

**OmicsArray™ Ganglioside and Phospholipid Antigen Array Antigen List - Catalog No. PA015**

<b>Antigen number</b>	<b>Antigen name</b>	<b>Description</b>	<b>Antigen Type</b>
1	CL	Cardiolipin (CL)	phospholipid
2	GD1A	Disialoganglioside GD1A	ganlioside
3	GD1B	Disialoganglioside GD1B	ganlioside
4	GD2	Disialoganglioside GD2	ganlioside
5	GD3	Disialoganglioside GD3	ganlioside
6	GM1	Monosialoganglioside GM1	ganlioside
7	GM2	Monosialoganglioside GM2	ganlioside
8	GM3	Monosialoganglioside GM3	ganlioside
9	GM4	Monosialoganglioside GM4	ganlioside
10	GQ1B	Ganglioside GQ1B	ganlioside
11	GT1A	Ganglioside GT1A	ganlioside
12	GT1B	Ganglioside GT1B	ganlioside
13	PA	phosphatidic acid (PA)	phospholipid
14	PC	Phosphatidylcholine (PC)	phospholipid
15	PE	Phosphatidylethanolamine (PE)	phospholipid
16	PG	Phosphatidylglycerols (PG)	phospholipid
17	PI	Phosphatidylinositol (PI)	phospholipid
18	PS	Phosphatidylserine (PS)	phospholipid
19	S-1-P	Sphingosine-1-phosphate (S-1-P)	phospholipid
20	SM	Sphingomyelin	phospholipid
21	Sulfatide	Sulfatides	ganlioside
22	IgC	Ig control	control
23	algC	anti-Ig Control	control