Rheumatoid Arthritis: Epigenetic Drug Targets

Adam P. Cribbs^{1,2}, Marc Feldmann² and Udo Oppermann^{2,3}

¹Computational Genomics and Training Centre, MRC Weatherall Institute of Molecular Medicine, University of Oxford, John Radcliffe Hospital, Headington, Oxford, OX3 9DS, UK ²Botnar Research Centre, University of Oxford, Windmill Road, Oxford, OX3 7LD, UK ³Structural Genomics Consortium, University of Oxford, Old Road Campus, Headington, Oxford, OX3 7DQ, UK

Autoimmune diseases, such as rheumatoid arthritis (RA) arise as a result of a breakdown in immune tolerance, for reasons that are as yet unknown. RA affects up to 1% of the population worldwide and has a lifetime risk of 4% for women and 2% for men. While a number of successful immunotherapies, such as anti-TNF, CTLA-4lg and anti-CD20, have revolutionized the treatment of RA in recent years, many patients still fail to respond sufficiently to treatment. Moreover, these costly therapies are not curative, so treatment must be continued for many years to manage disease. Although the etiology and pathogenesis of RA remains elusive, studies have revealed a genetic predisposition, with variants of several genes, including human leukocyte antigen DRB1 (HLA-DRB1), IRF5, PTPN22, CD28 and CTLA-4, having strong association with the condition. Despite this extensive genetic predisposition, the concordance of RA in monozygotic twins is low (12–20%), suggesting that environmental and epigenetic factors are important in disease development⁽¹⁾. The environmental factor most consistently associated with RA is cigarette smoke, with alcohol consumption seeming to have a mild protective effect.

RA Pathology

RA is a chronic destructive inflammatory disease that is characterized by severe swelling, pain and stiffness of the joints. In healthy individuals the synovial membrane, which lines the joint capsule, acts to lubricate the joint by producing synovial fluid. The synovial membrane contains two main immune cell types, macrophages and synovial fibroblasts (synoviocytes) that lie within the connective tissue. In RA, however, there is an influx of T cells, NK cells and B cells, in addition to other chronically inflamed cell types; the synoviocytes become activated and invade the surrounding cartilage, resulting in cartilage degradation and bone erosion.

RA Inflammation

Antigen-activated T cells produce pro-inflammatory cytokines, such as IL-17, that stimulate monocytes, macrophages and synovial fibroblasts to produce other pro-inflammatory cytokines, IL-1, IL-6 and TNF- α . This leads to recruitment of other immune cells, such as NK cells and B cells, which in turn secrete IFN-y and immunoglobulins such as rheumatoid factor, respectively. Overall, the combination of these proinflammatory cytokines contribute to the articular inflammation. The activation of macrophages, lymphocytes and fibroblasts can also induce angiogenesis, which may explain the increased vascularity found in the synovium of RA patients. Moreover, the endothelial cells in these newly formed vessels express high levels of adhesion markers leading to increased inflammator cell recruitment and infiltration. Activated CD4⁺ T cells express osteoprotegerin ligands that stimulate osteoclastogenesis similarly NK cells from the RA synovium express increased RANKL, which also stimulates osteoclastogenesis, leading to bone erosion.



Figure 1 Typical X-ray of hand with advanced rheumatoid arthritis. Bone erosion, cartilage degradation and bone displacement are classic features of this disease (from Bernd Brägelmann via Wikimedia Commons)



TOCRIS a **biotechne** brand

Neurochemicals | Signal Transduction Agents | Peptides | Biochemicals

www.tocris.com

Products available from Tocris

Aurora Kinases

Anacardic acid, PE 03814735. TC-A 2317, VX 680, ZM 447439 Bromodomains

(-)-JQ1, (+)-JQ1, BI 9564, I-BET 151, -CBP 112, Ischemin, PFI 1, SGC-CBP30

DNA Methyltransferases 5-Azacytidine, Decitabine, RG 108, SGI 1027, Zebularine

Histone Acetyltransferases

C 646, Garcinol, L002, NU 9056 **Histone Deacetylases**

CI 994, FK 228, MI 192, SAHA Scriptaid, Trichostatin A, UF 010 Valproic acid sodium salt, GSK J1 GSK J2, GSK J4, GSK J5, IOX 1 **JIB 04** JAK

AG 490, AZD 1480, CP 690550 Lysine Methyltransferases

3-Deazaneplanocin A, UNC 1999, UNC 2400

Sir2-like Family Deacetylases EX 527, Resveratrol, Sirtinol

Ubiquitin E3 Ligases GS 143, SMER 3, SZL P1-41

Calcium-Sensitive Proteases MG 132

Chemokine Receptors

AMD 3100, AMD 3465, SB 225002 Cvtokines

(D)-(+)-Neopterin, 4-IPP. CRID3. DMXAA, ISO 1, Pirfenidone DMARDs

Methotrexate

Indoleamine 2,3-dioxygenase 1-Methyl-D-tryptophan

Matrix Metalloprotease

Batimastat, GI 254023X, GM 6001, Marimastat

Nitric Oxide Signaling

1400W, BYK 191023, L-NAME, SNAP

p38

AMG 548, SB 202190, SB 203580, SB 239063, VX 745

Ac, acetylation DMARD, disease-modifying

antirheumatic drug IFN, interferon

IL, interleukin

NK, natural killer **PBMC**, peripheral blood mononuclear

RANKL, receptor activator of nuclear factor κ-B ligand

Th, T helper

TNF, tumor necrosis factor **T_{regs}, regulatory T cells**

- L. Svendsen *et al.* (2013) *PloS one* 8(2):e57304. . Wegner *et al.* (2010) Arthritis Rheum.
- Yang et al. (2007) Am.J.Physiol.Lung Cell.Mol.Physiol. 292(2):L567.
 Cribbs et al. (2014) Arthritis Rheum.
- 66(9):2344.
- Cribbs et al. (2015) Ther.Adv.
- *Musculoskelet.Dis.* 7(5):206. 5. Klein *et al.* (2016) *Ann.Rheum.Dis.*
- Oppermann (2013) Arthritis Res.Ther. 15(2):209.

For copies of this poster, please visit tocris.com © 2016 Tocris Cookson, Ltd. Tocris is a Bio-Techne brand