

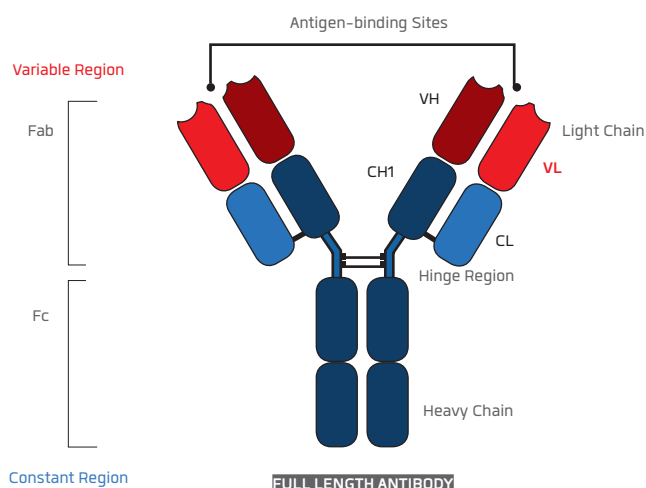
KANEKA KanCap™ L

- > High binding capacity for antibody molecules containing a κ light chain
- > Wide binding range for antibody molecules containing a κ light chain
- > High purification

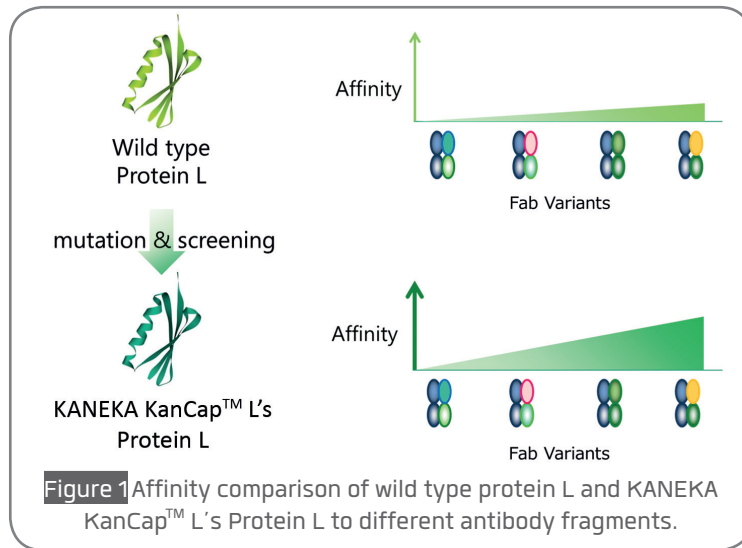


Kaneka proposes an antibody purification platform of antibody fragments using Kaneka's new Protein L resin.

KANEKA KanCap™ L is a newly developed cellulose-based Protein L resin designed for capture and purification of a wide range of antibody fragment formats such as scFv, Fab and F(ab')₂. It consists of a highly cross linked cellulose matrix and a newly developed Protein L ligand that exhibits an increased affinity to VL region of a κ light chain. It is also an alternative tool to purify full length antibody formats containing a κ light chain which poorly binds to Protein A resins.



Multiple Protein L mutants were screened to select the one with the widest binding spectra and the highest binding affinity to Fabs.



Excellent purification tool for antibody fragments

