

OmicsArray™ Bacterial Surface Antigen Array, 14-plex Antigen List - Catalog No. PA007

Antigen Number	Antigen ID	Bacterial Strain
1	NM	<i>Neisseria mucosa</i>
2	EY	<i>Eubacterium yurii</i>
3	AO	<i>Actinomyces oris</i>
4	RD	<i>Rothia dentocariosa</i>
5	NM2	<i>Neisseria macacae</i>
6	NS	<i>Neisseria sicca</i>
7	NE	<i>Neisseria elongata</i>
8	NF	<i>Neisseria flavescens</i>
9	AA	<i>Aggregatibacter aphrophilus</i>
10	SP	<i>Streptococcus parasanguinis</i>
11	LG	<i>Lactobacillus gasseri</i>
12	SS	<i>Streptococcus sanguinis</i>
13	PMSSI	<i>H. pylori</i> (PMSSI)
14	BC	<i>Blantia Coccoides</i>
15	Ig control_1	Ig control
16	Ig control_2	Ig control
17	Ig control_3	Ig control
18	Ig control_4	Ig control
19	anti-Ig control_1	anti-Ig control
20	anti-Ig control_2	anti-Ig control
21	anti-Ig control_3	anti-Ig control
22	anti-Ig control_4	anti-Ig control