



# High-Impact Publications in Stem Cell Research, 2022–2023

## ZE5 Cell Analyzer Publications

### **Aikins ME et al. (2022).**

Cancer stem cell antigen nanodisc cocktail elicits anti-tumor immune responses in melanoma.  
J Control Release 351, 872–882.

### **Augimeri G et al. (2023).**

A hybrid breast cancer/mesenchymal stem cell population enhances chemoresistance and metastasis.  
JCI Insight 8, e164216.

### **Bealer E et al. (2023).**

Extrahepatic transplantation of 3D cultured stem cell–derived islet organoids on microporous scaffolds.  
Biomater Sci 11, 3645–3655.

### **Chen YH et al. (2023).**

Rewilding of laboratory mice enhances granulopoiesis and immunity through intestinal fungal colonization.  
Sci Immunol 8, eadd6910.

### **Crees ZD et al. (2023).**

Motixafortide and G-CSF to mobilize hematopoietic stem cells for autologous transplantation in multiple myeloma:  
A randomized phase 3 trial.  
Nat Med 29, 869–879.

### **Harper TC et al. (2023).**

GATA1 deletion in human pluripotent stem cells increases differentiation yield and maturity of neutrophils.  
iScience 26, 107804.

### **Kang YA et al. (2023).**

Secretory MPP3 reinforce myeloid differentiation trajectory and amplify myeloid cell production.  
J Exp Med 220, e20230088.

### **Larouche JA et al. (2023).**

Spatiotemporal mapping of immune and stem cell dysregulation after volumetric muscle loss.  
JCI Insight 8, e162835.

**Leclerc K et al. (2023).**

Hox genes are crucial regulators of periosteal stem cell identity.  
Development 150, dev201391.

**Lee JS et al. (2022).**

The insulin and IGF signaling pathway sustains breast cancer stem cells by IRS2/PI3K-mediated regulation of MYC.  
Cell Rep 41, 111759.

**Li F et al. (2022).**

Fas ligand-modified scaffolds protect stem cell derived  $\beta$ -cells by modulating immune cell numbers and polarization.  
ACS Appl Mater Interfaces 15, 50549–50559.

**Lopez de Lapuente Portilla A et al. (2022).**

Genome-wide association study on 13,167 individuals identifies regulators of blood CD34+cell levels.  
Blood 139, 1659–1669.

**Otsuka T et al. (2022).**

Overexpression of NDST1 attenuates fibrotic response in murine adipose-derived stem cells.  
Stem Cells Dev 31, 787–798.

**Tolstova T et al. (2023).**

The effect of TLR3 priming conditions on MSC immunosuppressive properties.  
Stem Cell Res Ther 14, 344.

Visit [bio-rad.com/ZE5](https://www.bio-rad.com/ZE5) for more information.

BIO-RAD is a trademark of Bio-Rad Laboratories, Inc. All trademarks used herein are the property of their respective owner.  
© 2024 Bio-Rad Laboratories, Inc.



**Bio-Rad  
Laboratories, Inc.**

Life Science  
Group

**Website** [bio-rad.com](https://www.bio-rad.com) **USA** 1 800 424 6723 **Australia** 61 2 9914 2800 **Austria** 00 800 00 24 67 23 **Belgium** 00 800 00 24 67 23 **Brazil** 4003 0399  
**Canada** 1 905 364 3435 **China** 86 21 6169 8500 **Czech Republic** 00 800 00 24 67 23 **Denmark** 00 800 00 24 67 23 **Finland** 00 800 00 24 67 23  
**France** 00 800 00 24 67 23 **Germany** 00 800 00 24 67 23 **Hong Kong** 852 2789 3300 **Hungary** 00 800 00 24 67 23 **India** 91 124 4029300 **Israel** 0 3 9636050  
**Italy** 00 800 00 24 67 23 **Japan** 81 3 6361 7000 **Korea** 82 080 007 7373 **Luxembourg** 00 800 00 24 67 23 **Mexico** 52 555 488 7670  
**The Netherlands** 00 800 00 24 67 23 **New Zealand** 64 9 415 2280 **Norway** 00 800 00 24 67 23 **Poland** 00 800 00 24 67 23 **Portugal** 00 800 00 24 67 23  
**Russian Federation** 00 800 00 24 67 23 **Singapore** 65 6415 3188 **South Africa** 00 800 00 24 67 23 **Spain** 00 800 00 24 67 23 **Sweden** 00 800 00 24 67 23  
**Switzerland** 00 800 00 24 67 23 **Taiwan** 886 2 2578 7189 **Thailand** 66 2 651 8311 **United Arab Emirates** 36 1 459 6150 **United Kingdom** 00 800 00 24 67 23

