

Protein Production in Pichia: High yield and Low O-glycosylation Platform

Nathalie Pirlot², Scientist, New Technologies Development, Eurogentec Biologics Toru Watanabe¹, Biologics Research Groupe Biotechnology Research Laboratories, Kaneka

¹Kaneka Corporation, 1-8, Miyamae-cho, Takasago-cho, Takasago, Hyogo 676-8688, Japan; ²Kaneka Eurogentec S. A., 14, rue Bois Saint-Jean, 4102 Seraing, Belgium



Kaneka Eurogentec is a well-established, FDA inspected, full-service biopharmaceutical CDMO producing recombinant proteins by microbial fermentation for preclinical, clinical and commercial uses. Together with Kaneka Corp., the company is continuously improving its technological expertise for recombinant proteins production through the development of Pichia pastoris strains and our multicopy strategy offers a powerful expression platform for low O-glycosylated proteins (available for licensing).

Pool of 1 copy clones



Kguekg

Eurogentec at a glance

One Stop Shop From Tox/API to Commercial

Eurogentec is a well established, FDA inspected, full-service CDMO producing GMP plasmid DNA, RNA and recombinant proteins for preclinical, clinical and commercial supply.





up to 4 vectors



Isolated 2-copy clone colonies

O-Glycosylation reduction in V_HH production

Combination of our optimized strains screening and our multicopy approach:

Positive impact on O-Glycosylation level and secretion yield of the Pol



This research was supported by AMED under Grant Number JP17ae0101004h0005





Projects

Geographical breakdown





CONTACT US

For CDMO services, contact us at BIO.BD@eurogentec.com





