

Analysis of Histone H3R8me2a Antibody on the EpiTitan™ Histone Peptide Array

a-bG	70	21	79	8	74	25	75	61	13	78	22	66	17	82	a-bG
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	406	163	383	138	626	167	65	308	146	382	184	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	56	255	181	370	237	375	174
88	202	320	379	140	210	357	361	156	216	378	321	161	238	400	246
217	158	853	337	211	141	614	341	203	126	0	316	169	4	333	a-bG
0	0	317	330	0	0	342	615	0	0	338	854	0	a-Hs	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	a-bG	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
46	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	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
0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0
2	68	19	81	6	72	23	77	99	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	166	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-bG	82	5	71	22	78	88	10	75	25	62	14	79	21
17	66	26	99	13	61	30	103	74	8	83	35	70	a-bG	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	189	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-Hs	0	a-bG	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
0	0	0	0	0	0	0	0	0	0	0	0	0	0	0	0
a-bG	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 R8Me2a Antibody (Arg8 asymmetric 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
141	90297	5689	H3R8me2a
182	78458	15500	H3R8me2a + K9me1
180	70367	1146	H3R8me2a + K9me3
181	63333	9334	H3R8me2a + K9me2
96	50689	1980	H3K4me3 + R8me2a + K9me3
221	33511	5269	H3T6p + R8me2a + K9me3
224	21900	2724	H3R26me2a
159	21699	1631	H3R17me2a (1-25)
140	19088	6255	H3R8me1
149	16365	2406	H3R17me2a
178	15681	110	H3R8me1 + K9me3
200	14781	520	H3R26me2a + K27me1
179	11999	654	H3R8me1 + K9me2
198	9131	1700	H3R26me2a + K27me3
183	8748	5365	H3R8me2s + K9me3
155	7866	9556	H3K14me2
73	6652	1353	H4R3me2a
160	6410	2934	H3R17me2a + K18ac (1-25)
158	6356	4837	H3K18ac (1-25)
142	4794	486	H3R8me2s
156	4789	5328	H3K14me3
184	4722	1637	H3R8me2s + K9me2
61	4429	5897	H3R2me2s + K4me2
211	4375	5861	H3.3 75-94
11	3853	386	H3K14ac + K18ac
185	3815	1615	H3R8me2s + K9me1
241	3505	4660	H3K27ac
10	3469	377	H3K18ac
614	3404	4598	cnp1 (1-23)
12	3173	403	H3K9ac + K18ac
401	2738	3506	H2BK5me3
2	2716	260	H3K14ac
6	2552	87	H3K9ac + K14ac
47	2525	404	H3R2me2a
217	2409	1029	H3K14-N1e (1-20)
3	2258	705	H3K9ac
57	2210	901	H3T3p
246	2206	1979	H3 (1-20) N-ac
216	2056	2211	H3K9-N1e (1-20)
53	2051	990	H3Cit2