

| OmicsArray™ Rheumatoid Arthritis (RA) Antigen Array Antigen List |                                                      |                                                                                       | Catalog No. PA022                                                                        |                        |
|------------------------------------------------------------------|------------------------------------------------------|---------------------------------------------------------------------------------------|------------------------------------------------------------------------------------------|------------------------|
| Antigen Number                                                   | Antigen ID                                           | Antigen Name                                                                          | Gene symbol                                                                              | UniProt ID             |
| 1                                                                | CCP2                                                 | cyclic citrullinated peptide 2                                                        |                                                                                          |                        |
| 2                                                                | PAD2                                                 | Peptidylarginine deiminase II                                                         | PADI2                                                                                    | <a href="#">Q9Y2I8</a> |
| 3                                                                | PAD4                                                 | Peptidylarginine deiminase IV                                                         | PADI4                                                                                    | <a href="#">Q9UM07</a> |
| 4                                                                | Collagen type II                                     | Collagen Alpha-1 type II                                                              | COL2A1                                                                                   | <a href="#">P02458</a> |
| 5                                                                | H2A                                                  | Histone H2A                                                                           | H2AC7                                                                                    | <a href="#">P20671</a> |
| 6                                                                | H2B                                                  | Histone H2B                                                                           | H2BC3                                                                                    | <a href="#">P33778</a> |
| 7                                                                | H3                                                   | Histone H3                                                                            | H3Y1                                                                                     | <a href="#">P0DPK2</a> |
| 8                                                                | H4                                                   | Histone H4                                                                            | H4C1; H4C2; H4C3; H4C4; H4C5; H4C6; H4C8; H4C9; H4C11; H4C12; H4C13; H4C14; H4C15; H4C16 | <a href="#">P62805</a> |
| 9                                                                | Ro52 (SSA-52)                                        | Ro/SSA 52KD, TRIM21                                                                   | TRIM21                                                                                   | <a href="#">P19474</a> |
| 10                                                               | Ro60 (SSA-60)                                        | Ro/SSA 60KD, TROVE2                                                                   | RO60                                                                                     | <a href="#">P10155</a> |
| 11                                                               | Centromere B                                         | Centromere protein B (CENP-B)                                                         | CENPB                                                                                    | <a href="#">P07199</a> |
| 12                                                               | IGFBP2                                               | Insulin-like Growth Factor Binding Protein 2                                          | IGFBP2                                                                                   | <a href="#">P18065</a> |
| 13                                                               | Fibrinogen                                           | Fibrinogen                                                                            | FGB                                                                                      | <a href="#">P02675</a> |
| 14                                                               | Cit-Fibrinogen                                       | Human Fibrinogen (PAD2 Citrullinated)                                                 | FGB                                                                                      | <a href="#">P02676</a> |
| 15                                                               | α-Enolase                                            | Alpha-enolase (ENO1)                                                                  | ENO1                                                                                     | <a href="#">P06733</a> |
| 16                                                               | Cit-α-Enolase (CEP1)                                 | Citrullinated α-Enolase (human)                                                       | ENO1                                                                                     | <a href="#">P06733</a> |
| 17                                                               | Cit-H2A                                              | Citrullinated Histone H2A Type 1 (human)                                              | H2AC7                                                                                    | <a href="#">P20671</a> |
| 18                                                               | Cit-H2B                                              | Citrullinated Histone H2B                                                             | H2BC3                                                                                    | <a href="#">P33778</a> |
| 19                                                               | Cit-H3                                               | Citrullinated Histone H3                                                              | H3Y1                                                                                     | <a href="#">P0DPK2</a> |
| 20                                                               | Cit-H4                                               | Citrullinated Histone H4 (human)                                                      | H4C1; H4C2; H4C3; H4C4; H4C5; H4C6; H4C8; H4C9; H4C11; H4C12; H4C13; H4C14; H4C15; H4C16 | <a href="#">P62805</a> |
| 21                                                               | Vimentin                                             | Vimentin                                                                              | VIM                                                                                      | <a href="#">P08670</a> |
| 22                                                               | Cit-Vimentin (Sa)                                    | Citrullinated Vimentin                                                                | VIM                                                                                      | <a href="#">P08670</a> |
| 23                                                               | cit-LL37                                             | Citrullinated LL37                                                                    |                                                                                          |                        |
| 24                                                               | Cit-FLG                                              | citrullinated Filaggrin                                                               | FLG                                                                                      | <a href="#">P20930</a> |
| 25                                                               | GNPTG                                                | N-acetylglucosamine -1-phosphotransferase subunit gamma                               | GNPTG                                                                                    | <a href="#">Q9UJJ9</a> |
| 26                                                               | RF                                                   | Rheumatoid factor (RF)                                                                |                                                                                          |                        |
| 27                                                               | BRAF                                                 | Serine/threonine-protein kinase B-raf, v-Raf murine sarcoma viral oncogene homolog B1 | BRAF                                                                                     | <a href="#">P15056</a> |
| 28                                                               | Acetylated ornithine (AcOrn)                         | Nα-Acetyl-L-ornithine                                                                 |                                                                                          |                        |
| 29                                                               | N-acetylglucosamine -1-phosphatase transferase (OGT) | Human O-GlcNAcTransferase/OGT Protein                                                 | OGT                                                                                      | <a href="#">O15294</a> |
| 30                                                               | Malondialdehyde-acetaldehyde (MAA) modified protein  | MAA modified albumin (MAA-Alb)                                                        |                                                                                          |                        |
| 31                                                               | Carbamylated Albumin                                 | Carbamylated Albumin (Car-Alb)                                                        |                                                                                          |                        |

|    |                         |                                   |  |  |
|----|-------------------------|-----------------------------------|--|--|
| 32 | Carbamylated Fibrinogen | Carbamylated fibrinogen (Car-Fib) |  |  |
| 33 | Ig control 1            | Internal control                  |  |  |
| 34 | Ig control 2            | Internal control                  |  |  |
| 35 | Ig control 3            | Internal control                  |  |  |
| 36 | Ig control 4            | Internal control                  |  |  |
| 37 | anti-Ig control 1       | positive control                  |  |  |
| 38 | anti-Ig control 2       | positive control                  |  |  |
| 39 | anti-Ig control 3       | positive control                  |  |  |
| 40 | anti-Ig control 4       | positive control                  |  |  |