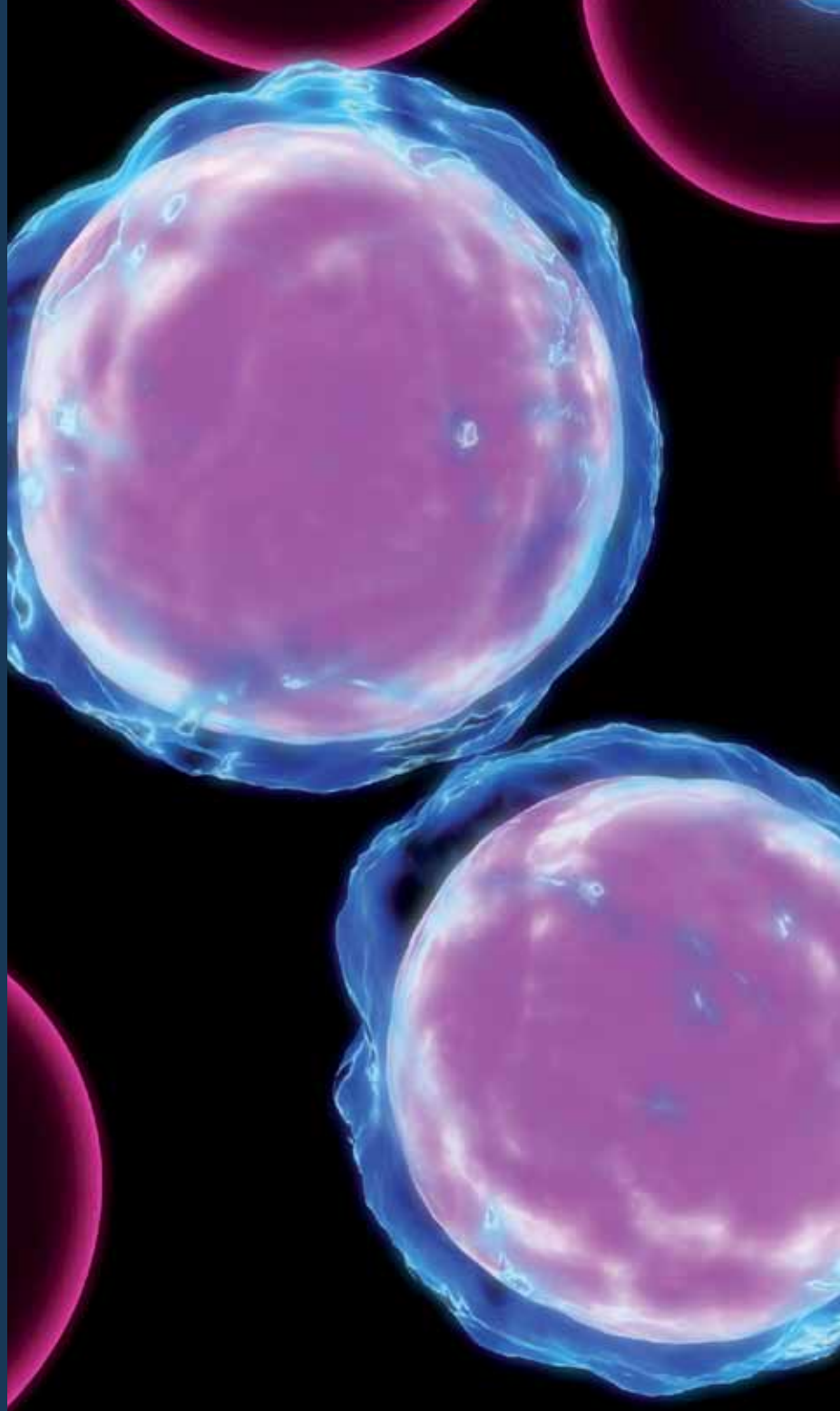


3

3.9 Custom Protein Expression



Custom Protein Expression

Whether you are interested in crystallography, NMR, protein-protein interactions, or activity assays, Eurogentec now offers a protein expression service.

You choose your protein of interest, we will produce it for you according to your specifications:

- ▣ Expression in *E. coli*
- ▣ Expression in HEK or CHO cells
- ▣ Choice of different protein tags
- ▣ Choice of purification levels from standard purity (85 %) to high purity (> 90 %)

Confidentiality

Eurogentec maintains all projects in the strictest confidentiality. Should you require a material transfer agreement please e-mail us at info@eurogentec.com

How to start

Fill in the order form on our website or send us on e-mail mentioning: protein.expression@eurogentec.com

- ▣ Your gene/protein sequence
- ▣ The host
- ▣ The tag
- ▣ The required amount of protein to be expressed
- ▣ The purification level

You will receive a sequence analysis report and a quotation by e-mail within the next 5 business days.

Project description

1. Each project starts with the codon-optimized gene synthesis
2. We clone it (optionally with the tag) into the best-suited vector for your chosen host.
3. We produce a pilot or small-scale reactor to determine the gene's expression level.
4. Based on previous results we move to the large-scale protein expression phase.

Guarantee

It may happen that your protein is too difficult to express. You will only be invoiced for the services completed. You are also free to exit the protocol at any step.

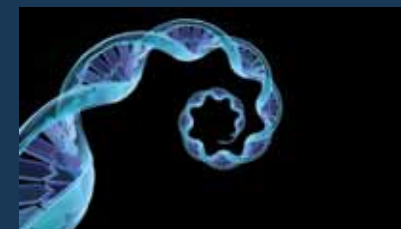
Protein expression in *E. Coli*

Description	References
Gene synthesis & cloning	GL-GENE-SYNTH
Small scale protein expression	GL-PEXP-PREP
Large scale protein expression >10 mg	GL-PEXP-10MG

Protein expression in CHO or HEK cells

Description	References
Gene synthesis & cloning	VB-GENE-SYNTH
Pilot scale protein expression	VB-PEXP-PREP
Large scale protein expression	VB-PEXP-PROD

Custom Genes



- ❖ Any Gene in Any Vector
- ❖ Up to 50 000 bp
- ❖ Gene Optimization
- ❖ Synthesis of Complex Genes
- ❖ 100 % Guaranteed Sequence