



Antibody internalization assays for cancer drug discovery

N. Bevan, T. Dale & D. Trezise (sponsor Daniel Martinez Molina) Essen BioScience, Welwyn Garden City, AL7 3AX, UK

Summary & Impact

- Monoclonal antibodies (mAb) and antibody-drug conjugates are widely used biological therapeutics.
- A key property is the extent and rate of internalization, governing efficacy, safety and PD profile, thus
- An antibody-binding fragment coupled to a pH-sensitive dye (FabFluor-pH) was used to label and test mAbs using a single-step, no-wash protocol.
- As expected, an increase in fluorescence signal was

Continuous Live-Cell Analysis: Methodology







- quantifying and comparing Ab internalization is a critical step in the biopharmaceutical process.
- Here, we describe an automated, novel and enabling cell-based Ab internalization assay that is turnkey, medium throughput and geared toward industrial biologics discovery.
- Internalization measurements are made over time on 96well microplates using live-cell analysis (IncuCyte[®])
- observed as the mAb complex was internalized into the acidic lysosome.
- This approach has been validated for use in pharmacological and temporal characterisation of mAb.
- Proof of concept as part of a screening cascade to compare Ab characteristics and specificity of signal across cell lines and epitopes.





- BT-474 Her2-positive breast cancer cells (10K/well) were treated with IncuCyte FabFluor-pH labeled-Herceptin or IgG, isotype control.
- Red fluorescence and phase images were taken (10x) every 30 min over 48 h.
- Quantification of the red fluorescence area, normalized for phase area (confluence) demonstrates a rapid, concentration-dependent increase over time with labeled Hercpetin but no signal with IgG.



- Raji cells (30K/well) were treated with either IncuCyte FabFluor-pH ____ labeled-Rituxan or IgG, isotype control.
- Red fluorescence and phase images were taken (20x) every 30 min over 36 h.
- Quantification of the red fluorescence, normalized for phase area (confluence) demonstrates a rapid, concentration dependent increase in signal over time with labeled Rituxan but no signal with IgG.

Drug discovery screening application





- Multiple anti-CD71 Abs were labeled and internalization properties compared (triplicate wells, CRC) in HT1080 fibrosarcoma cells (8K/well).

	Seed cells (50 µL/well, 5,000-	IncuCyte [®] FabFluor Reagent	labeled Antibody Addition
	30,000 cells/well), into 96-well plate and leave to adhere (2-24 h, depending on cell type).	Mix antibody and FabFluor	Add antibody-FabFluor mix (50 µL/well) to cell plate.
		Reagent at a molar ratio of	
		1:3 in media, 2x final assay	
		concentration. Incubate for 15	
		minutes to allow conjugation.	





- When the signal area was normalized

Capture images every 15-30

IncuCvte[®] for 24-48 hours Analyze using integrated

minutes (10x or 20x) in

software.



- Positive control wells (α -CD71) display a marked increase in fluorescence, whereas little or no signal was observed in the negative controls (IgG isotype); Z' values of 0.75 & 0.87.
- Of the 6 Abs, 3 (Ab 1a, 2 & 1b) produced a large signal and were detected at low concentrations (<0.05 μ g/mL).
- Ab 1a & 1b were the Ab clone from different suppliers and as expected produced similar responses.
- Abs 3, 4 & 5 were internalized only at higher test concentrations.



- Various FabFluor-pH-Abs were added to Jurkat or Raji cells (4 µg/mL).

con

- Both cell lines internalized CD71 and CD45 (general lymphocyte markers) but only Jurkat internalized CD3 (T cell marker) and not CD20, with the converse in Raji cells (B cells).
- Data correlates with known CD surface marker expression in these cell lines providing strong confidence in the signal specificity and generic utility of the method.

