

Analysis of anti-H3R17Me2a 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	145	382	164	352	150	365	151
147	309	180	264	139	305	184	353	260	133	198	220	400	89	374	226
55	224	221	371	360	129	185	195	364	56	255	181	370	237	375	174
68	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-GST
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	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	26	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	263	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	180	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
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	359	120	183	259	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	616	342	202	88	330	317
a-His	0	a-GST	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

Data Specificity:

EpiTitan™ Histone Peptide Array was used to analyze binding specificity of the H3R17me2a 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).

Result:

Anti-H3R17me2a only recognizes H3R17me2a peptides on the array.

Peptide #	Signal	STDEV	Peptide Name
160	343669	178002	H3R17me2a + K18ac (1-25)
159	273965	96587	H3R17me2a (1-25)
149	185544	60487	H3R17me2a
151	14390	11443	H3R17me1
179	8466	11409	H3R8me1 + K9me2
218	5823	7678	H3K18-Nle (11-26)
411	4343	5750	H2BK20ac
60	3902	480	H3R2me2a + K4me2
219	3556	1335	H3K23-Nle (15-31)
351	3258	4563	H4 (1-23) (no N-ac)
141	3256	133	H3R8me2a
59	3000	4242	H4K5ac + K8ac + K12ac + K16ac
75	2845	3639	H4R3me1
47	2395	2355	H3R2me2a
412	2028	1997	H2BK12ac + K15ac + K16ac + K20ac
147	1979	661	H3K9me1 + S10p
2	1791	2402	H3K14ac
403	1508	2132	H2BK5me1
48	1268	1329	H3R2me2a + K4ac + K9ac + K14ac + K18ac
129	1085	666	H3K9me2 + K27me2
11	914	1293	H3K14ac + K18ac
260	857	860	H3K4me1 + K9me2
28	856	215	H3R2me2a + T3p + K4me3 + K9ac + K14ac + K18ac
43	845	1086	H3K4ac + K9me3 + K14ac + K18ac
343	837	853	H4K12acK16acK20me3 (1-25)
18	728	531	H3K4me3
65	710	1004	H3K4N3
335	683	735	H3K36ac (21-44)
361	665	790	H4K5ac + K20ac
614	642	411	cnp1 (1-23)
331	632	825	H3 (21-44)
178	582	340	H3R8me1 + K9me3
302	559	790	H2AK5ac
30	513	726	H3R2me2a + K4me3
332	512	724	H3K36me1 (21-44)
55	486	648	H3Cit2 + K4me3 + K9ac + K14ac + K18ac
408	483	249	H2BK12ac
195	481	622	H3K27me3

51	480	679	H3R2me1 + K4me3
68	467	661	H4K12ac
376	459	433	H4K5me1 + K8me1 + K12me1
140	454	643	H3R8me1
242	428	606	H3K27ac + S28p
174	428	100	H3R2me2s + K4me3 + K9ac + K14ac + K18ac
135	427	493	K3K4me1 + K18ac
402	416	588	H2BK5me2
44	407	575	H3K4me2 + K9ac + K18ac
340	406	575	H2B (108-125)
357	385	544	H4S1p (1-23)
258	356	235	H3K9me2 (1-15)
50	349	432	H3R2me2a + K4me3 + K9ac + K14ac + K18ac
208	348	492	H3 105-124
330	348	64	H3Cit8 (1-21)
166	345	488	H3K4me3 + T6p + K9ac + K14ac + K18ac
62	341	483	H3R2me1 + K4me2
180	340	422	H3R8me2a + K9me3
134	330	29	H3K9me1
104	329	465	H3 (74-84)
255	327	416	H3K56me3 (52-61)
360	323	247	H4K5ac + K16ac
77	318	450	H4S1p + R3me2s
26	314	443	H3T3p + K4me3 + K9ac + K14ac + K18ac
359	308	435	H4K5ac + K8ac
211	296	418	H3.3 75-94
197	291	223	H3K27me1
123	290	410	H3K36ac (27-45)
150	287	111	H3R17me2s
790	283	400	H3S31p + K36me3
215	280	395	H3K23me1 15-34
198	276	391	H3R26me2a + K27me3
88	274	193	H4K12ac + K16ac
162	270	382	H3T6p
307	263	372	H2AS1p + Cit3 + K5ac
259	262	51	H3K4me2 + K9me2
63	258	365	H3Cit2 + K4me2
181	248	296	H3R8me2a + K9me2
126	248	221	H3.3K36me1
381	241	341	H4K5me1 + K8ac + K12ac + K16ac
350	230	325	H4R3me2a + K5ac
185	230	316	H3R8me2s + K9me1
352	228	322	H4K20ac

254	222	313	H3K56ac (52-61)
187	216	134	H3K4ac + K9me1 + K14ac + K18ac
626	212	300	H2A.X K5ac
13	211	299	H3K4ac + K18ac
90	207	120	H3 (15-41)
85	206	291	H4K12ac + K16ac + K20me3
210	196	277	H3.3 30-49
52	195	276	H3R2me1 + K4me3 + K9ac + K14ac + K18ac
138	194	274	H3K18me2
220	188	265	H3T6p + K9me3
42	184	261	H3K9me3
35	184	260	H3K4me1 + K9ac + K14ac + K18ac
4	179	253	H3K4ac
38	173	244	H3K4me3 + S10p
80	168	238	H4R3me2s + K5ac + K8ac + K12ac + K16ac + K20ac
370	168	64	H4K12ac + KQ5,8,16,20
86	164	232	H4K12ac + K16ac + K20me2
382	163	68	H4K5ac + K8me1 + K12ac + K16ac
136	162	230	H3T11p
322	157	99	H2A.Z K4ac + K8ac + K12ac (1-19)
89	151	214	H3K4me3 + R8me2s + K9me3
156	147	208	H3K14me3
364	145	205	H4K12ac + K16ac
305	142	201	H2AS1p + R3me2a + K5ac
401	141	181	H2BK5me3
306	140	198	H2ACit3 + K5ac
146	140	56	H3K9me2 + S10p
341	132	187	H4 (1-25)
53	132	153	H3Cit2
661	130	184	H3K4me3 (1-11)
171	129	183	H3K23ac
17	129	183	H3K4ac + K9ac + K14ac + K18ac
336	127	179	H3K27acK36me1 (21-44)
54	124	176	H3Cit2 + K4me3
338	123	174	H3K27acK36me3 (21-44)
84	115	163	H4K20me1
184	115	163	H3R8me2s + K9me2
36	113	160	H3S10p
40	112	158	H3R2me2a + K4me3 + S10p
56	110	155	H3Cit2 + K4ac + K9ac + K14ac + K18ac
93	109	154	H3K36me3
1	108	153	H3 (1-20)
125	102	144	H3T3p

209	100	142	H3.3 15-34
27	97	44	H3T3p + K4me3
70	94	134	H4K5ac + K12ac
23	92	130	H3K4me3 + K9ac + K18ac
789	88	124	H3K36me3
165	87	124	H3T6p + K9ac + K14ac + K18ac
155	87	106	H3K14me2
95	83	75	H3K18me3 + K36me3
225	80	2	H3K27me3 + S28p
121	80	113	H3K36me2 (27-45)
342	79	112	H4K20me3 (1-23)
120	77	109	H3K36me3 (27-45)
15	75	2	H3K4ac + K14ac + K18ac
301	74	104	H2AK5ac + K9ac + K13ac + K15ac
300	68	96	H2A (1-17)
304	64	91	H2AR3me2a + K5ac
45	63	89	H3K4me1 + K9ac + K18ac
362	57	81	H4K8ac + K12ac
61	56	80	H3R2me2s + K4me2
124	56	80	H3 (27-45)
144	48	68	H3K9ac + S10p
99	44	62	H4 (11-27)
186	42	43	H3K4ac + K9me2 + K14ac + K18ac
7	42	59	H3K4ac + K9ac
145	40	57	H3K9me3 + S10p
353	37	52	H4S1p + K5ac + K8ac + K12ac + K16ac
87	34	48	H4K12ac + K16ac + K20me1
8	33	47	H3K4ac + K9ac + K14ac
303	33	4	H2AS1p + K5ac
71	30	42	H4K8ac + K16ac
323	29	41	H2A.Z K4ac + K8ac + K12ac (1-19) N-ac
5	27	39	H3K4ac + K14ac
372	27	39	H4K8me1
237	27	38	H3K9me2 (1-32)
58	20	28	H4 (1-23)
400	20	28	H2B (1-24)
321	20	28	H2A.Z (1-19) N-ac
196	17	12	H3K27me2
6	16	23	H3K9ac + K14ac
164	16	9	H3K4me2 + T6p
308	16	23	H2AS1p + K5ac + K9ac + K13ac + K15ac
182	10	14	H3R8me2a + K9me1
625	10	14	H2A.X (1-17)

34	7	10	H3K4me1
253	7	10	H3 (52-61)
83	7	9	H4K20me2
20	5	7	H3K4me3 + K14ac
67	1	1	H4K8ac