

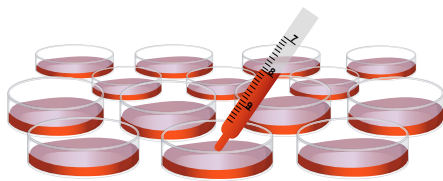
# OncoFlow-Profiler

Elevate your ADC antigen discovery with our rapid, high-throughput flow cytometry screening platform

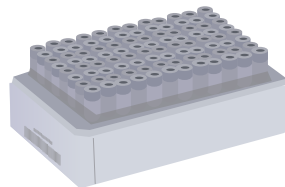
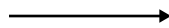


Introducing OncoFlow-Profiler - our state-of-the-art high-throughput screening solution crafted to fast-track the detection of your target antigen expression in a well-established panel of 160 cancer cell lines. Within just 2 weeks, you can obtain high-quality flow cytometry data, facilitating early informed decisions in the ADC drug discovery process.

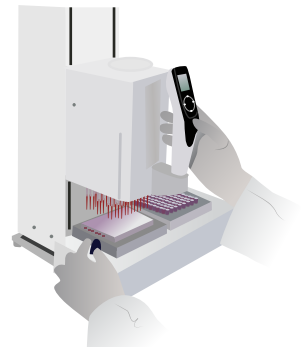
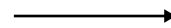
## Experimental Procedure



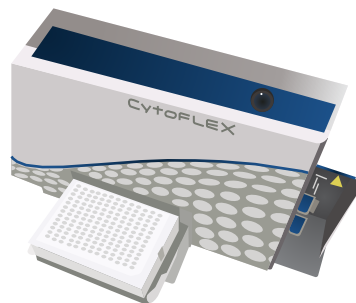
Culture & expansion of 160 cell lines



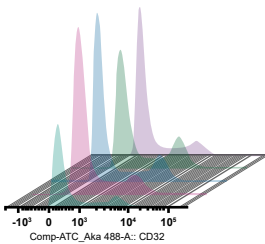
Viability stain, fixation, cryopreservation



Thawing, washing, staining



Data acquisition



Data analysis

## Benefits

- Ⓜ High-throughput screening platform
- Ⓜ Gain access to 160 well-established cancer cell lines
- Ⓜ Quick turnaround; results in 2 weeks
- Ⓜ Capable of testing up to 4 ADCs simultaneously within a single experimental batch
- Ⓜ Detects carbohydrate-based antigens as well
- Ⓜ Streamlined experimental staining procedure with analysis of 4 markers in 160 cell lines within a single day



Take the guesswork out of target antigen expression. Reach out to explore the power of OncoFlow-Profiler now.