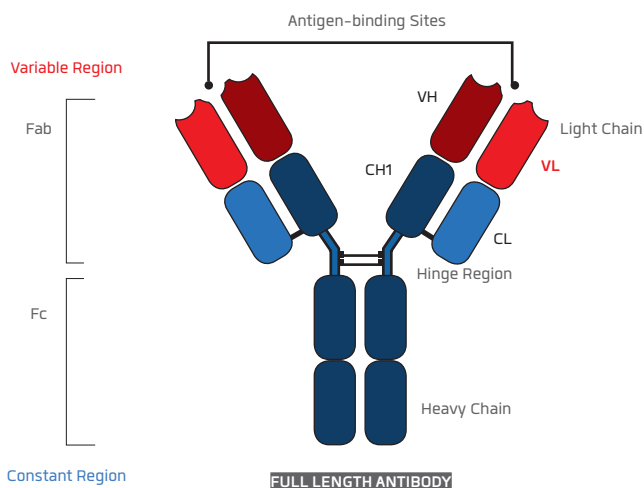


Protein A Affinity Chromatography Resin

KANEKA KanCapA™

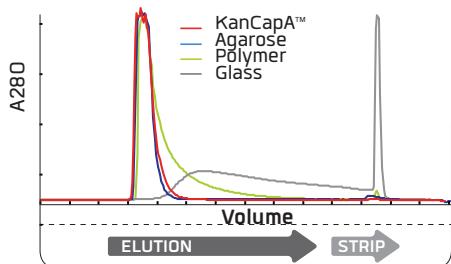
- > Mild acidic elution conditions for all mAbs
- > High flow rate operation and easy scale up from pilot to process
- > Track record of successful mAb GMP manufacturing for clinical trial



KANEKA KanCapA™ resin is composed of a proprietary recombinant alkaline stable Protein A ligand and an innovative highly cross-linked cellulose base matrix.

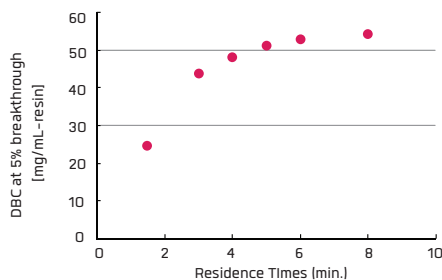
KANEKA KanCapA™ is designed for industrial-scale purification of monoclonal antibodies. The resin is manufactured under a validated GMP process to ensure a stable and reliable supply chain.

Mild Acidic Elution

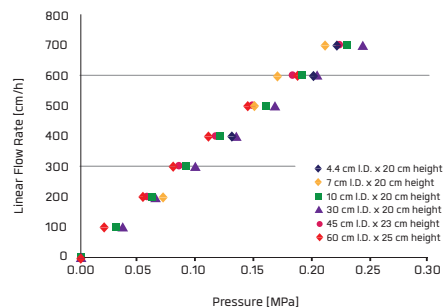


mAb A (VH3 subfamily) elution pH 3.5

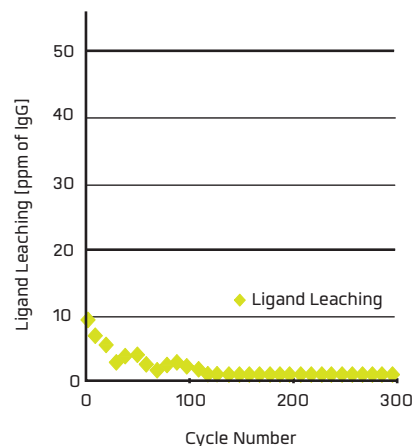
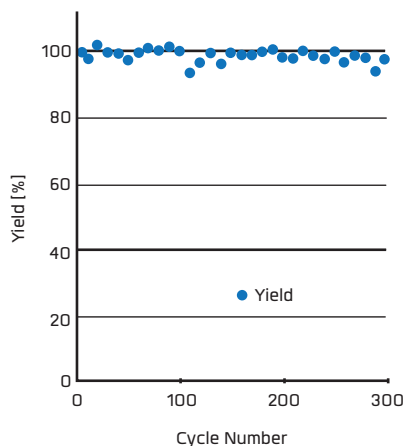
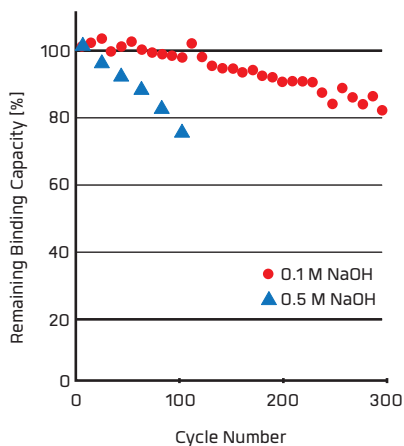
High Binding Capacity



High Flow Rate Operation and Easy Scale Up from Pilot to Process



Alkaline Stability and Long Lifetime



Contact information

Japan

Kaneka Corporation
1-12-32, Akasaka, Minato-ku,
Tokyo 107-6028, Japan
bioproducts@kaneka.co.jp

USA

Kaneka US Innovation Center
7979 Gateway Blvd., Suite 220
Newark, CA 94560
bioproducts@kaneka.com

Europe

Kaneka Eurogentec S.A.
Rue du bois Saint-Jean, 14
4102 Seraing, Belgium
bioseparation@eurogentec.com

