

Protein A Affinity Chromatography Resin

KANEKA KanCapA™ 3G

Resin with advanced performance

- > Enhanced binding capacity
- > Excellent elution profile
- > Unique impurity removal properties



KANEKA KanCapA™ 3G Formats

KANEKA KanCapA™ 3G represents a new cellulose-based protein A resin with a high dynamic binding capacity and an outstanding impurity removal characteristics.

KANEKA KanCapA™ 3G is the best choice for purifying monoclonal antibodies and their Fc fusion derivatives under milder elution conditions with increased purity.

Enhanced binding capacity

KANEKA KanCapA™ 3G shows enhanced binding capacity, mainly at longer residence times, which makes it suitable for monoclonal antibody purification from high titer feedstocks.

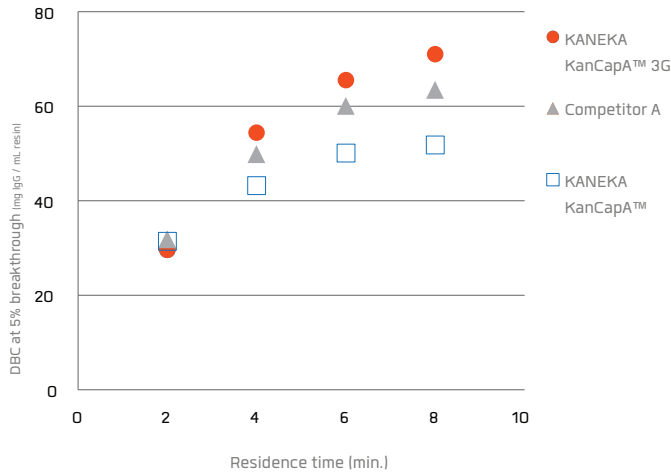


Figure 1 Residence time vs DBC (Polyclonal Human IgG)

Mild elution pH

KANEKA KanCapA™ 3G's ligand was designed for efficient elution of monoclonal antibodies and their Fc fusion derivatives under mild acidic conditions.

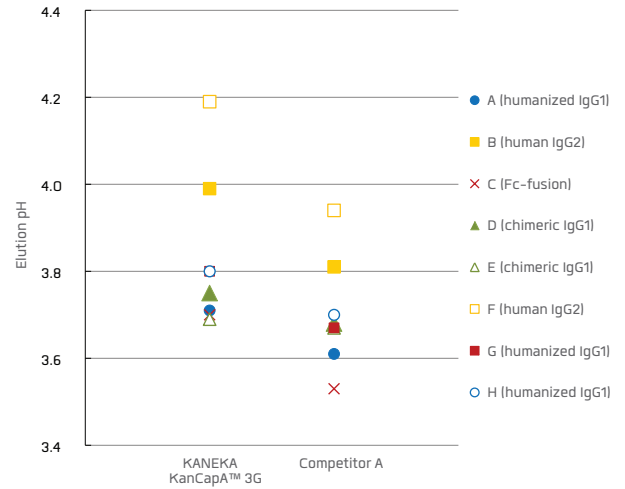


Figure 2 Comparison of the elution pH of KANEKA KanCapA™3G to competitor A. Elution pH was determined by a linear pH gradient elution using citrate buffer from pH 6 to pH 3.

Unique impurity removal properties

KANEKA KanCapA™ 3G has a better potential for impurity removal compared to current resins available on the market.

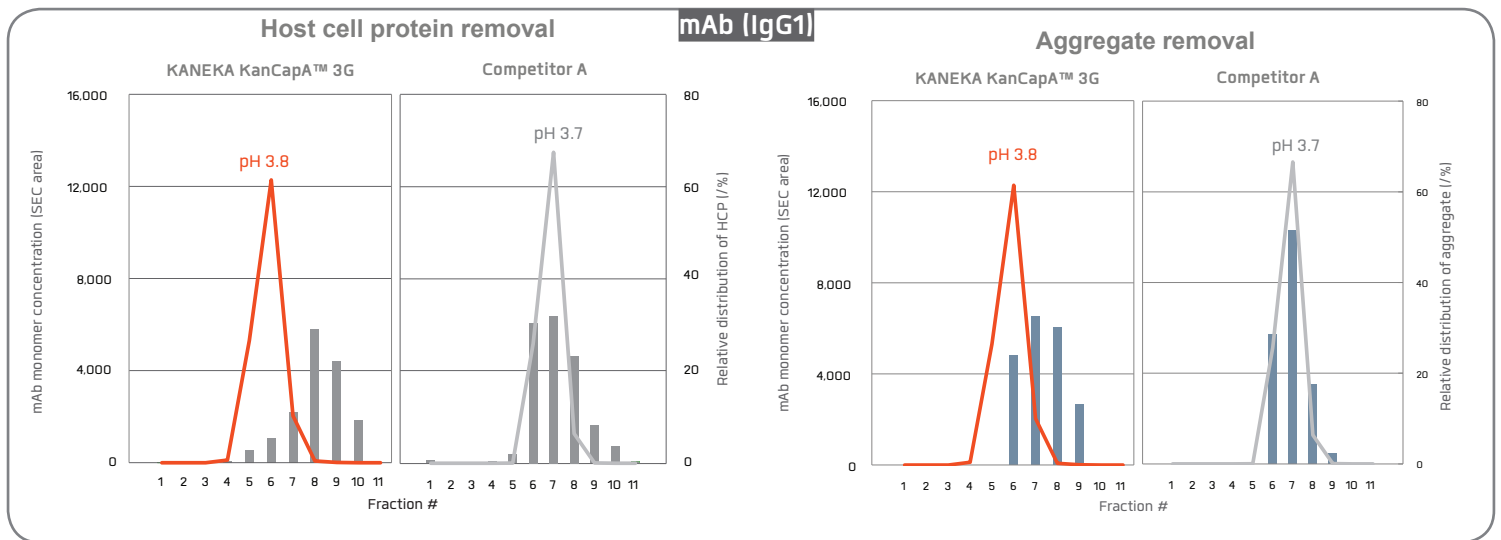


Figure 3 Efficient HCP (left) and aggregates (right) separation during mAb elution by linear pH gradient.

■ Monomer Peak ■ Host cell protein (HCP) ■ Aggregate **Molecules:** IgG1 (CCCF), Load: 5 mg/mL-resin **Elution:** 50 mM Citrate buffer, (pH 4.5→3, gradient, 10 CV)



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

