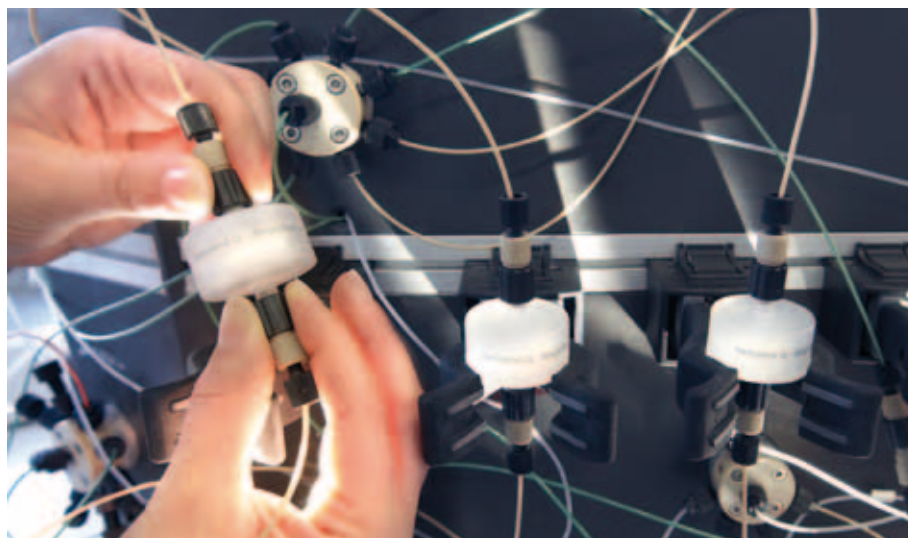




Host cell protein removal

A comparison between Sartobind STIC® PA and Sartobind Q



Materials

Membrane	1) Sartobind STIC® PA 2) Sartobind Q
Membrane ligand	1) Primary Amine (PA) 2) Quaternary ammonium (Q)
Membrane volume (area)	0.41 mL (15 cm ²)
Number of membrane Layers	3
Load material (HCP) concentration	500 ppm (500 ng HCP/mg MAb)

Introduction

In anion exchange flow-through (FT-AIEX) applications, membrane chromatography offers higher throughput, less buffer consumption, and more convenient handling than traditional bead columns. Conventional quaternary amine based chemistries – although established as a standard FT polishing step in Monoclonal Antibody (MAb) purification processes – typically require low feed conductivity. This often involves dilution of feedstreams and can result in facility fit limitations when high titer processes are accommodated in existing plants¹. Furthermore, high impurity levels limit load densities and would require larger membrane adsorber volumes and lead to increased production cost. To address these limitations and facilitate a wider design space for FT-AIEX membrane chromatography at commercial scale, Sartobind STIC® (Salt Tolerant Interaction Chromatography), a novel membrane concept based on weak anion exchange chemistry was developed². The Sartobind STIC® PA (primary amine) anion-exchange membrane is composed of a primary amine ligand that is attached to a cross-linked, regenerated macroporous cellulose base matrix.

The superior performance of Sartobind STIC® PA membrane at higher conductivities can be attributed to the ligand and ligand density. It is important to note that while the working pH of quaternary amine chemistries are ~pH 8, the weak anion exchange ligand of Sartobind STIC® PA may require operating at a lower pH for optimal impurity clearance.

Method

A MAb feedstream produced in CHO (Chinese Hamster Ovary) cell culture was processed through affinity and cation exchange chromatography steps before being processed through either Sartobind STIC® PA or Sartobind® Q under flow-through conditions. Sartobind® Q was loaded at pH 8 and a conductivity of 7 mS/cm, while Sartobind STIC® PA was loaded at pH 8 and at conductivities of 7, 10 and 15 mS/cm. The MAb feedstream had a host cell protein (HCP) concentration of 500 ppm (500 ng per mg of MAb), and was loaded at a rate of 10 membrane volumes/minute, up to a load density of 10 kg/liter of membrane volume³. HCP levels were measured in the filtrate at various intervals.

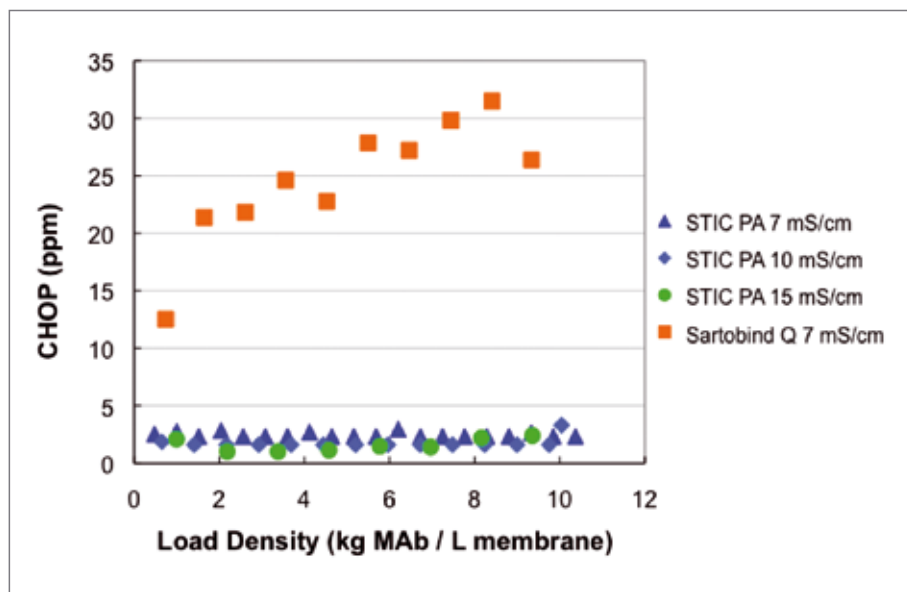


Fig. 1: HCP concentration in Sartobind STIC PA and Sartobind Q pools as a function of load density at various conductivities. Sartobind STIC PA shows higher HCP clearance than Sartobind Q at conductivities up to 15 mS/cm.

Results

Figure 1 shows the superior HCP clearance with Sartobind STIC® PA as compared to Sartobind® Q. At a load condition of pH 8, 7 mS/cm, Sartobind® Q showed increasing levels of HCP in the pool through the entire duration of load while Sartobind STIC® PA cleared HCP to less than detectable at a load density up to 10 kg/L. At higher conductivity conditions, 10 and 15 mS/cm, Sartobind STIC® PA also cleared HCP to less than detectable at a load density up to 10 kg/L. This exemplifies the ability of Sartobind STIC® PA to operate at high conductivity conditions without adverse impact to HCP clearance.

Summary

Sartobind STIC® PA is able to overcome the limitations of HCP removal at high conductivities by utilizing a primary amine ligand that is attached to a cross-linked, regenerated macroporous cellulose base matrix. This novel combination results in higher binding capacities of impurities at high conductivities, which may eliminate facility fit limitations caused by feedstream dilution when using adsorbers that necessitate operating at low conductivities.

References

1. Faber, R., Yang, Y. & Gottschalk, U. Salt tolerant interaction chromatography for large-scale polishing with convective media. *BioPharm Int. Suppl.* 11–14 (2 October 2009).
2. Fraud, N., Shomglin, K., Faber, R., van Reis, R., Gottschalk, U. & Mehta, A. New membrane adsorber for polishing at high salt concentration. *Recovery of Biological Products XIV Conference Lake Tahoe, California, USA, 1–6 August 2010*
3. Fraud, N., Faber, R., Mehta, A., Tully, T., Ultra small membrane adsorbers for process development & screening of new membrane adsorbers. *238th ACS Meeting, Washington August 16–20, 2009.*

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen, Germany

Phone +49.551.308.0
Fax +49.551.308.3289
www.sartorius-stedim.com

USA Toll-Free +1.800.368.7178
UK +44.1372.737159
France +33.442.845600
Italy +39.055.63.40.41
Spain +34.90.2110935
Japan +81.3.3740.5407

Sartobind®, and Sartobind STIC® are trademarks of Sartorius Stedim Biotech GmbH
Technical data are subject to change without notice.
Printed in Germany on paper that has been bleached without any use of chlorine.

First published:
May 2011

Publication No.: SL-4060-e11051
Order No.: 85032-540-18
Ver. 05 | 2011