

**OmicsArray™ ANA/ENA Array Panel - 32-plex antigen list -
Catlog# PA008**

| Antigen# | Antigen ID |
|-----------------|-----------------------------|
| 1 | dsDNA |
| 2 | ssDNA |
| 3 | Histone Core |
| 4 | Histone H1 |
| 5 | Histone H2B |
| 6 | Histone H2A |
| 7 | Histone H3 |
| 8 | Histone H4 |
| 9 | SSA-Ro52 |
| 10 | SSA-Ro60 |
| 11 | SSB-La |
| 12 | U1-sn-RNP A |
| 13 | U1-sn-RNP BB |
| 14 | U1-sn-RNP C |
| 15 | U1-sn-RNP 70 |
| 16 | RNP/Sm |
| 17 | SmD |
| 18 | SmD1 |
| 19 | SmD2 |
| 20 | SmD3 |
| 21 | SmD4 |
| 22 | Ribo Phosphaprotein P0 |
| 23 | Ribo Phosphaprotein P1 |
| 24 | Ribo Phosphaprotein P2 |
| 25 | Nucleosome |
| 26 | Scl-70/Topoisomerase I |
| 27 | Jo-1 |
| 28 | CENP-A |
| 29 | CENP-B |
| 30 | Ku 70/80 |
| 31 | Th/To (RPP25) |
| 32 | RNA polymerase III (POLR3D) |
| 33 | Ig control 1 |
| 34 | Ig control 2 |
| 35 | Ig control 3 |
| 36 | Ig control 4 |
| 37 | anti-Ig control 1 |
| 38 | anti-Ig control 2 |
| 39 | anti-Ig control 3 |
| 40 | anti-Ig control 4 |