

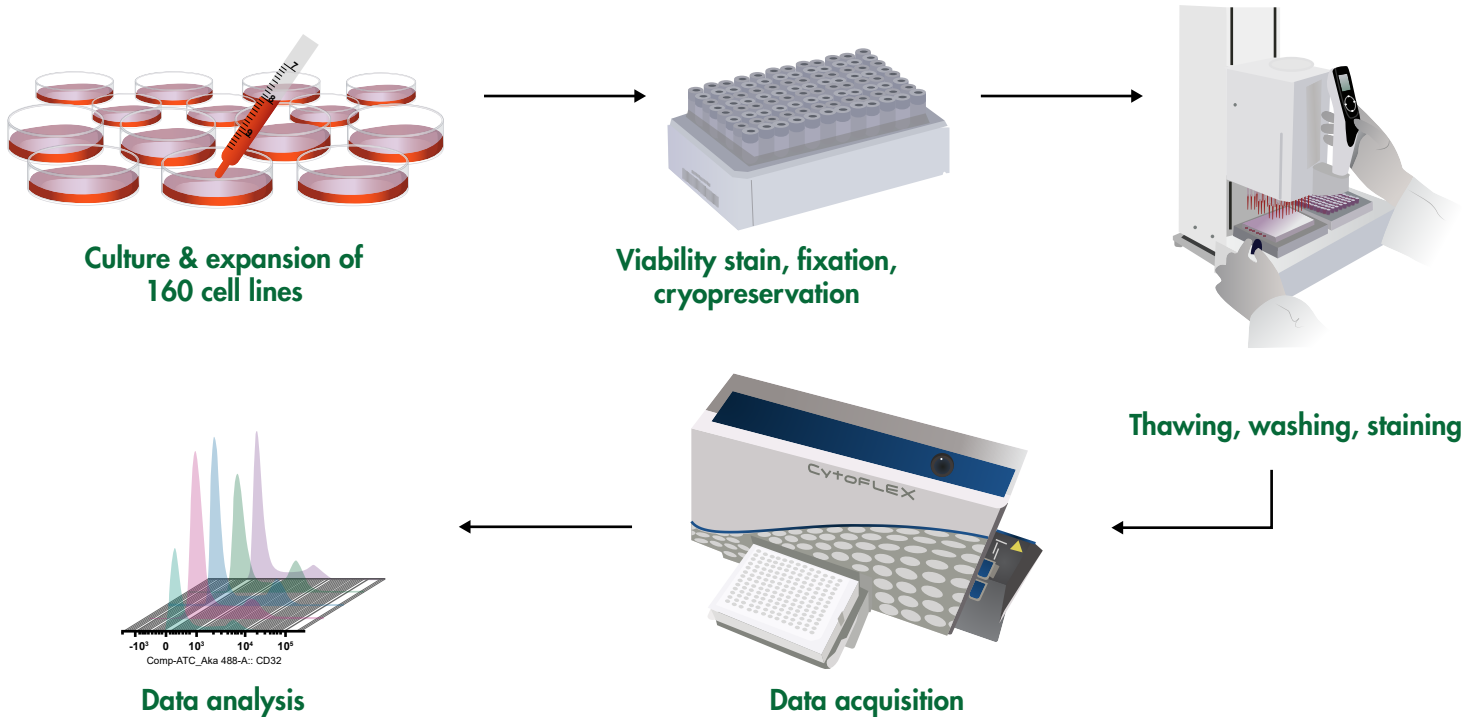
OncoFlow-Profiler

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



Benefits

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