

Evaluating CAR-T Cell Dynamics in a 3D Tumor Microenvironment Using Confocal Imaging

Evaluation of the immune effect of CAR-T (Chimeric Antigen Receptor T) cell therapy is usually performed using a model organism, which is costly and time consuming. This application note introduces an example of building a simple 3D immune cell-mediated killing assay using AIM Biotech's 3D cell culture chips, and measuring the immune effects of T cells by *in vitro* imaging. The 3D assay makes it easy to probe different conditions *in vitro* such as the cancer microenvironment and T cell regulation, and it can be customized in various ways according to the purpose of the research. This assay reproduces the more spatiotemporal dynamics of cells *in vitro* and enables the analysis of immune cell-mediated killing under more physiological conditions as compared to 2D models.

Experiment Overview

We evaluated the killing efficiency of antitumor effector T cells (CAR-T cells) by constructing a 3D microenvironment that mimics a tumor using AIM Biotech's idenTx 9 3D cell culture chip and observing and quantitatively analyzing images using AX Confocal Microscope from Nikon.

The AIM Biotech idenTx 9 system used in this experiment consists of 3 fluid channels (Fig. 1A). Firstly, GFP-labeled cancer cells were mixed with Collagen Type I derived from rat tails, and seeded in the central hydrogel channel. The 3 channels can be compartmentalized by having polymerized hydrogel in the center (Fig. 1B, left).

Next, CAR-T cells were added to the media channel adjacent to the hydrogel channel (Fig. 1B, center). Then, the infiltration of CAR-T cells into the cancer cell-populated hydrogel channel was observed in time series (Fig. 1B, right). In this experiment, 3 types of samples, "Cancer cells only," "Cancer cells + CD133 specific CAR-T cells" and "Cancer cells + Mock transfected T cells," were each prepared and fixed after 24 hours and 120 hours. The infiltration of CAR-T cells was evaluated and the rate of apoptosis of cancer cells was quantified based on the 3D fluorescent images (Fig. 2 and Fig. 3).

Fig.1: Creating a 3D assay model using the idneTx 9 system

(A) Top view and cross section of a chip in the idneTx 9 system

a: Site, b: Media channel, c: Gel channel, d: Media inlet, e: Gel inlet, f: Port, g: Trough

(B) Cell seeding procedures using the idneTx 9 system

Cancer cells: GFP labeled Hep3B cells (CD133 positive hepatocellular carcinoma),

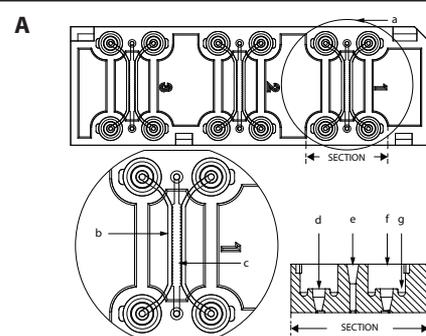
Effector T cells: mCherry labeled CD133 specific CAR-T cells

Control effector cells: mCherry labeled mock transfected T cells,

Primary antibody: Cleaved Caspase-3 (Asp175) antibody; Cell Signaling Technology; cat #9661,

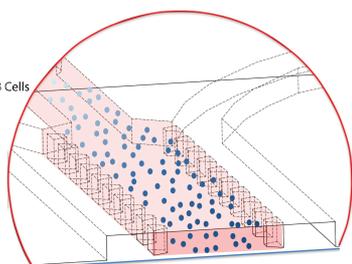
Secondary antibody: Goat anti-Rabbit IgG (H+L) Highly Cross-Adsorbed Secondary Antibody, Alexa 633;

Thermo Fisher; cat # A-21071



B

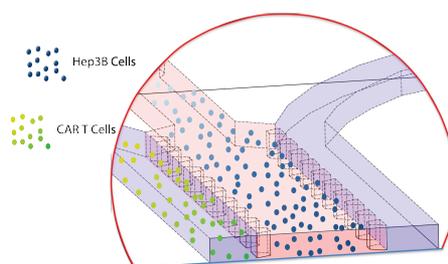
Hep3B Cells



Collagen gel that was mixed with Hep3B cells was loaded into idenTx 9

Hep3B Cells

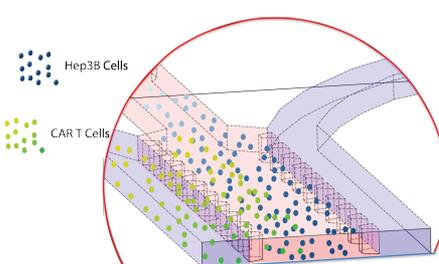
CAR T Cells



CAR-T cells were seeded in the media channel

Hep3B Cells

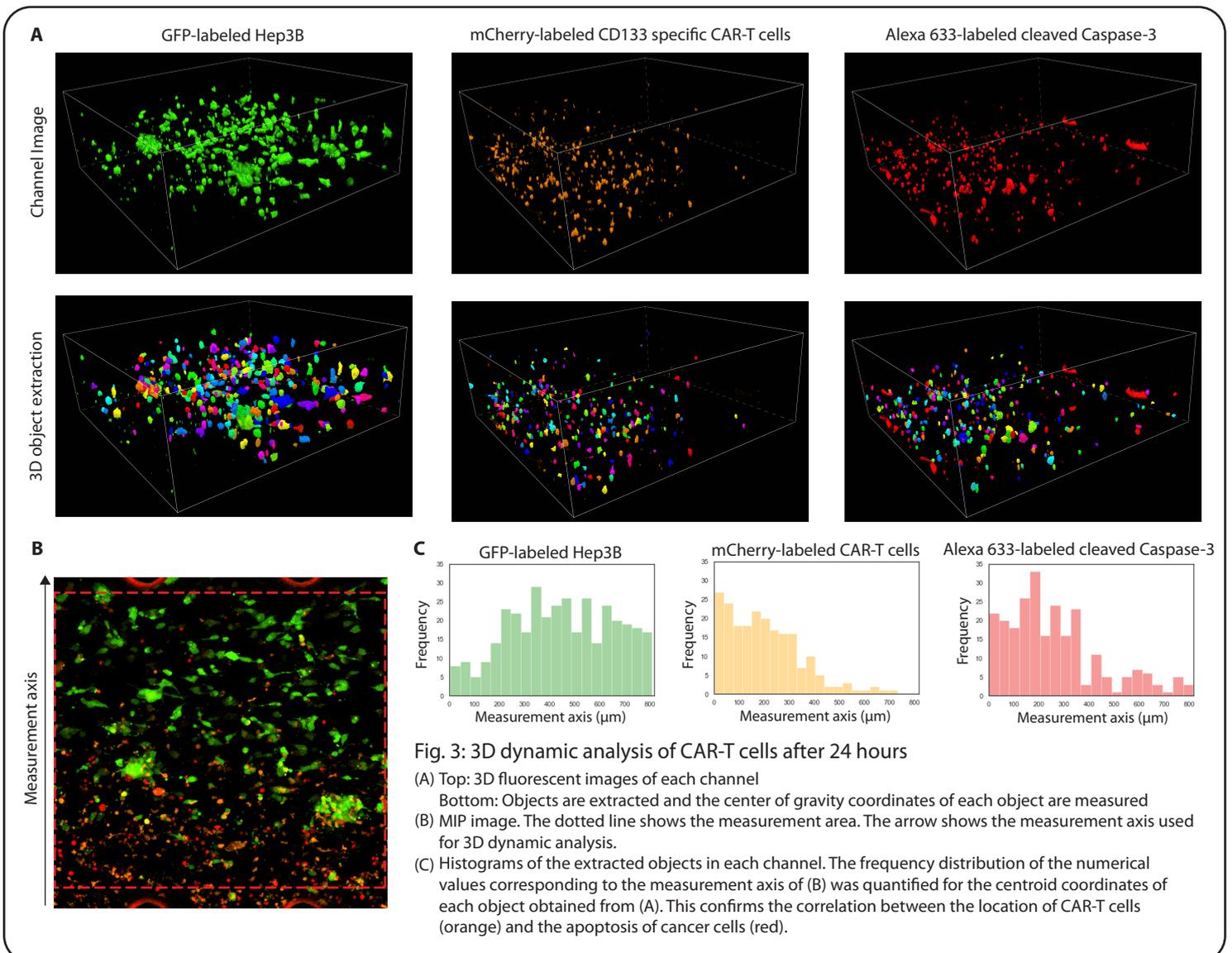
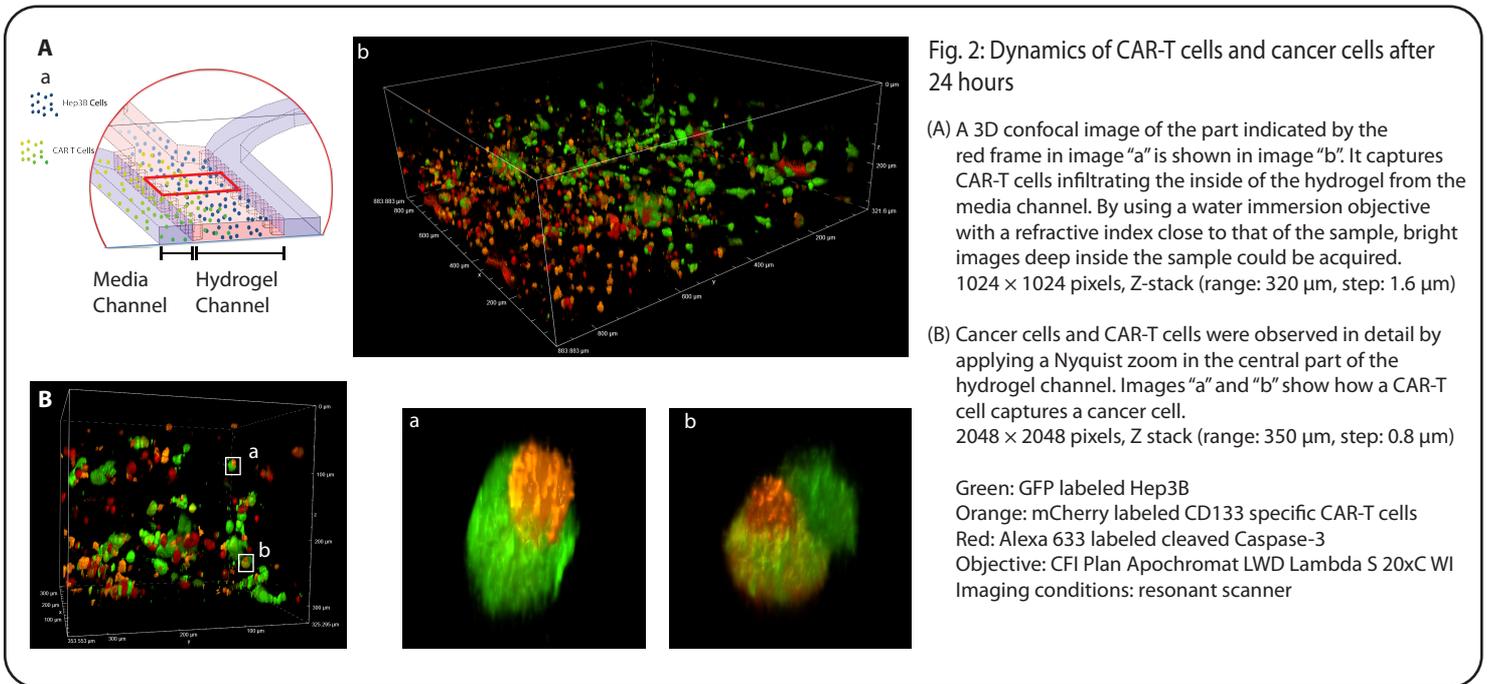
CAR T Cells



Infiltration of CAR-T cells and their killing ability against target cells were monitored

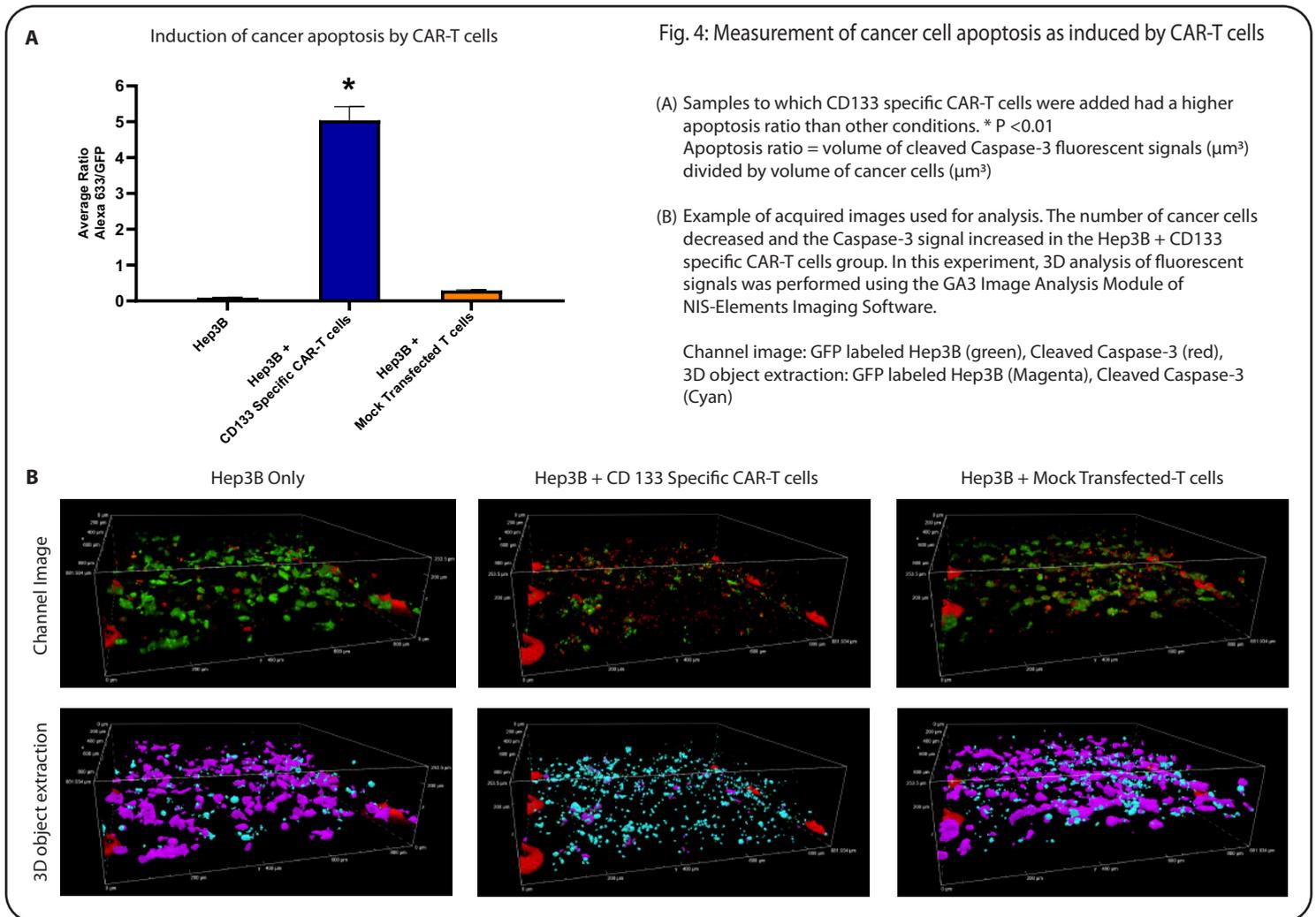
Infiltration of CAR-T cells induced the apoptosis of cancer cells

CD133 specific CAR-T cells, the effector cells, infiltrated the tumor microenvironment spontaneously within 24 hours and came into contact with the cancer cells (Fig. 2). Also, the dynamics of the fluorescent signals of cancer cells, CAR-T cells, and cleaved Caspase-3 were analyzed, and a correlation was found in the position of their centers of gravity (Fig. 3). In other words, this suggests that CAR-T cells moved towards the cancer cells, stayed near them and induced their apoptosis.



CD133 specific CAR-T cells have a high killing ability against cancer cells

Next, a comparison of the induction of cancer cell apoptosis by CAR-T cells was performed using 3 types of samples: "Cancer cells only", "Cancer cells + CD133 specific CAR-T cells", and "Cancer cells + Mock transfected T cells". All were fixed after 120 hours. The results revealed that the ratio (volume ratio) of fluorescent signals of cleaved Caspase-3 per cancer cell was significantly higher in the group with CD133 specific CAR-T cells than in the group with mock transfected T cells (Fig. 4). The "Cancer cells only" control group had the lowest apoptosis ratio, as expected.



Product Information

idenTx 9 Plate



The idenTx 9 fully integrates the capacity of three individual idenTx 3 chips into a standard SBS plate format, enabling 9 simultaneous experiments on a single plate, and rapid scaling up for higher throughput drug screening models. The idenTx 9 retains the same ease-of-use as the idenTx 3 and is a seamless way to progress through feasibility while maintaining the common handling features and laboratory automation compatibility that comes with the standard lab plate format.

AX Confocal Microscope (Nikon)

Achieves high resolution images of 8192 x 8192 pixels, which are four times that of conventional models. With a large diagonal field of view of 25 mm, wide areas of samples can be acquired at once, reducing phototoxicity. An automatic shading correction function enables acquisition of images with uniform brightness.

