

A Novel CAR-T Cell Therapy Approach Using Fluorescence-Activated Cell Sorting and Stem Cell Transplantation

By Mariko Alexander

Successful chimeric antigen receptor- (CAR-) T cell therapy requires a target antigen that is unique to cancer cells. But what happens when there are no unique antigens? Researchers at Columbia University Medical Center addressed this problem by replacing healthy non-target cells with genetically modified versions lacking the CAR-T cell target (Borot et al. 2019). Their results, published in PNAS, may provide a new avenue for treatment of some types of cancer.

CAR-T Cell Therapy and the Unique Antigen Problem

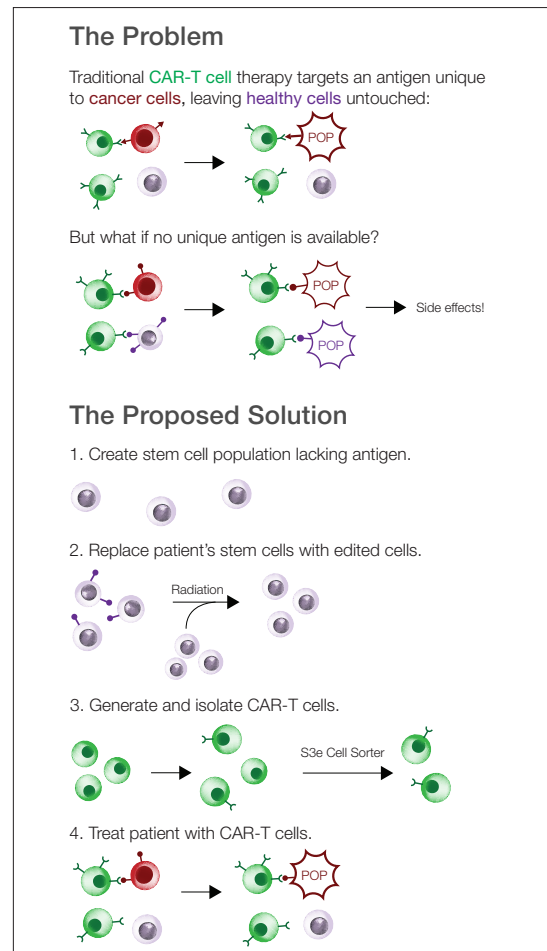
CAR-T cell therapy, an emerging cancer treatment strategy, uses T cells modified to express a CAR — a receptor engineered to recognize an antigen found on cancer cells. CAR-T cell therapy has proven highly successful for some treatment-resistant leukemias and initial studies extending the strategy to other types of cancer have been promising.

However, one of the major drawbacks for this approach is that CAR-T cells will target any cell displaying the target antigen, whether healthy or cancerous. Damage to healthy cells can result in severe side effects, such as immune cell depletion or organ damage, and even death.

Unfortunately, there are many types of cancer for which no unique antigen is known, leaving clinicians and researchers to decide whether the risk of side effects outweighs the potential benefits. This is the case for acute myeloid leukemia (AML), a type of bone marrow cancer for which new treatments are desperately needed. Immunotherapy targeting CD33, an antigen present in more than 80% of cases, has been successful for treatment of some patients with AML. However, the CD33 antigen is also found in certain immune system progenitor cells, causing severe complications in some cases.

A New Approach for CAR-T Cell Therapy

To get around the unique antigen problem, researchers Florence Borot and colleagues tested a creative solution. Rather than searching for a new antigen, they used gene editing to generate a population of healthy cells immune to the effects of CAR-T cell therapy targeting CD33.



To do this, researchers started with CD34-expressing hematopoietic stem cells isolated from bone marrow or cord blood. Using CRISPR-Cas9, expression of CD33 was ablated. Transplantation of the edited cells into mice effectively replaced the native population of bone marrow cells and their progenitors with the knockout cells, permitting safe treatment with CAR-T cells targeting CD33.

Results in vitro and in mouse models were promising, showing that the CAR-T cells successfully targeted cancer cells without affecting the edited healthy cells. These data provide hope that CAR-T cell therapy may yet be successful not only for AML, but also for other cancers faced with a lack of unique antigens.



Isolating T Cells with the S3e Cell Sorter

Researchers are increasingly incorporating a cell sorting step when generating modified cells. In this study, Borot and colleagues used a Bio-Rad S3e Cell Sorter to isolate CAR-expressing T cells after

transfection. The S3e Cell Sorter is particularly well suited for experiments like these due to its ease of use and its ability to produce highly pure cell populations while preserving viability.

The S3e Cell Sorter makes cell sorting approachable for novices and experts alike. Because it has a shallow learning curve, unlike many other cell sorters, it can easily be incorporated into laboratories that do not regularly perform flow cytometry. The simple setup also reduces the risk of mistakes introduced by unnecessarily complicated equipment, saving researchers' time and materials.

In this study, T cells isolated with the S3e Cell Sorter were injected into mice. For experiments like this one, the purity of the cells after sorting is critical. The S3e Cell Sorter produces highly pure cell populations, ensuring that contaminating cell types do not confound results or cause side effects in treated organisms. Poor purity may also lead to results that are not reproducible, wasting time and resources on downstream experiments and clinical trials.

The cell sorter must also be gentle enough to work without compromising viability. Samples that do not survive the duration of the experiment are useless regardless of purity. Stressed cells may also behave differently even if they are still viable, affecting reproducibility and confounding results. The S3e Cell Sorter was designed to handle cells gently, preserving their health and allowing researchers to experiment without worry.

Visit bio-rad.com/S3e for more information about the S3e Cell Sorter.

References

- Borot F et al. (2019). Gene-edited stem cells enable CD33-directed immune therapy for myeloid malignancies. *Proc Natl Acad Sci USA* 116, 11,978–11,987.
- Borot F et al. (2019). Correction for Borot et al., Gene-edited stem cells enable CD33-directed immune therapy for myeloid malignancies. *Proc Natl Acad Sci USA* 116, 14,780–14,781.

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

Bioradiations.com is an online magazine that offers scientific professionals around the globe timely tips, techniques, and topics related to Bio-Rad products and services. The site offers new product information, technical and application notes describing experiments performed using Bio-Rad products, interviews with renowned researchers in the field, and product demonstrations in the form of system tours, videos, podcasts etc. Visit us at bioradiations.com.



**Bio-Rad
Laboratories, Inc.**

Life Science
Group

Website 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 2 3473 4460 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

