

Increasing Dynamic Binding Capacity of Oligo dT Using CIM 96 Well Oligo dT Plates

Nina Mencin, Andreja Krušič, Jure Ličen, Sebastijan Peljhan, Jana Vidič, Urh Černigoj, Tomas Kostelec, Aleš Štrancar, Rok Sekirnik*

BIA Separations d.o.o., A Sartorius Company, Mirce 21, 5270 Ajdovščina, Slovenia
* Corresponding author: rok.sekirnik@biaseparations.com

Affinity Purification of mRNA

Affinity-based chromatographic isolation of mRNA is robust and simple, lending itself as a useful industrial platform. mRNA constructs typically contain a 3' polyA tail to increase stability in vivo, thereby affording the possibility of affinity purification using oligo deoxythymidinic acid (Oligo dT) probes covalently coupled to a solid support. Polyadenylated mRNA forms a stable hybrid with Oligo dT under high-salt conditions which is destabilized when the salt is removed, allowing mRNA to be released (Figure 1).

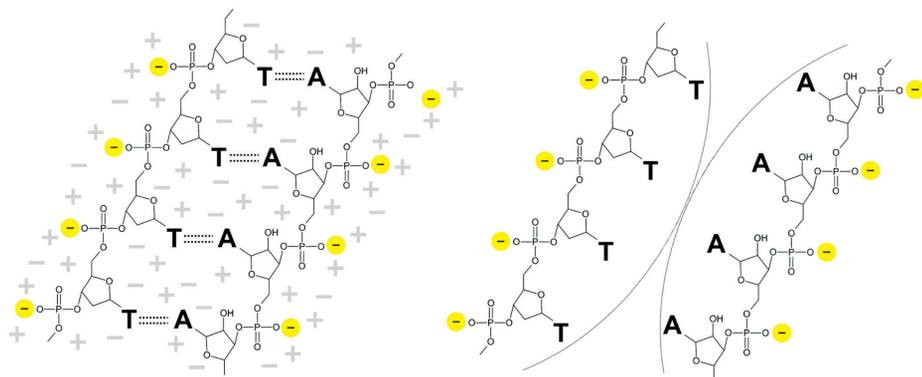


Figure 1: Binding Interactions Between mRNA Molecules and Immobilized Oligo dT.

Note: Left: High-salt conditions (binding), Right: Low-salt conditions (elution).

Typical dynamic binding capacity (DBC) of CIMmultus® Oligo dT for mRNA is 2-4 mg/mL. Higher IVT productivity will require higher binding capacities. Screening experiments to elucidate factors affecting CIMmultus® Oligo dT binding capacity for mRNA were performed in CIM® 96 well Oligo dT format (Figure 2). A simplified model identified NaCl, guanidine hydrochloride (Gu-HCl) and MgCl₂ concentration as the key factors contributing to DBC. Buffer chemistry, buffer pH, salt type and mRNA concentration had little or no effect on DBC.

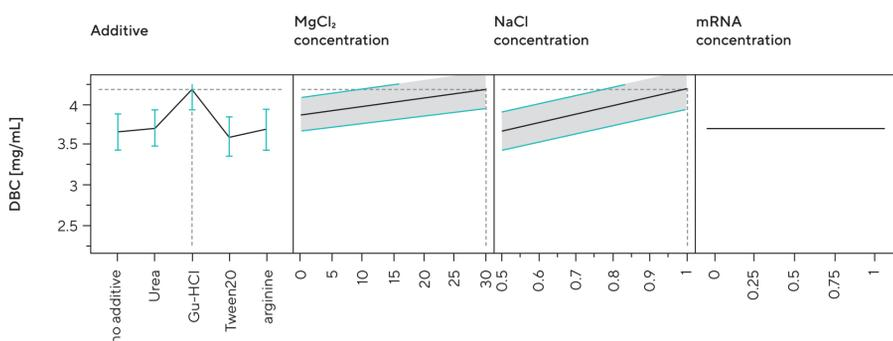


Figure 2: Prediction Graphs of Factors With Significant Contribution to Oligo dT DBC.

Note: Experimental results were fitted by a multiple regression model and refined with a backward selection approach with a p-value threshold of <0.05.

Conclusions

- CIM 96 well Oligo dT plates were used for optimization of DBC
- The main contributing factors to DBC were identified as NaCl, Gu-HCl, MgCl₂. DBC of >6 mg/mL can be achieved with Gu-HCl
- Due to higher chaotropicity of Gu-HCl compared to NaCl, stronger binding is achieved by reducing the hydration shell around, and thus minimizing repulsive interactions between, mRNA and Oligo dT

Guanidinium, Mg²⁺ and NaCl Effects on DBC

Combinatorial effect of Gu-HCl/NaCl and MgCl₂/NaCl was evaluated in 96-well format. Gu-HCl demonstrated higher impact than NaCl; DBC of 6 mg/mL was reached (Figure 3 a). Above 1 M Gu-HCl/1 M NaCl, precipitation of mRNA was observed. Effect of Mg²⁺ was also positive, though less pronounced, reaching DBC of 4.5 mg/mL (Figure 3 b). Gu-HCl alone was titrated as a loading salt, resulting in dose-response between 0.1-1 M Gu-HCl with DBC 6.4 mg/mL for 1 kb mRNA (Figure 3 c) and 3.8 mg/mL for 4 kb mRNA (Figure 3 d).

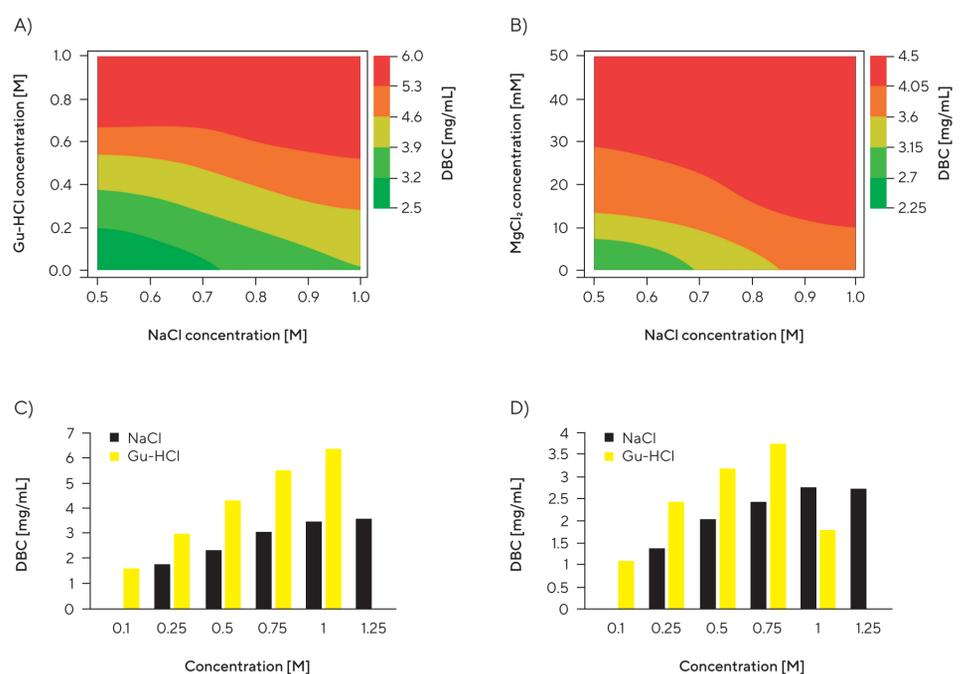


Figure 3: Contour Plots of Oligo dT Dynamic Binding Capacity When mRNA Is Loaded In a) Gu-HCl/NaCl, b) MgCl₂/NaCl Combination. Titration of Gu-HCl as Loading Salt for Binding Of c) eGFP (1000 nt) or d) mRNA Encoding a Proprietary Sequence (4000 nt) to Oligo dT 96 Well Plates.

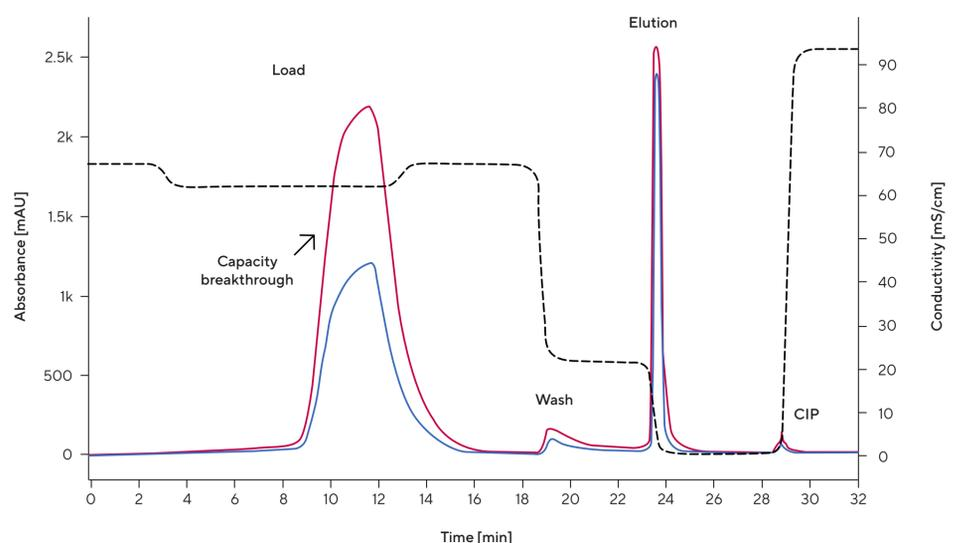


Figure 4: Chromatogram of mRNA DBC Determination on CIM® Oligo dT (0.1 mL) Column in Presence of 0.75 M Gu-HCl. Red: UV 260 nm; Blue: UV 280 nm

Effect of Gu-HCl was then transferred to a chromatographic separation mode using CIM® Oligo dT column (0.1 mL) confirming DBC of 5.5 mg/mL (Figure 4). A scale up to CIMmultus® Oligo dT (1 mL) column confirmed DBC of 5.5 mg/mL at 0.75 M Gu-HCl.