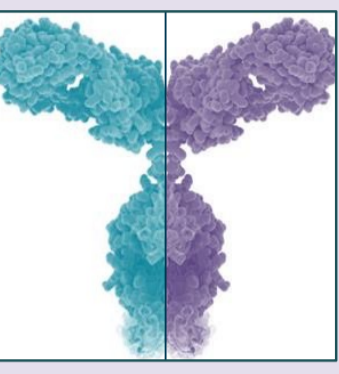


Charge variant microheterogeneity explored via icIEF and offline MS analysis of infliximab biosimilars

Nora Crushell¹, Anna Mulligan^{1,2}, Sara Carillo¹, Zarnab Yasmeen³, Jean-Francois Bellec³, Jonathan Bones^{1,2}

1. National Institute for Bioprocessing Research & Training, Fosters Avenue, Mount Merrion, Blackrock, A94 X099, Co. Dublin, Ireland
 2. School of Chemical and Bioprocess Engineering, University College Dublin, Belfield, Dublin 4, D04 V1W8, Ireland.
 3. Bio-Techne Ltd., 19 Barton Lane, Abingdon Science Park, Abingdon, OX14 3NB, United Kingdom



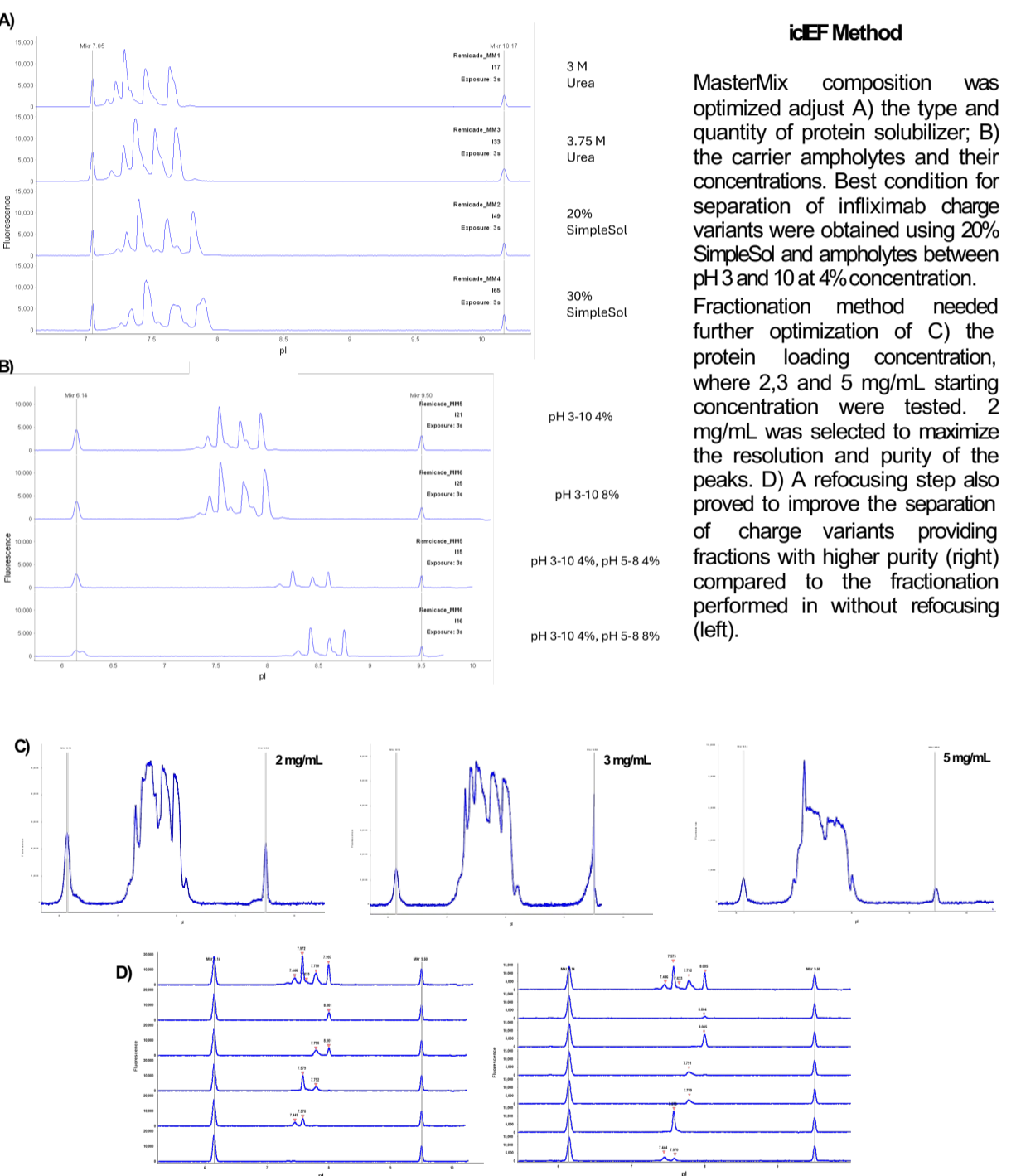
Introduction

Biosimilarity assessment requires a plethora of analytical assay to decipher the structural microheterogeneity of both target originator and candidate biosimilar. A compulsory assay performed in quality control laboratories is the charge variant analysis, based on the separation of proteoforms with different pI values. The gold standard for charge variant profiling is image-capillary isoelectric focusing (icIEF) although detection methods paired with this technology normally prevent a direct identification by mass spectrometry, due to matrix incompatibility or low recovery of each species. In this study we will develop a complete workflow using icIEF fractionation and analysis of each fraction by intact analysis using microchip CE-MS technology or bottom-up analysis by nano LC-MS/MS workflow using innovator (Remicade®) and biosimilar (Flixabi®) of infliximab as a model analytes¹.

About infliximab

- Infliximab is an anti-TNFα chimeric monoclonal antibody used against inflammatory diseases (Crohn's disease)
- MoA: it binds to TNFα inhibiting its binding to the receptor of immune cells
- Remicade (Janssen) was first approved by FDA in 1998 and by EMA in 1999
- Biosimilars started to appear on the market in 2013.
- 8 biosimilars now present on the market.

icIEF Method Optimization



Master Mix
 0.35% MC, 4% pH 3-10 ampholytes, 50 mM Arg
 20% SimpleSol
 1.5% pI markers 6.14, 9.50

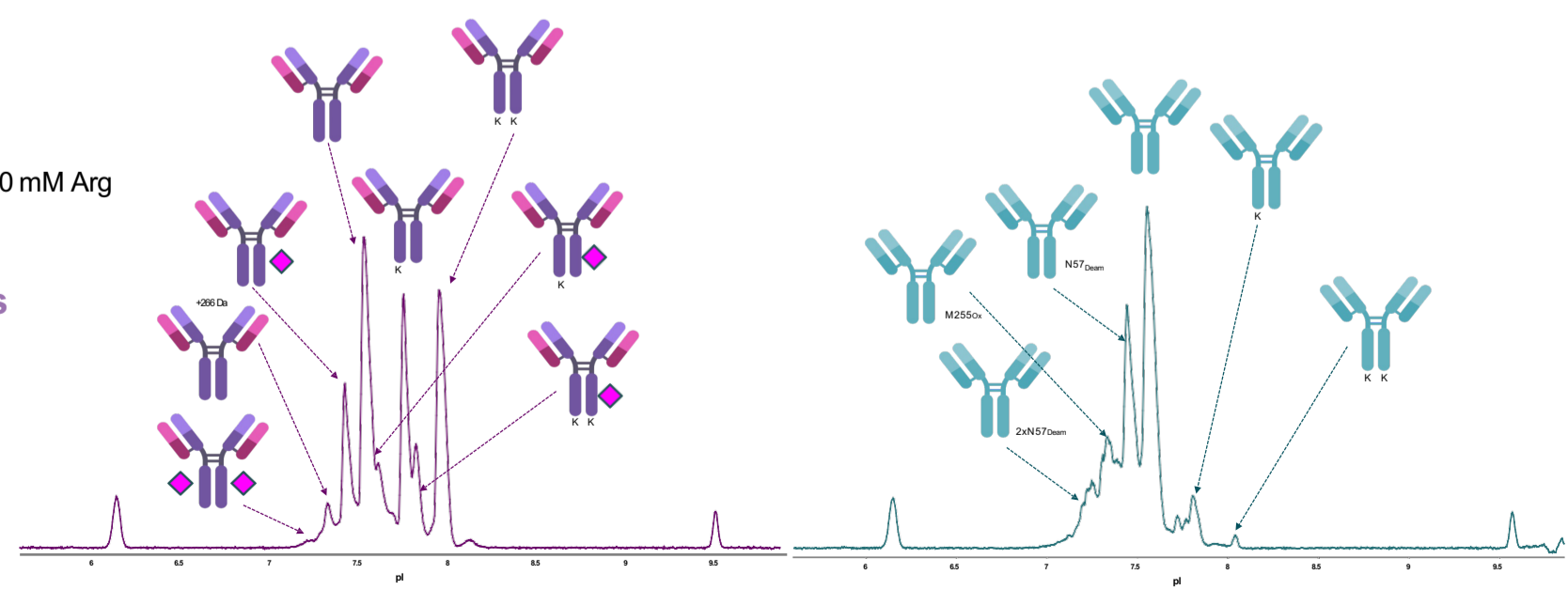
icIEF – Optimized Conditions
 Separation: 10 min 500 V, 10 min 1,000 V, 25 min 1,500 V
 Fluorescence detection: 0.2 sec
 Sample Load: 20 sec
 Mobilisation: 30 min 1,500 V
 Refocus: 5 min 1,500 V
 Fractionation 25 s, 1,500 V
 pI Markers 6.14, 9.50



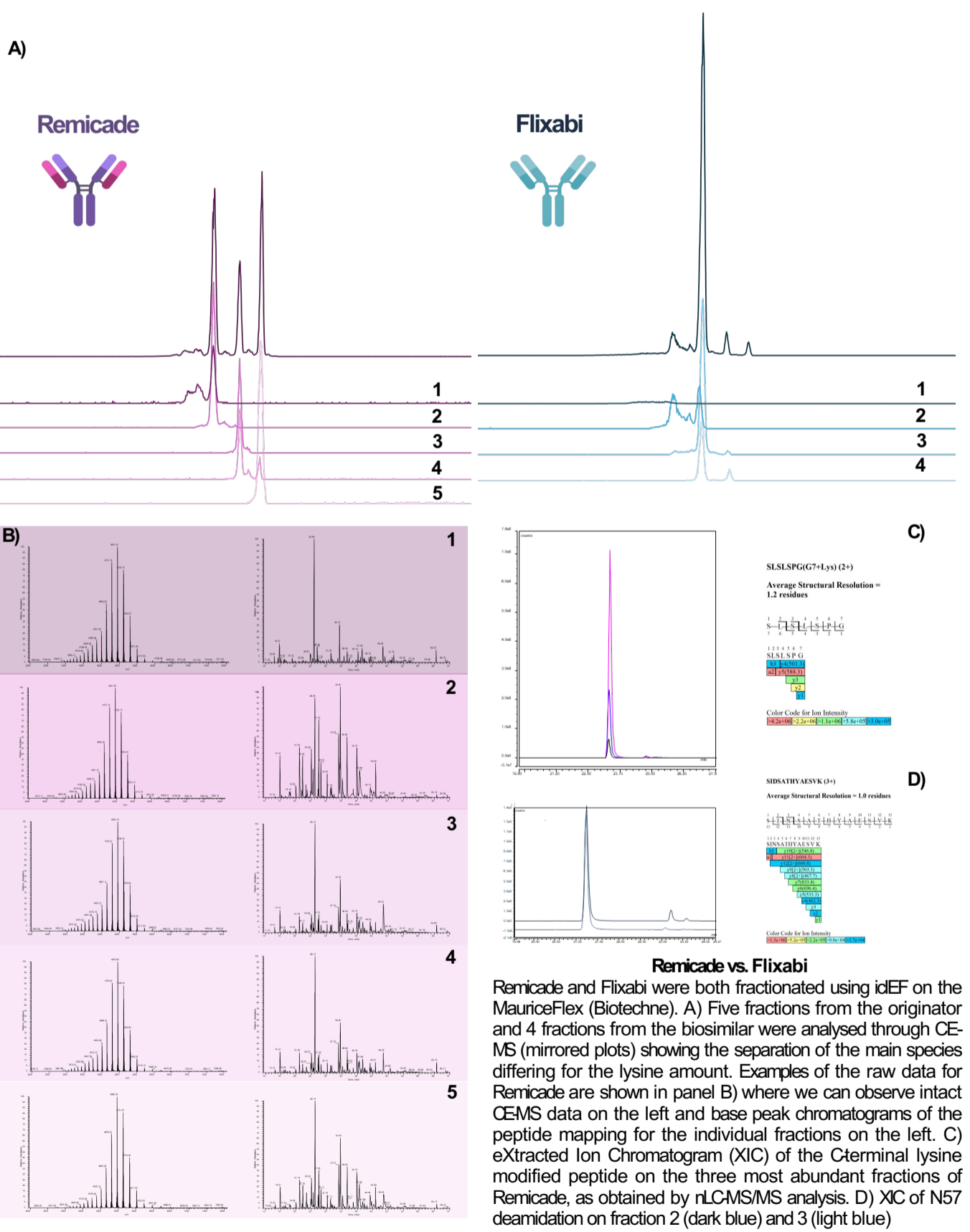
Offline fractions analysis



Separation and MS Settings	
CE-MS	CE sample transfer 10 µL, injection 2 nL, field strength 500 V/cm, pressure assist start time 0.5 min, replicate delay 20 s. MS: Application Mode: Intact Protein, High Pressure Tune Settings - Sheat gas: 2 arbitrary units (au), ion transfer tube: 300 °C Scan Parameters - Resolution: 30K (at m/z 200), Scan range: 4,000-8,000 m/z, RF lens 200%, acquisition gain control (AGC) 100%, Maximum injection time: 20 ms, microscan: 3, in-source CID: 135 eV
nLC-MS/MS	nLC: A – 0.1% formic acid in MS grade water, B – 0.1% formic acid in MS grade acetonitrile Flow rate 0.250 µL/min. Gradient 2% - 35% B in 55 minutes, 40%-45% B in 5 minutes. Column wash 80% B for 5 min Column temperature 45 °C MS: Application Mode – Peptide Scan Parameters – Resolution 120K (at m/z 200), Scan range: 200-2,000 m/z, RF lens %70, AGC 100%, Maximum injection time 100 ms, microscan 1 MS/MS: DDA Top 5. Isolation window 1.2 m/z. Collision energy (HCD) 28%, Resolution 15K (at 200 m/z), AGC 100%, maximum injection time 250 ms, intensity threshold 5e3, dynamic exclusion 7s, charge states 2-7



Results



Conclusions

- Charge variant analysis remains a key analytical tool to solve the complex structural heterogeneity of biotherapeutics.
- icIEF is a robust and reliable platform for charge variant analysis and the fractionation feature of the new MauriceFlex allows direct offline MS analysis using both intact and bottom-up approach.
- Although the differences highlighted in infliximab biosimilars are not clinically meaningful, the icIEF fractionation allowed to confidently distinguish acidic variants generated from sialylated N-glycans and acidic variants derived from deamidation events in the CDR region.
- The fractionation of species not allowing easy identification with commonly known modifications allows further studies to be carried for complete analysis.

1) Hong J. MAbs. 2017 Feb/Mar;9(2):364-382. doi: 10.1080/19420862.2016.1264550.
 2) Carillo S. J Pharm Biomed Anal. 2020 Jun 5;185:113218. doi: 10.1016/j.jpba.2020.113218.
 3) Smith J. Eur J Pharm Biopharm. 2023 Aug;189:276-280. doi: 10.1016/j.ejpb.2023.06.019.

