

Analysis of Histone H3R8me1 Antibody on the EpiTitan™ Histone Peptide Array

a-IgG	70	21	79	8	74	25	75	61	13	78	22	66	17	82	a-IgG
14	62	29	102	10	58	34	121	71	5	86	38	67	1	95	42
47	301	39	90	52	312	35	83	300	48	103	30	304	43	99	26
132	403	163	383	138	626	167	65	308	146	382	164	352	150	366	151
147	309	180	254	139	305	184	353	260	133	198	220	400	89	374	226
55	224	221	371	360	129	185	195	364	66	255	181	370	237	375	174
88	202	320	379	140	210	357	361	156	216	378	321	161	238	409	246
217	158	853	337	211	141	614	341	203	126	0	316	169	4	333	a-IgG
0	0	317	330	0	0	342	615	0	0	338	854	0	a-His	334	410
0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0
0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0
0	0	0	0	0	0	0	0	0	0	0	0	0	a-IgG	0	0
3	69	20	80	7	73	24	76	60	12	77	23	243	16	81	19
15	63	28	101	11	59	33	311	72	6	85	37	68	2	93	41
45	302	40	91	51	123	36	84	124	50	104	32	303	44	100	27
125	402	162	789	137	625	166	350	307	145	381	165	351	149	790	157
148	310	179	253	144	306	183	120	264	134	197	187	401	96	373	225
54	259	229	372	359	362	186	196	200	57	241	182	242	53	376	178
87	171	265	380	136	209	323	363	155	215	365	322	160	219	408	258
218	159	412	336	213	142	587	340	208	135	662	315	170	18	332	319
0	0	318	331	0	0	343	661	0	0	339	586	0	0	335	411
0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0
0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0
0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0
2	68	19	81	6	72	23	77	59	11	76	24	63	15	80	20
16	243	27	100	12	60	32	104	73	7	84	36	69	3	91	40
44	303	41	93	50	124	37	85	123	51	311	33	302	45	101	28
96	401	157	790	134	264	165	381	306	144	350	166	310	148	789	162
149	351	178	376	145	307	182	241	625	137	196	186	402	125	372	229
53	242	225	373	57	200	187	197	362	369	120	183	239	54	253	179
18	170	258	408	135	208	322	365	142	213	363	323	159	218	380	265
219	160	411	335	215	155	586	339	209	136	661	343	171	87	331	318
0	0	319	332	0	0	315	662	0	0	340	587	0	0	336	412
0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0
0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0
0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0
0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0
1	67	a-IgG	82	5	71	22	78	58	10	75	25	62	14	79	21
17	66	26	99	13	61	30	103	74	8	83	35	70	a-IgG	90	39
43	304	42	95	48	300	38	86	312	52	121	34	301	47	102	29
89	400	151	366	133	260	164	382	305	139	65	167	309	147	383	163
150	352	174	375	146	308	181	255	626	138	195	185	403	132	371	221
237	370	226	374	56	364	220	198	129	360	353	184	224	55	254	180
4	169	246	409	126	203	321	378	141	211	361	357	158	217	379	320
238	161	410	334	216	156	854	338	210	140	615	342	202	88	330	317
a-His	0	a-IgG	333	0	0	316	0	0	0	341	614	0	0	337	853
0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0
0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0
a-IgG	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0

Antibody Specificity:

EpiTitan™ Histone Peptide Array was used to analyze binding specificity of the Histone H3 R8Me1 Antibody (Arg8 monomethyl) 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
140	34782	9185	H3R8me1
178	33375	4946	H3R8me1 + K9me3
179	28414	8438	H3R8me1 + K9me2
11	9117	251	H3K14ac + K18ac
160	8400	4959	H3R17me2a + K18ac (1-25)
158	8090	4813	H3K18ac (1-25)
2	5973	1307	H3K14ac
14	4827	2572	H3K9ac + K14ac + K18ac
159	4627	2031	H3R17me2a (1-25)
24	4448	1081	H3K4me3 + K14ac + K18ac
10	4269	1387	H3K18ac
53	4196	3532	H3Cit2
61	3720	3046	H3R2me2s + K4me2
217	3573	1604	H3K14-Nle (1-20)
149	3302	1992	H3R17me2a
12	3247	901	H3K9ac + K18ac
1	3159	650	H3 (1-20)
63	3134	31	H3Cit2 + K4me2
60	3115	1979	H3R2me2a + K4me2
6	3107	595	H3K9ac + K14ac
238	3006	3343	H3K9me2 + K27me1 (1-32)
135	2964	319	H3K3K4me1 + K18ac
210	2950	1768	H3.3 30-49
258	2749	313	H3K9me2 (1-15)
260	2566	638	H3K4me1 + K9me2
23	2541	15	H3K4me3 + K9ac + K18ac
22	2526	671	H3K4me3 + K9ac + K14ac
156	2513	2585	H3K14me3
161	2407	1877	H3 (1-25)
3	2392	1283	H3K9ac
33	2378	563	H3K4me2 + K9ac + K14ac + K18ac
19	2374	149	H3K4me3 + K9ac
13	2295	1558	H3K4ac + K18ac
62	2273	1013	H3R2me1 + K4me2
21	2229	985	H3K4me3 + K18ac
25	2113	19	H3K4me3 + K9ac + K14ac + K18ac
150	2067	2018	H3R17me2s
133	2055	858	H3K9me2