

Design of Experiments for early phase process development of Antibody Drug Conjugates

Xavier Despinoy & Erik Johansson
06 May 2021

The logo for Sartorius, consisting of the word "SARTORIUS" in a bold, black, sans-serif font, centered within a bright yellow rectangular background.



SARTORIUS

- Antibody
- DOE software MODDE expertise



 **Piramal**
Pharma Solutions

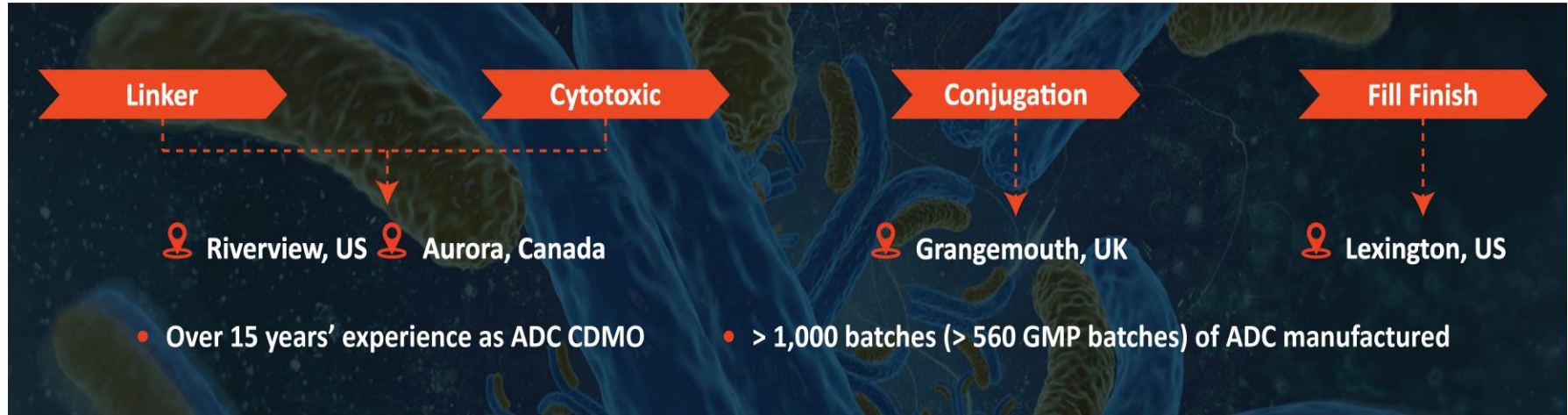
- Drug Linker
- Bioconjugation expertise

- Introduction to Piramal bioconjugation services
- Introduction to ADC Development
- Case Study: Process development of an ADC Reactive stage
 - Process
 - Considerations for DOE
 - Parameters/responses
- Design Of Experiments

- Piramal Grangemouth is a world leading Bio-conjugation CDMO
- >15 years experience in Antibody Drug Conjugation and Bio-conjugation
- Support Proof of Concept (milligram scale) through to Bulk Drug Substance Commercial manufacturing
- Highly skilled workforce across Development, Manufacturing and Quality Units



- Supply chain simplification for ADCs & bioconjugates



- Piramal GMP manufacture:
 - 40 distinct conjugates
 - 2 commercial products



Introduction to ADC Development

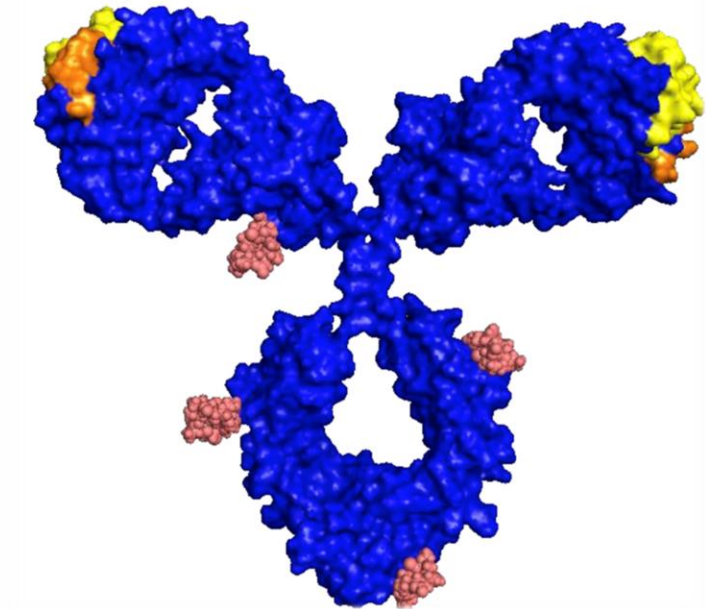
- Develop scientifically sound analytical methods suitable to support pre-clinical and ultimately, clinical release and stability testing of ADC
- Develop process conditions to meet key quality attributes for the ADC
- Have sufficient understanding of process robustness to enable safe scale-up
- Establish control strategy



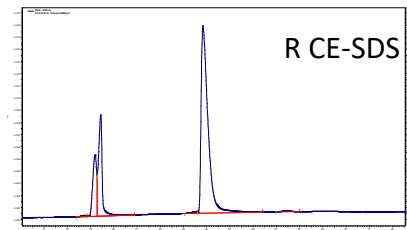
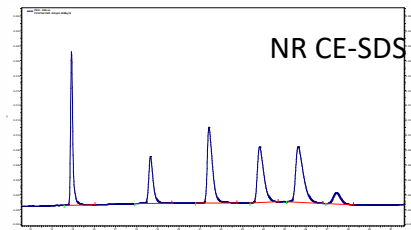
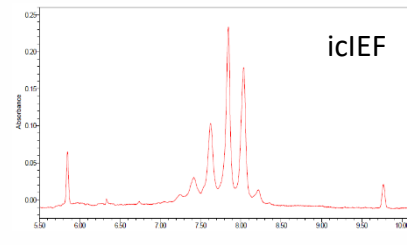
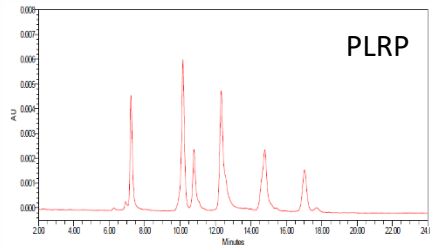
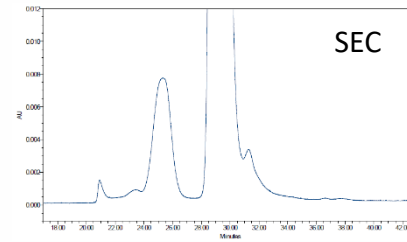
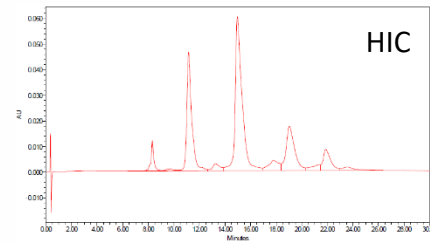
key components:

- Monoclonal antibody highly specific to target cell antigen
- Anticancer drug (payload) highly potent for cell killing activity
- Linker to covalently join payload to antibody

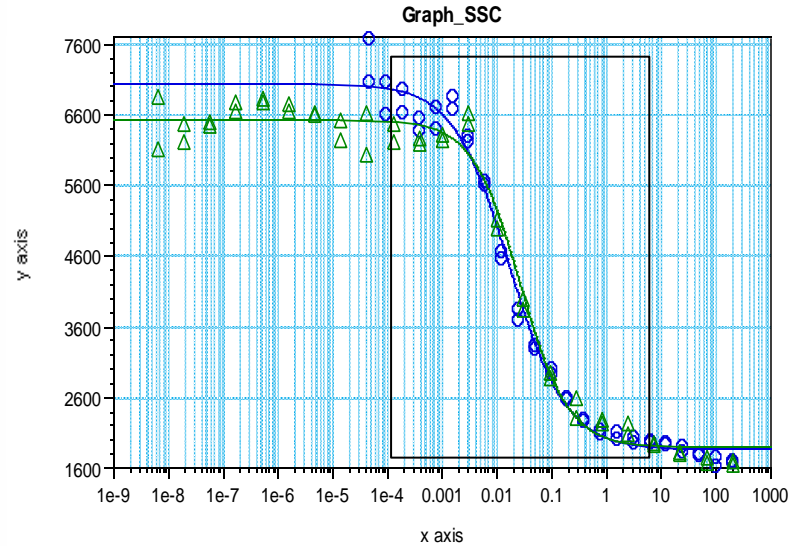
“ ADCs Bring together the best features of
Antibodies &
Cytotoxic drugs “



- Analytical complexity: antibody + payload + conjugate
- Methods developed immediately for key quality attributes: SEC, DAR and distribution (HIC, PLRP) and icIEF to support quick Process Development start
- Free drug and CE-SDS (R+NR) also initiated at early stage



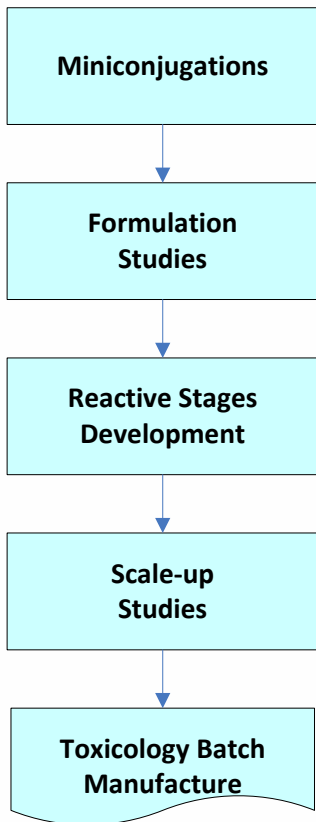
- Antibody Functional assays
 - ELISA (binding)
 - Potency (cell killing assay)
 - Effector functions
- Significant development required
 - Methods development starts early



4-P Fit: $y = (A - D) / (1 + (x/C)^B) + D$:

	A	B	C	D	R ²
○ std (Standards: conc vs OD)	7.04e+03	0.842	0.0168	1.87e+03	0.99
△ Test1 (Test1: conc vs OD)	6.53e+03	0.987	0.0261	1.9e+03	0.988

Curve Fit Option - Fixed Weight Value



- 5 Stages of development to go to GMP
- Approach designed to speed up overall program
 - (Analytical Development throughout Process Development)
 - Formulation Studies & Reactive Stages run in parallel
- Specific set of activities for each stage
- Final process with purification tested at scale prior to GMP manufacture

Case Study :

Use of DOE to develop the reactive stages of an ADC process

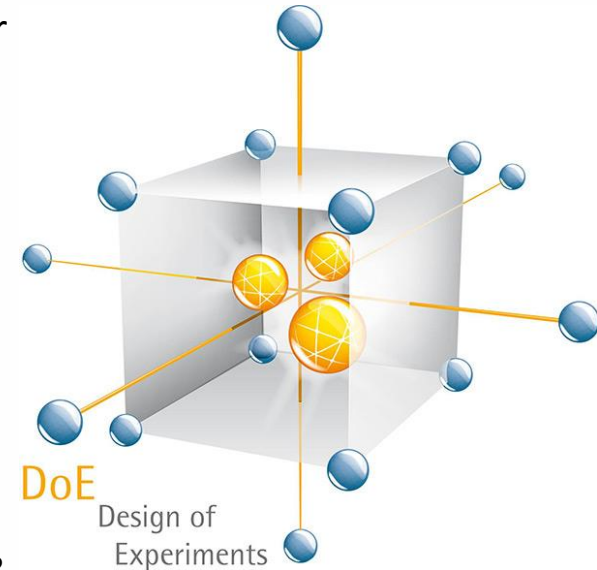


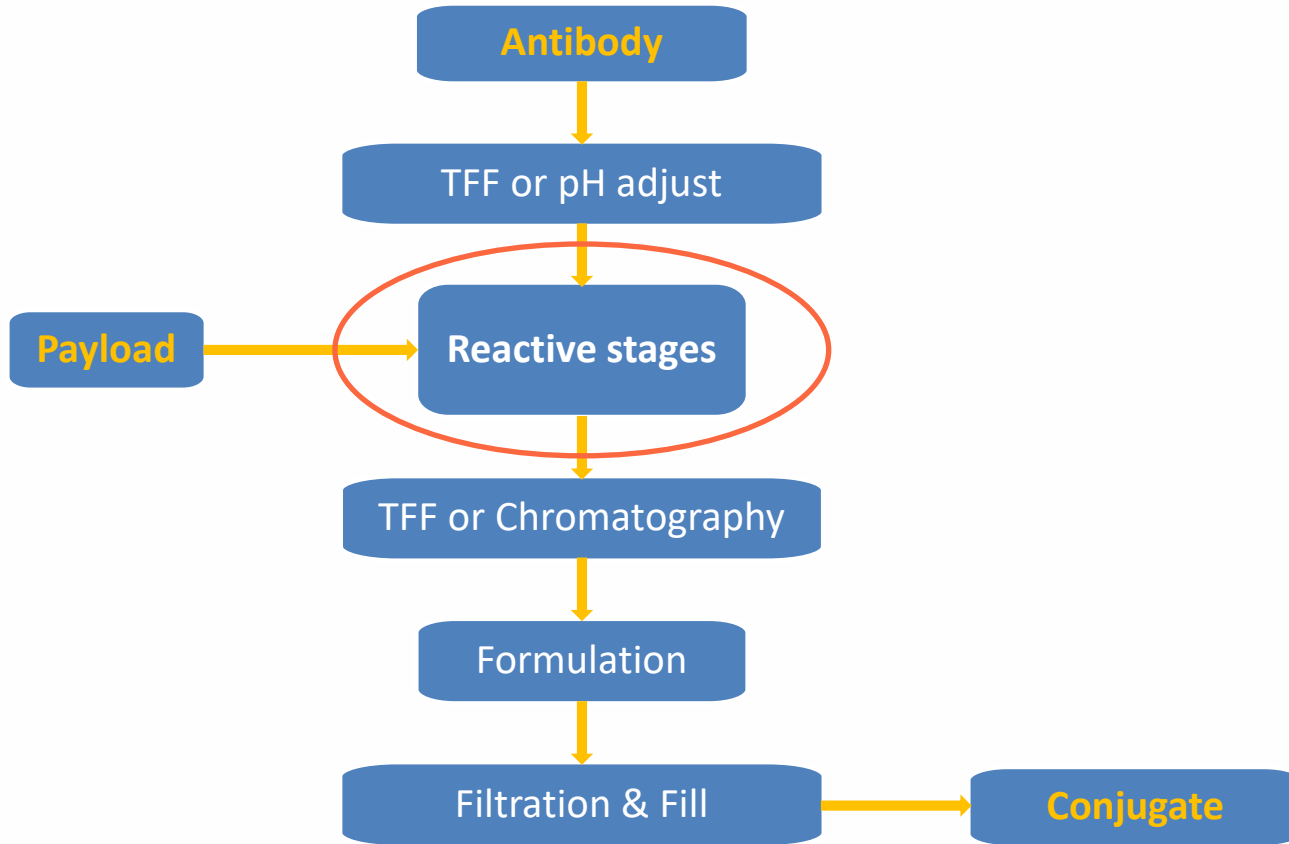
Why DoE?

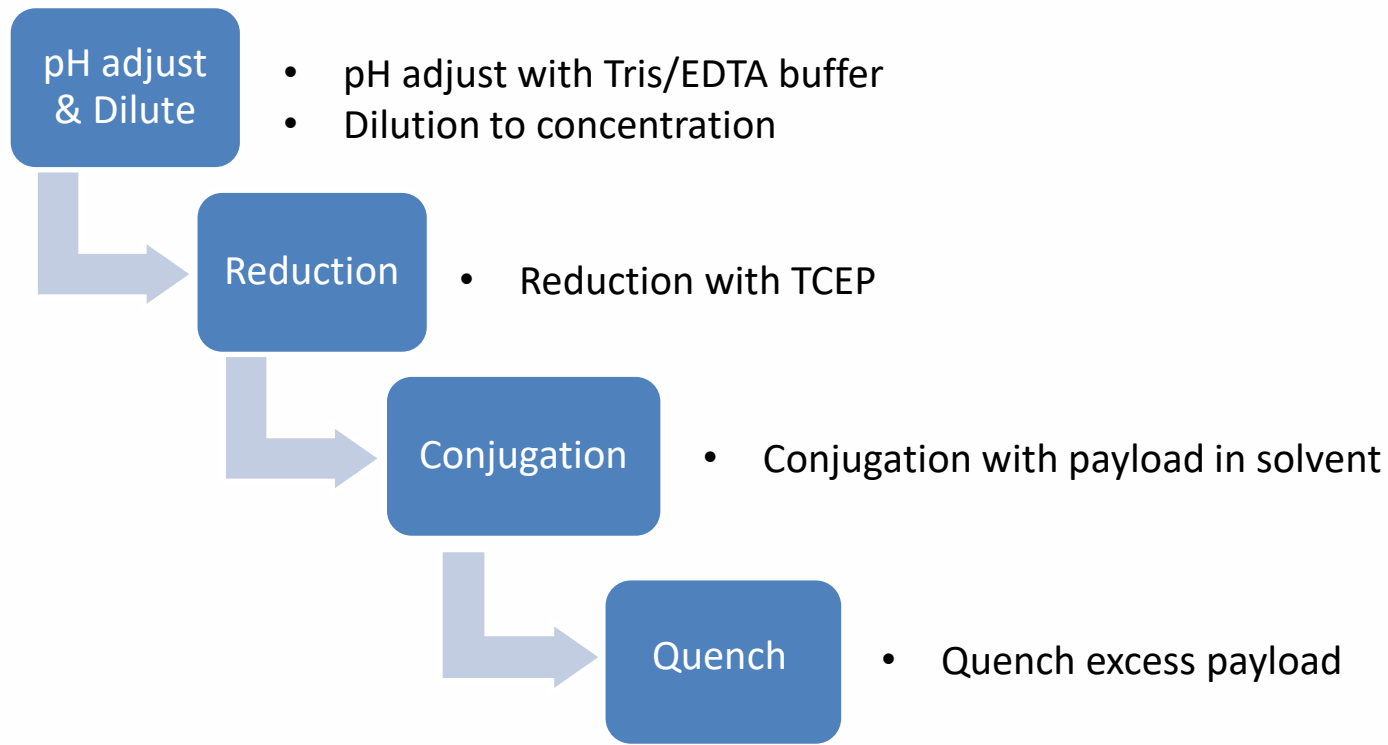
- Maximizes the information content while keeping the number of experiments low
- Allows systematic assessment of effect of multiple Factors (and their interactions) on Responses
 - Factor = Process Parameter
 - Response = Critical Quality Attribute
- Allows definition of a “Design Space” – safe operating conditions with CQAs meeting targets/ranges.

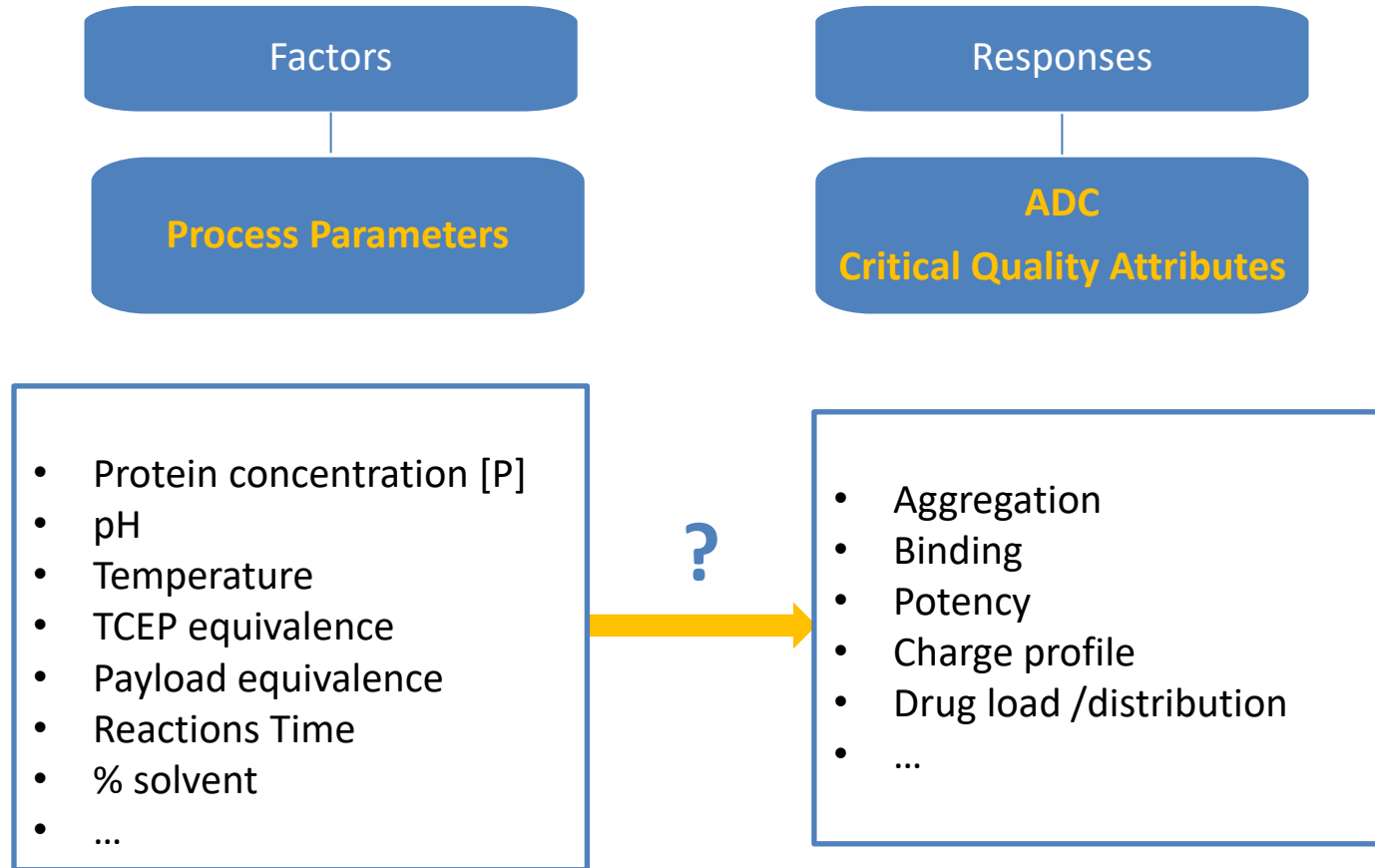
When and where ?

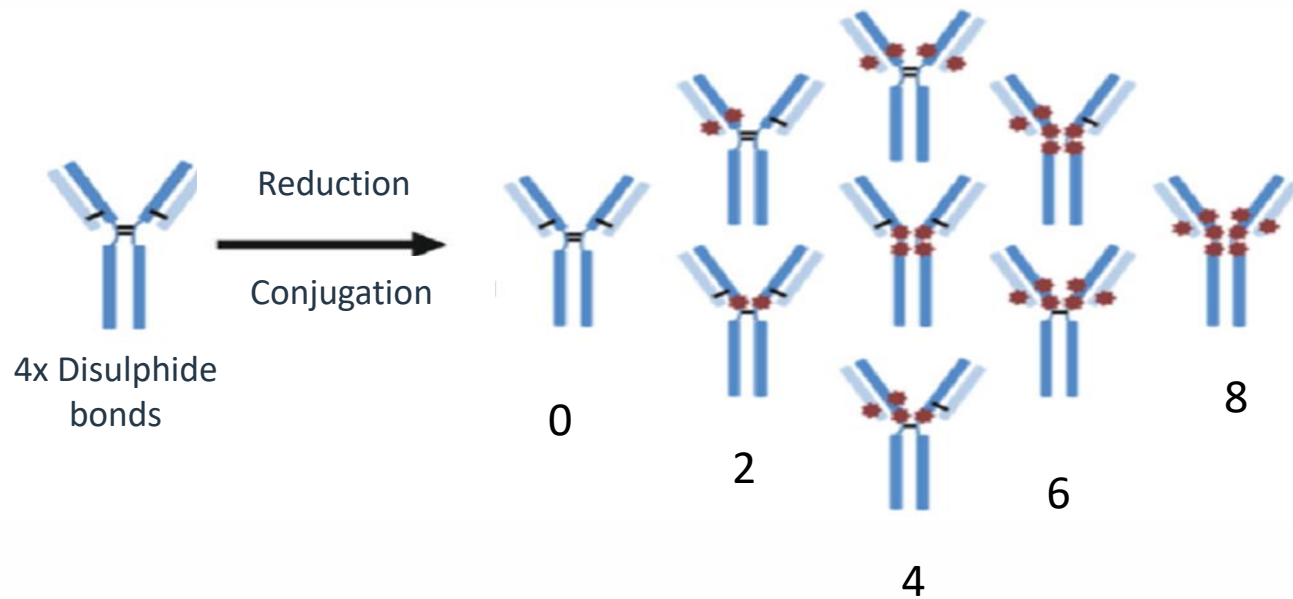
- DoE as an early Development tool to support ADC analytical & process development
- DOE for later phases to define the Design Space for parameters during pre-commercial Process Characterisation studies





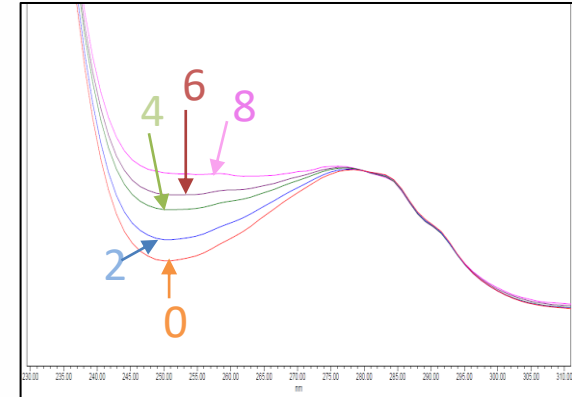
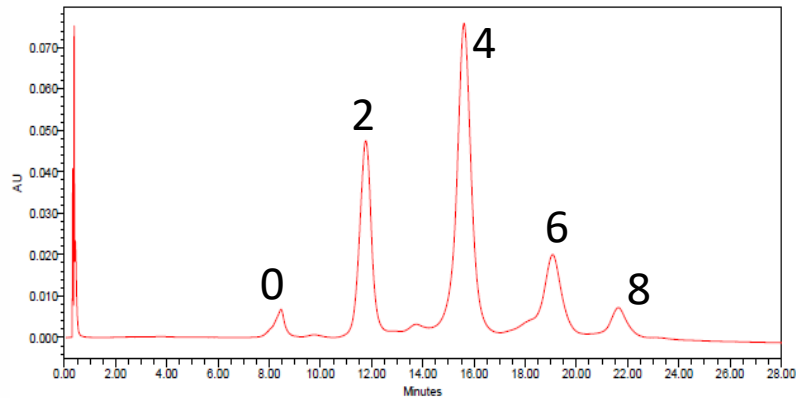






- Heterogeneous mixture of species
- Different number of “Drug or Payload” per molecule of Antibody

- Hydrophobic interaction (HIC) spectrum :
 - Distribution
 - Unconjugated antibody
 - Average DAR



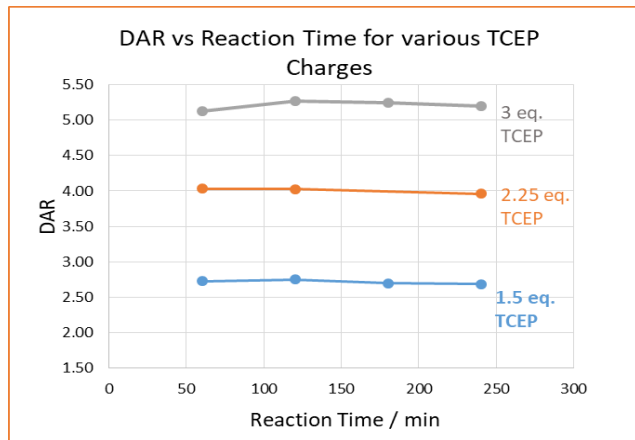
Average DAR – response for our study

Considerations prior to DOE

- Good analytics
- Parameters selection and range
 - Prior knowledge
 - Scouting experiments
 - Manufacture fit
- Statistical Design selection
- Preparation of input materials to design (eg pH)
- Use of scale-down model

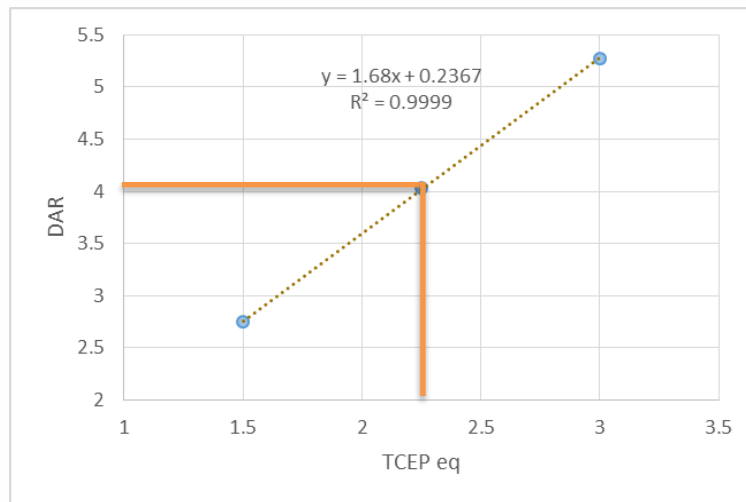
TCEP equ.	Time (h)	DAR	Time (h)	DAR
1.5	1	2.73	2	2.75
	3	2.70	4	2.69
2.25	1	4.03	2	4.03
	3	2.21*	4	3.96
3	1	5.12	2	5.27
	3	5.25	4	5.19

* undercharge of payload



- Factor evaluation :

- TCEP
- time



TCEP-DAR relationships

- Proposed DOE Factors and ranges :

Factors				
Name	Units	Low	High	Control Range (\pm)
Protein Conc.	mg/mL	5	15	1
Temperature	$^{\circ}$ C	16	26	2
pH		6.8	7.8	0.2
Reduction Time	min	60	180	30
Response				
Name	Abbrev.	Min	Target	Max
Drug Antibody Ratio	DAR	3.4	3.9	4.4

- Checked 'extremes' of proposed ranges :
 - All 'LOW' versus All 'HIGH'
 - Significant DAR variation 2.9 & 4.1 😊