

## OmicsArray™ Allergen Antigen Array Antigen List - Catalog No. PA006

Antigen number	Antigen Name
1	a Bet v 1-homologous panallergen, belonging to the PR-10 protein family
2	a seed storage protein from <i>Arachis hypogaea</i> (peanut) found in the cotyledon.
3	a seed storage protein, 11 S globulin and trypsin inhibitor from peanut
4	Alfalfa_ <i>Medicago sativa</i>
5	allergen for plant polysensitization
6	<i>Alternaria alternata</i> / <i>Alternaria tenuis</i>
7	<i>Anisakis</i> / <i>Anisakis simplex</i>
8	Ant, Fire_ <i>Solenopsis invicta</i>
9	Ash, White_ <i>Fraxinus americana</i>
10	<i>Aspergillus flavus</i>
11	Bahia_ <i>Paspalum notatum</i>
12	Banana_ <i>Musa paradisiaca</i> var. <i>paradisiaca</i>
13	Barley Grain Dust
14	Barley Smut_ <i>Ustilago nuda</i>
15	Beef_ <i>Bos taurus</i>
16	Bermuda Grass Smut_ <i>Ustilago cynodontis</i>
17	Bermuda_ <i>Cynodon dactylon</i>
18	birch pollen allergen in apple, pear, orange, mango, lychee and other plants
19	Birch, Spring_ <i>Betula occidentalis</i> ( <i>fontinalis</i> )
20	calcium-binding pollen allergen <i>Aln g 4</i>
21	Cat Dander/Antigen_ <i>Felis catus</i> (domesticus)
22	Cat Epithelia_ <i>Felis catus</i> (domesticus)
23	Cat Hair_ <i>Felis catus</i> (domesticus)
24	Catfish_ <i>Ictalurus</i> spp.
25	Cedar, Red_ <i>Juniperus virginiana</i>
26	Cockroach, American_ <i>Periplaneta americana</i>
27	conglutin, belonging to the 2S albumin protein family
28	Corn Grain Dust
29	Crab_ <i>Callinectes</i> spp.

30	Dog Dander, Mixed-Breed_ <i>Canis familiaris</i>
31	Dog Epithelia, Mixed-Breed_ <i>Canis familiaris</i>
32	Dog Hair_ <i>Canis familiaris</i>
33	Glycine max seed storage protein
34	grass pollen allergen
35	group II/III of grass pollen allergen
36	group V allergen-starch granules of <i>Phleum pratense</i> pollens
37	hazelnut allergen from <i>Corylus avellana</i> .
38	Honey Bee_ <i>Apis mellifera</i>
39	House Dust
40	Iodine Bush_ <i>Allenrolfea occidentalis</i>
41	Lamb_ <i>Ovis aries</i>
42	Lobster_ <i>Homarus americanus</i>
43	Major apple ( <i>Malus domestica</i> ) allergen Mal d 1, a ribonuclease
44	major pollen allergen
45	Mattress Dust
46	Milk, Bovine_ <i>Bos taurus</i>
47	Mite, Food/Storage_ <i>Acarus siro</i>
48	Mite, Food/Storage_ <i>Lepidoglyphus destructor</i>
49	Mite, House Dust_ <i>Blomia tropicalis</i>
50	Mite, House Dust_ <i>Dermatophagoides farinae</i>
51	Mite, House Dust_ <i>Dermatophagoides pteronyssinus</i>
52	Mite, Storage_ <i>Chortoglyphus arcuatus</i>
53	Mixed Hornet_ <i>Dolichovespula</i> spp.
54	Oak, Western White/Garry_ <i>Quercus garryana</i>
55	Oat Grain Dust
56	Oat, Cultivated_ <i>Avena sativa</i>
57	peanut allergen
58	Peanut_ <i>Arachis hypogaea</i>
59	Plantain, English_ <i>Plantago lanceolata</i>
60	pollen allergen from <i>Corylus avellana</i>
61	pollen allergen of <i>Carpinus betulus</i>
62	pollen allergen of <i>Phleum pratense</i>
63	pollen allergen of <i>Phleum pratense</i>

64	pollen of <i>Alnus glutinosa</i>
65	profilin-1
66	Salmon_ <i>Salmo salar</i>
67	Shrimp_ Penacidae
68	soybean protein
69	The lipid transfer protein (LTP)
70	Tree allergen Ole e 1-like protein family
71	Upholstery Dust
72	Weed allergen of mugwort pollen
73	Wheat Grain Dust
74	wheat protein
75	Wheat, Whole_ <i>Triticum aestivum</i>
76	Ig control
77	Ig control
78	Ig control
79	Ig control
80	anti-Ig control
81	anti-Ig control
82	anti-Ig control
83	anti-Ig control