

<b>OmicsArray™ Allergen Antigen Array Antigen List - Catalog No. PA006</b>			
<b>Allergen#</b>	<b>ID</b>	<b>Allergen Description</b>	<b>Category</b>
1	Cat Dander	Cat Dander	animal Allergen
2	Cat Epi	Cat Epithelia Felis catus (domesticus)	animal Allergen
3	Dog Dander	Dog Dander	animal Allergen
4	Dog EPI	Dog Epithelia	animal Allergen
5	Dog Hair Canis lupus familiaris	Dog Hair Canis lupus familiaris	animal Allergen
6	Mite, Food/Storage Acarus siro	Mite, Food/Storage Acarus siro	environment allergen
7	Mite, Food/Storage Lepidoglyphus destructor	Mite, Food/Storage Lepidoglyphus destructor	environment allergen
8	Mite, House Dust Glycyphagus domesticus	Mite, House Dust Glycyphagus domesticus	environment allergen
9	Ara h 1.0101	Arachis hypogaea (peanut)	food allergen
10	Ara h 2.0201	peanut allergen	food allergen
11	Ara h 3.0101	11 S globulin from peanut	food allergen
12	Ara h 6.0101	conglutin, 2S albumin protein family	food allergen
13	Ara h 8.0101	a Bet v 1-homologous panallergen	food allergen
14	Ara h 9.0101	The lipid transfer protein (LTP)	food allergen
15	Banana Musa acuminata	Banana Musa acuminata	food allergen
16	Beef taurus	Beef, Bos taurus	food Allergen
17	Cor a 1.0401	allergen from Corylus avellana	food allergen
18	Crab Callinectes sapidus	Crab Callinectes sapidus	food allergen
19	CSN1	Casein (milk protein) / CSN1	food allergen
20	Gliadin	Gliadin (recombinant; deamidated)	food allergen
21	Gly m 4.0101	soybean protein	food allergen
22	Gly m 5.0101	Glycine max seed	food allergen
23	HEL	Egg-white Lysozyme (HEL)	food allergen
24	Honey Bee	Honey Bee_Apis mellifera	food allergen
25	Lamb Ovis aries	Lamb Ovis aries	food allergen
26	Lobster Homarus americanus	Lobster Homarus americanus	food allergen
27	Mal d 1.0108	Major apple allergen Mal d 1	food allergen
28	Milk, Bovine Bos taurus	Milk, Bovine Bos taurus	food allergen
29	NEW Tri a 14.0101	a major wheat (Triticum aestivum) allergen	food allergen
30	Oats, Common/Cultivated Avena sativa	Oats, Common/Cultivated Avena sativa	food allergen
31	OVA	Ovalbumin (OVA)	food allergen
32	Peanut	Peanut_Arachis hypogaea	food allergen
33	Pen a 1.0101	allergen in brown shrimp Penaeus aztecus	food allergen
34	Pork Sus scrofa	Pork Sus scrofa	food allergen
35	Salmon Salmo salar	Salmon Salmo salar	food allergen
36	Shrimp	Shrimp_Penacidae	food allergen
37	Alternaria alternata, Alternaria	Alternaria alternata Alternaria tenuis	fungus
38	Aspergillus flavus	Aspergillus flavus	fungus
39	Art v 1.0101	Weed allergen of mugwort pollen	Grass allergen
40	Bahia Paspalum notatum	Bahia Paspalum notatum	Grass allergen
41	Bermuda	Bermuda Cynodon dactylon	Grass allergen

42	Phl p 1.0102	pollen allergen of Phleum pratense	Grass allergen
43	Phl p 12.0101	grass pollen allergen	Grass allergen
44	Phl p 2.0101	group II/III of grass pollen allergen	Grass allergen
45	Phl p 5.0203	group V allergen-starch granules of Phleum pratense pollens	Grass allergen
46	Phl p 6.0101	pollen allergen of Phleum pratense	Grass allergen
47	Phl p 7.0101	a minor, non-glycosylated, allergen of timothy grass pollen	Grass allergen
48	Ant, Fire Solenopsis invicta	Ant, Fire Solenopsis invicta	insect Allergen
49	CockRoACH	Cockroach, American Periplaneta	insect Allergen
50	Dermatophagoides farinae	Mite, Dermatophagoides farinae	mite allergen
51	Mite, Dermatophagoides	Dermatophagoides pteronyssinus	mite allergen
52	NEW Cor a 9 (non recombinant)	an 11S legumin-like seed storage globulin and a major allergen from Corylus avellana	plant allergen
53	Cor a 14.0101	Seed storage proteins, a major allergens in several other tree nuts and peanut	plant food allergen
54	Wheat	Wheat, Whole Triticum aestivum	plant food allergen
55	Mal d 3.0101	major apple (Malus domestica) allergen	plant food allergen
56	NEW Cor a 8.0101	allergens in fruits and tree	plant food allergen
57	Aln g 1.0101	pollen of Alnus glutinosa	Tree allergen
58	Aln g 4.0101	polcalcin Aln g 4, calcium-binding pollen allergen Aln g 4	Tree allergen
59	Ash, White Fraxinus americana	Ash, White Fraxinus americana	Tree allergen
60	Bet v 1.0101	major pollen allergen	Tree allergen
61	Bet v 2.0101	profilin-1, pollen of Betula verrucosa	Tree allergen
62	Bet v 4.0101	allergen for plant polysensitization	Tree allergen
63	Bet v 6.0102	birch pollen allergen	Tree allergen
64	Car b 1.0109	pollen allergen of Carpinus betulus	Tree allergen
65	Cedar	Cedar, Red Juniperus virginiana	tree Allergen
66	Cor a 1.0103	pollen allergen from Corylus avellana	Tree allergen
67	Fra e 1.0101	Tree allergen Ole e 1-like protein family	Tree allergen
68	Oak, Western White Quercus garryan	Oak, Western White Quercus garryana	Tree allergen
69	Plantain	Plantain, English Plantago lanceolata	weeds allergen
70	C. Albicans	Candida Albicans (wild strain)	fungus
71	anti-Ig control 1	control	internal control
72	anti-Ig control 2	control	internal control
73	anti-Ig control 3	control	internal control
74	anti-Ig control 4	control	internal control
75	Ig control 1	control	internal control
76	Ig control 2	control	internal control
77	Ig control 3	control	internal control
78	Ig control 4	control	internal control