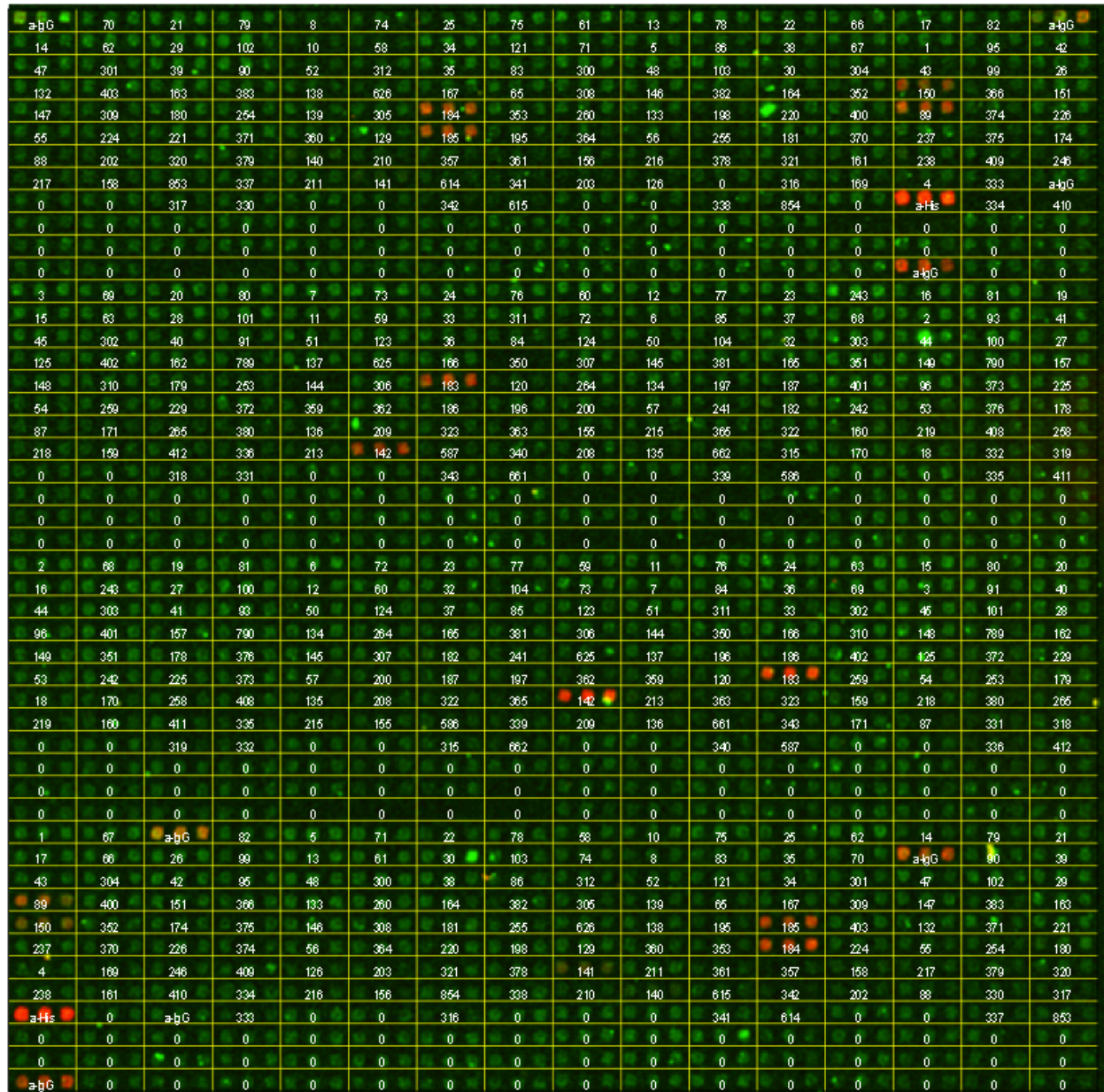


Analysis of Histone H3R8me2s Antibody on the EpiTitan™ Histone Peptide Array



Antibody Specificity:

EpiTitan™ Histone Peptide Array was used to analyze binding specificity of the Histone H3 R8Me2s Antibody (Arg8 symmetric dimethyl) antibody at 1:5,000. Antibody signal appears in red, while the spotting tracer appears in green. The array image is overlaid with a frame containing numbers corresponding to the peptide number (“Peptide #”) in the raw data list below. Raw data “Signal” is an average of the antibody signal intensities of all 6 instances of the corresponding peptide on the array (2 instances of three peptide spots in a row).

Peptide #	Signal	STDEV	Peptide Name
142	66654	25177	H3R8me2s
183	66009	11545	H3R8me2s + K9me3
184	59167	4254	H3R8me2s + K9me2
185	50324	2091	H3R8me2s + K9me1
89	48129	9765	H3K4me3 + R8me2s + K9me3
150	35690	4424	H3R17me2s
90	20422	28299	H3 (15-41)
141	12906	11122	H3R8me2a
140	6546	703	H3R8me1
182	5112	1164	H3R8me2a + K9me1
178	2656	2219	H3R8me1 + K9me3
159	2577	1898	H3R17me2a (1-25)
180	2250	1927	H3R8me2a + K9me3
179	2092	468	H3R8me1 + K9me2
11	2068	1255	H3K14ac + K18ac
160	1950	829	H3R17me2a + K18ac (1-25)
181	1917	949	H3R8me2a + K9me2
213	1864	2201	H3 120-135
61	1772	131	H3R2me2s + K4me2
32	1717	1435	H3K4me2
74	1687	634	H4R3me2s
158	1628	1076	H3K18ac (1-25)
2	1437	510	H3K14ac
19	1377	1454	H3K4me3 + K9ac
13	1329	595	H3K4ac + K18ac
238	1314	1538	H3K9me2 + K27me1 (1-32)
40	1305	284	H3R2me2a + K4me3 + S10p
24	1287	482	H3K4me3 + K14ac + K18ac
151	1261	115	H3R17me1
365	1259	1628	H4K12ac + K20ac
63	1188	1162	H3Cit2 + K4me2
10	1184	645	H3K18ac
353	1132	870	H4S1p + K5ac + K8ac + K12ac + K16ac
125	1126	1056	H3T3p
5	1104	386	H3K4ac + K14ac
149	1085	223	H3R17me2a
1	1079	405	H3 (1-20)
38	1073	663	H3K4me3 + S10p
60	1022	668	H3R2me2a + K4me2