48186042	ab179273 - H3 K56me1
ab179274	24.49968125	ab179274 - H3 K56me2
ab179275	19.52060208	ab179275 - H3 K56me3
ab179276	23.75555417	ab179276 - H3 R63me1
ab179277	20.41583646	ab179277 - H3 R63me2s
ab179278	23.1545	ab179278 - H3 R63me2a
ab179279	20.30383438	ab179279 - H3 K64ac
ab179280	20.56258021	ab179280 - H3 K64me2
ab179281	26.17547188	ab179281 - H3 K64me3
ab179282	23.12322292	ab179282 - H3 unmod(73-103)
ab179283	21.73566771	ab179283 - H3 K79ac
ab179284	18.15834792	ab179284 - H3 K79fo
ab179285	20.77502708	ab179285 - H3 T80p
ab179286	24.44528542	ab179286 - H3 K79me3 + T80p
ab179287	22.18361979	ab179287 - H3 unmod(99-135)
ab179288	23.80768854	ab179288 - H3 T118O-ac
ab179289	22.01921146	ab179289 - H3 K122fo
ab179290	18.98959063	ab179290 - H3 K122me1
ab179291	22.65434688	ab179291 - H3 R128me1
ab179292	23.95919271	ab179292 - H3 R128me2s
ab179293	24.41060208	ab179293 - H3.3 unmod(10-40)
ab179294	21.18154688	ab179294 - H3.3 S28p
ab179295	23.77723646	ab179295 - H3.3 S28p + S31p
ab179296	24.20695313	ab179296 - H3.3 S31p
ab179297	27.82645938	ab179297 - H4 unmod(1-30)
ab179298	23.84069063	ab179298 - H4 S1p + R3me1

ab179299	22.51855	ab179299 - H4 S1p + R3me2s
ab179300	25.55649688	ab179300 - H4 S1p + R3me2a
ab179301	22.70941771	ab179301 - H4 S1p + K5ac + K8ac + K12ac + K16ac
ab179302	21.50264167	ab179302 - H4 R3cit
ab179303	23.55241979	ab179303 - H4 R3me1 + K5ac + K8ac + K12ac + K16ac + K20ac
ab179304	26.19058854	ab179304 - H4 R3me2s
ab179305	24.62583021	ab179305 - H4 R3me2s + K5ac + K8ac + K12ac + K16ac + K20ac
ab179306	25.07174896	ab179306 - H4 R3me2a
ab179307	21.95333125	ab179307 - H4 R3me2a + K5ac
ab179308	26.60351667	ab179308 - H4 R3me2a + K5ac + K8ac + K12ac + K16ac + K20ac
ab179309	21.79237396	ab179309 - H4 K5ac
ab179310	25.22160938	ab179310 - H4 K5ac + K8ac
ab179311	21.64852083	ab179311 - H4 K5ac + K12ac
ab179312	30.02351563	ab179312 - H4 K5ac + K16ac
ab179313	20.65655938	ab179313 - H4 K5ac + K8ac + K12ac
ab179314	33.51526979	ab179314 - H4 K5ac + K8me1 + K12ac
ab179315	27.10750625	ab179315 - H4 K5ac + K8ac + K12ac + K16ac
ab179316	27.93181146	ab179316 - H4 K5ac + K8me1 + K12ac + K16ac
ab179317	31.61876458	ab179317 - H4 K5ac + K8ac + K12me1 + K16ac
ab179318	23.06190938	ab179318 - H4 K5me1
ab179319	23.47719479	ab179319 - H4 K5me1 + K8ac + K12me1
ab179320	25.12428021	ab179320 - H4 K5me1 + K8me1 + K12me1
ab179321	24.74863333	ab179321 - H4 K5me1 + K8ac + K12ac + K16ac
ab179322	25.52731458	ab179322 - H4 K5me2
ab179323	23.23993958	ab179323 - H4 K5me3
ab179324	24.15354167	ab179324 - H4 K5cr
ab179325	22.32519167	ab179325 - H4 K8ac + K12ac
ab179326	22.63799375	ab179326 - H4 K8ac + K16ac
ab179327	25.3257625	ab179327 - H4 K8cr
ab179328	21.14718021	ab179328 - H4 K8me1
ab179329	22.75163021	ab179329 - H4 K12ac
ab179330	22.9851375	ab179330 - H4 K12ac + K16ac
ab179331	26.46762917	ab179331 - H4 K12ac + K16ac + K20me2
ab179332	24.16223125	ab179332 - H4 K12ac + K16ac + K20me3
ab179333	25.75001771	ab179333 - H4 K5mutQ + K8mutQ + K12ac + K16mutQ + K20mutQ
ab179334	24.30459688	ab179334 - H4 K12cr
ab179335	25.60109688	ab179335 - H4 K12fo
ab179336	21.51937396	ab179336 - H4 K12me1
ab179337	23.39520313	ab179337 - H4 K12me2
ab179338	22.20632813	ab179338 - H4 K12me3
ab179339	23.87375625	ab179339 - H4 K16ac
ab179340	20.92573542	ab179340 - H4 K16ac + K20ac
ab179341	21.95750729	ab179341 - H4 K16cr
ab179342	24.66406354	ab179342 - H4 K16me1
ab179343	18.58925625	ab179343 - H4 K16me3
ab179344	24.69743646	ab179344 - H4 R17me1
ab179345	26.46142396	ab179345 - H4 R17me2s
ab179346	28.64511354	ab179346 - H4 R17me2a
ab179347	23.61065833	ab179347 - H4 R19me1
ab179348	21.07713646	ab179348 - H4 R19me2s
ab179349	25.27621875	ab179349 - H4 R19me2a
ab179350	23.34781979	ab179350 - H4 K20ac
ab179351	22.53856042	ab179351 - H4 R23me1
ab179352	23.10151042	ab179352 - H4 R23me2s
ab179353	21.819925	ab179353 - H4 R23me2a
ab179354	23.10463229	ab179354 - H4 unmod(24-54)
ab179355	23.97440938	ab179355 - H4 K31fo
ab179356	20.92620625	ab179356 - H4 K31me1
ab179357	24.04422917	ab179357 - H4 K31me2
ab179358	19.51737292	ab179358 - H4 K31me3
ab179359	27.98632083	ab179359 - H4 P32hyd
ab179360	22.40721563	ab179360 - H4 R35me1
ab179361	26.01731146	ab179361 - H4 R35me2s
ab179362	24.10889688	ab179362 - H4 R35me2a
ab179363	24.73289375	ab179363 - H4 S47O-ac
ab179364	24.03245417	ab179364 - H4 S47GlcNAc
ab179365	23.5614125	ab179365 - H4 S47p
ab179366	25.34489063	ab179366 - H4 unmod(45-75)
ab179367	24.64602813	ab179367 - H4 Y51hyd
ab179368	29.3032375	ab179368 - H4 R55me1
ab179369	29.79409479	ab179369 - H4 R55me2s
ab179370	22.27138021	ab179370 - H4 R55me2a
ab179371	26.37703438	ab179371 - H4 K59fo
ab179372	20.78781354	ab179372 - H4 K59me1

ab179373	19.68897917	ab179373 - H4 K59me2
ab179374	22.74469167	ab179374 - H4 R67me1
ab179375	21.06574688	ab179375 - H4 R67me2s
ab179376	19.47133542	ab179376 - H4 R67me2a
ab179377	20.46855417	ab179377 - H4 unmod(71-102)
ab179378	21.70653229	ab179378 - H4 K77me1
ab179379	20.78488125	ab179379 - H4 K77me2
ab179380	22.4091875	ab179380 - H4 K77me3
ab179381	24.76293542	ab179381 - H4 K79fo
ab179382	23.42055104	ab179382 - H4 K79me1
ab179383	22.54787708	ab179383 - H4 K79me3
ab179384	24.43902708	ab179384 - H4 T82O-ac
ab179385	26.19659063	ab179385 - H4 Y88hyd
ab179386	27.97651042	ab179386 - H4 Y88nitro
ab179387	24.29722188	ab179387 - H4 K91fo
ab179403	21.23151563	ab179403 - H2A K5ac
ab179404	24.14152396	ab179404 - H2A K9me1
ab18504	18.724775	ab18504 - H2B S32p
ab19828	20.30288125	ab19828 - ScH2A S129p
ab20631	22.06046354	ab20631 - H3 R26cit
ab20635	19.58199688	ab20635 - H3 T3p + K4me3
ab21043	21.52949688	ab21043 - H3 K37me1
ab21998	19.11721042	ab21998 - H1.4 K25me2
ab22079	18.61682292	ab22079 - H4 K59me3
ab22080	19.15758854	ab22080 - H3 K122me2
ab22082	21.46463333	ab22082 - H3 K122me3
ab22214	22.53802917	ab22214 - H3 K37me2
ab22398	30.92825313	ab22398 - H2A R3me2a
ab22399	30.36653646	ab22399 - H2A R3me2s
ab23539	20.34239896	ab23539 - H3 T22p
ab24003	19.69046354	ab24003 - H3 K18ac
ab24404	19.23385104	ab24404 - H3 K27ac
ab24417	19.38628542	ab24417 - H3 K37me3
ab24444	21.93716458	ab24444 - H3 T11p
ab24659	18.71306563	ab24659 - H2A K127me3
ab24660	20.57188021	ab24660 - H2A K125me3
ab26307	21.6656375	ab26307 - H3 T45p
ab26342	21.49418958	ab26342 - H2B K5me2
ab26739	24.51748021	ab26739 - H4 T30p
ab27896	27.01396563	ab27896 - H4 T96p
ab27898	22.14634271	ab27898 - H4 T80p
ab28756	29.06278125	ab28756 - H2A R3me1
ab32876	24.591075	ab32876 - H3 R2cit + R8cit + R17cit
ab33047	38.66283229	ab33047 - H2A K127me2
ab33505	22.82052188	ab33505 - H3 T118p
ab33594	24.91965208	ab33594 - H3 R128me2a
ab34466	23.21849479	ab34466 - H3 K122ac
ab36927	18.62439063	ab36927 - H2A S122p
ab4555	21.86023438	ab4555 - H3 K79me1
ab4556	23.00338333	ab4556 - H3 K79me2
ab4557	21.52213958	ab4557 - H3 K79me3
ab4560	22.03589479	ab4560 - H4 K79me2
ab46854	16.85520208	ab46854 - H3 K18me3
ab48359	19.37908125	ab48359 - H3 K23ac
ab53244	22.19743542	ab53244 - H3 K64me1
ab54016	41.72969063	ab54016 - H2A K9ac
ab5499	16.90144375	ab5499 - H3 S28p
ab7768	21.98645729	ab7768 - H3 K4me2
ab89018	27.12613854	ab89018 - H1.4 S26p
ab92374	23.11838438	ab92374 - H3 K4me3 (HEX1)
ab95444	25.3893625	ab95444 - H4 K16me2