

OmicsArray™ Allergen Antigen Array - Catalog No. PA006-V2				
Serial#	Allergen Name	Allergen Code	Category	Description
1	D. farinae	D2	In-door_House dust mite	House dust extract
2	Blomia tropicalis	D201	In-door_House dust mite	House dust mite (Blomia tropicalis)
3	Acarus siro	D70	In-door_Flour mite	Mite, Food/Storage Acarus siro
4	Lepidoglyphus destructor	D71	In-door_Food/Storage Mite	Mite, Food/Storage Lepidoglyphus destructor
5	Glycyphagus domesticus	D73	In-door_House dust mite	Mite, House Dust Glycyphagus domesticus (Acarus domesticus)
6	Cat dander	E1	In-door_Pets allergen	Cat dander (e1)
7	Cat Epithelia	E1	In-door_Pets allergen	Cat Epithelia Felis catus (domesticus)
8	Dog Hair, canis familiaris	E5	In-door_Pets allergen	Dog Hair Canis lupus familiaris
9	Dog dander (skin flaks/extract)	E5	In-door_Pets allergen	dog skin flakes/extract
10	Dog dander (epithelia extract)	E5	In-door_Pets allergen	Mixed-Breed Dog Epithelia
11	Milk	F2	Food allergen	Milk, Bovine Bos taurus
12	Cod fish	F3	Food allergen	Codfish
13	Wheat	F4	Food allergen	Wheat, Whole Triticum aestivum
14	Rye	F5	Food allergen	Rye_Cereal
15	Barley	F6	Food allergen	Barley
16	Maize/Corn	F8	Food allergen	Corn
17	Sesame seed	F10	Food allergen	Sesame seed
18	Peanut	F13	Food allergen	Peanut
19	Soybean	F14	Food allergen	Soybean (f14)
20	Almond	F20	Food allergen	Almond
21	Tomato	F25	Food allergen	Tomato
22	Pork	F26	Food allergen	Pork Sus scrofa
23	Beef	F27	food allergen	Beef extract
24	Strawberry	F44	Food allergen	Strawbeery extract
25	Apple (Mal d 3)	F49	Food allergen	Apple (Mal d 3.0101)
26	Egg yolk	F75	Food allergen	Egg yolk (f75)
27	Milk casein	F78	Food allergen	Milk casein CSN1
28	Chicken	F83	Food allergen	Chicken
29	Lamb	F88	Food allergen	Lamb Ovis aries
30	Banana	F92	Food allergen	Banana Musa acuminata
31	Wheat gliadin	F98	Food allergen	Gliadin (recombinant; deamidated)
32	Oats	F154	Food allergen	Oats, Common/Cultivated Avena sativa
33	Rice	F200	Food allergen	Rice
34	Pecan nut	F201	Food allergen	Pecan nuts extract
35	Cashew nut	F202	Food allergen	Cashew Nut
36	Pistachio Nut	F203	Food allergen	Pistachio Nut
37	Clam	F207	Food allergen	Clam
38	Ovalbumin	F232	Food allergen	Egg white (albumin)
39	Egg white lysozyme	F245	Food allergen	Egg white (lysozyme)
40	Walnut	F256	Food allergen	Walnut extract
41	Blueberry	F288	Food allergen	Blueberry extract
42	Flaxseed/Linseed	F333	Food allergen	Flaxseed (Linum usitatissimum)

43	Peanuts (Ara h 8)	F352	Food allergen	Major peanut allergen Ara h 8
44	Soybean (Gly m 4)	F353	Food allergen	Soybean Glycine max (Gly m) 4.0101 protein
45	Peanuts (Ara h 1)	F422	Food allergen	Major peanut allergen Ara h 1
46	Peanut (Ara h 2)	F423	Food allergen	Major peanut allergen Ara h 2
47	Peanuts (Ara h 3)	F424	Food allergen	Major peanut allergen Ara h 3
48	Hazel nut (Cor a 8)	F425	Food allergen	Hazelnut (Cor a 8.0101)
49	Peanuts (Ara h 9)	F427	Food allergen	Major peanut allergen Ara h 9
50	Hazelnut (Cor a 1.0401)	F428	Food allergen	Bet v 1-related major hazelnut allergen Cor a 1.0401
51	Hazel pollen (Cor a 1.0103)	F428	Food allergen	PR-10 protein, a major pollen in hazel plants (Corylus avellana)
52	Soybean (Gly m 5)	F431	Food allergen	Soybean (Glycine max) allergen
53	Wheat (Tri a 14)	F433	Food allergen	Wheat (Tri a 14.0101)
54	Apple (Mal d 1)	F434	Food allergen	Major apple protein Mal d 1.
55	Hazelnut (Cor a 14)	F439	Food allergen	Hazelnut (Cor a 14.0101)
56	Hazelnut (Cor a 9)	F440	Food allergen	Hazelnut (Cor a 9)
57	Peanuts (Ara h 6)	F447	Food allergen	Major peanut allergen Ara h 6
58	Crab	F23	Food allergen_seafood	Crab Callinectes sapidus
59	Shrimp	F24	Food allergen_seafood	Shrimp
60	Tuna	F40	Food allergen_seafood	Tuna fish
61	Salmon	F41	Food allergen_seafood	Salmon Salmo salar
62	Lobster	F80	Food allergen_seafood	Lobster Homarus americanus
63	Scallops	F338	Food allergen_seafood	Scallops (f338)
64	Shrimp (Pen a 1)	F351	Food allergen_seafood	Shrimp (Pen a 1.0101)
65	Penicillium chrysogenum	M1	Mold (fungus)	Penicillium chrysogenum
66	Aspergillus flavus	M228	Mold (fungus)	Aspergillus flavus
67	Aspergillus fumigatus	M3	Mold (fungus)	Aspergillus fumigatus (m3)
68	Candida Albicans	M5	Mold (fungus)	Candida Albicans (wild strain)
69	Alternaria alternata	M6	Mold (fungus)	Alternaria alternata, Alternaria tenuis
70	Honey Bee	I1	Insect allergen	Honey Bee_Apis mellifera
71	Cockroach	I6	Insect Allergen	Cockroach,American_Periplaneta americana
72	Fire Ant	I70	Insect Allergen	Ant, Fire Solenopsis invicta
73	Johnson grass	G10	Out-door_Grass allergen	Johnson grass
74	Bahia grass	G17	Out-door_Grass allergen	Bahia Paspalum notatum
75	Bermuda grass	G2	Out-door_Grass allergen	Bermuda grass pollen extract (Cynodon dactylon)
76	Timothy grass (Phl p 1)	G205	Out-door_Grass allergen	Allergen derived from timothy grass (Phelum pratense) pollen
77	Timothy grass (Phl p 2)	G206	Out-door_Grass allergen	Major allergen from pollen of Timothy grass (Phleum pratense).
78	Timothy grass (Phl p 6)	G209	Out-door_Grass allergen	Phl p 6 allergen from Timothy grass (Phleum pratense) pollen
79	Timothy grass (Phl p 7)	G210	Out-door_Grass allergen	Timothy grass (Phl p 7.0101)
80	Timothy grass (Phl p 12)	G212	Out-door_Grass allergen	Minor allergen in timothy grass (Phleum pratense) pollen
81	Timothy grass (Phl p 5)	G215	Out-door_Grass allergen	Major Timothy grass (Phleum pratense) pollen allergen
82	Perennial ryegrass	G5	Out-door_Grass allergen	Perennial ryegrass
83	Timothy grass	G6	Out-door_Grass allergen	Timothy grass (g6)
84	Plantain	W9	Out-door_grass allergen	Plantain, English_Plantago lanceolata
85	Common ragweed	W1	Out-door_weeds allergen	Ragweed, Western, Ambrosia psilostachya
86	Rough pigweed	W14	Out-door_weeds allergen	Pigweed

87	Sheep sorrel	W18	Out-door_weeds allergen	Sorrel, Sheep/Red, Rumex acetosella
88	Nettle	W20	Out-door_weeds allergen	Nettle
89	Mugwort (Art v 1)	W231	Out-door_weeds allergen	Common mugwort (Artemisia vulgaris) pollen Art v 1
90	Walnut	T10	Out-door_Tree allergen	Walnut, Juglans regia
91	Maple leaf sycamore	T11	Out-door_Tree allergen	Sycamore
92	White ash	T15	Out-door_Tree allergen	Ash, White Fraxinus americana
93	Alder (Aln g 1)	T2	Out-door_Tree allergen	A primary pollen allergen of the European alder tree (Alnus glutinosa) Aln g 1.0101
94	Alder (Aln g 4)	T2	Out-door_Tree allergen	Aln g 4, a calcium-binding polcalcin (panallergen) protein from the pollen of European alder tree (Alnus glutinosa).
95	Horn beam tree (Car b 1)	T209	Out-door_Tree allergen	Car b 1, a major pollen allergen in the European hornbeam tree (Carpinus betulus).
96	Birch (Bet v 1)	T215	Out-door_Tree allergen	Major birch pollen allergen Bet v 1 (Betula verrucosa)
97	Birch (Bet v 2)	T216	Out-door_Tree allergen	Minor allergen in the pollen of the European white birch (Bet v 2 or Profilin-1)
98	Pecan/Hickory	T22	Out-door_Tree allergen	Pecan, Carya illinoensis
99	Birch (Bet v 4)	T220	Out-door_Tree allergen	Bet v 4, a minor allergen and calcium-binding protein (polcalcin) found in mature birch (Betula verrucosa) pollen
100	Birch (Bet v 6)	T225	Out-door_Tree allergen	A minor birch pollen allergen Bet v 6.
101	European ash (Fra e 1)	T25	Out-door_Tree allergen	A major pollen allergen in the European ash tree (Fraxinus excelsior)
102	Birch	T3	Out-door_Tree allergen	European white birch
103	Mountain cedar	T6	Out-door_Tree allergen	Cedar, Red_Juniperus virginiana
104	Oak	T7	Out-door_Tree allergen	Oak, Western White Quercus garryan
105	Mulberry pollen	T70	Out-door_Tree allergen	Mulberry
106	Olive pollen	T9	Out-door_Tree allergen	Olive
107	Positive control		internal control	
108	Negative control		internal control	