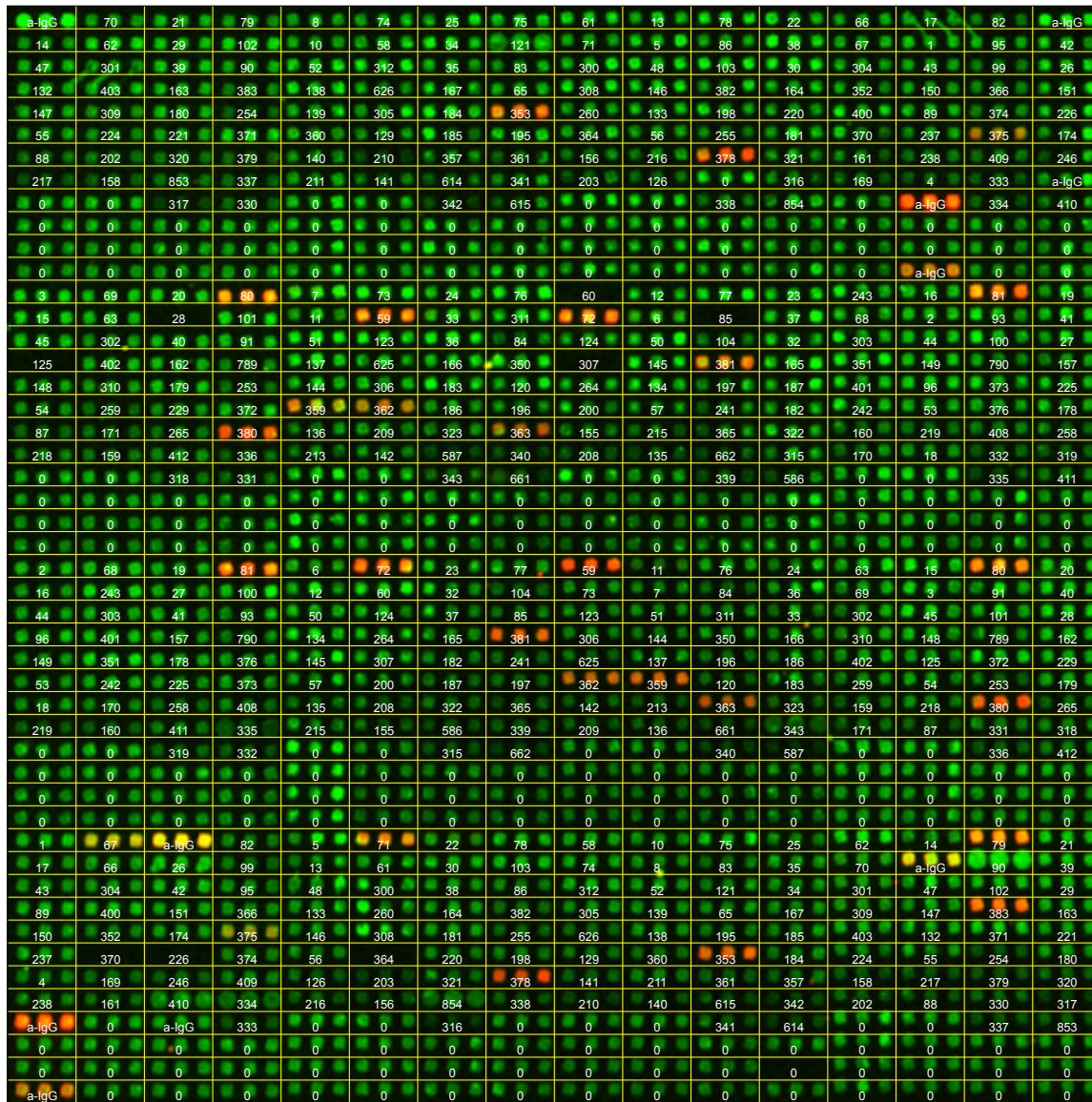


Analysis of anti-H4K8Ac (Cat. No. 13-0021) on the *EpiTitan™* Histone Peptide Array



Data Specificity:

EpiTitan™ Histone Peptide Array was used to analyze binding specificity of the Histone H4K8Ac Antibody (Catalog No. 13-0021) 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” shown in Table 1 is an average of the antibody signal intensities of all instances of the corresponding peptide on the array. Only peptides exhibiting signals that can be visualized on the array were listed in Table 1.

Result:

Anti-H4K8ac (**E1615**) recognizes all the peptides containing H4K8ac on the array.

Peptide #	Signal	STDEV	Peptide Name
81	198895	1342	H4R3me1 + K5ac + K8ac + K12ac + K16ac + K20ac
72	164189	3083	H4K5ac + K8ac + K12ac
59	162096	6468	H4K5ac + K8ac + K12ac + K16ac
80	149138	21192	H4R3me2s + K5ac + K8ac + K12ac + K16ac + K20ac
380	115995	8118	H4K5ac + K8ac + K16ac
79	109705	152523	H4R3me2a + K5ac + K8ac + K12ac + K16ac + K20ac
378	108402	20183	H4K8ac + K12ac + K16ac
353	103450	29830	H4S1p + K5ac + K8ac + K12ac + K16ac
381	99796	10051	H4K5me1 + K8ac + K12ac + K16ac
67	83557	1682	H4K8ac
359	82639	1797	H4K5ac + K8ac
383	82234	114829	H4K5ac + K8ac + K12me1 + K16ac
362	75718	2589	H4K8ac + K12ac
363	61702	8261	H4K8ac + K20ac
375	56205	5051	H4K5me1 + K8ac + K12me1
71	53036	73398	H4K8ac + K16ac