

2

2.2 Custom genes

113 Introduction

113 Case Study

114 Process Overview



Custom Genes

Introduction

Custom genes are synthetic genes optimized to face any hindrance encounter with classical PCR amplified nucleotide fragments.

Custom Genes synthesis considerably reduces time spending on oligonucleotides design, cloning process and sequencing. Moreover, the quality of genes coming from artificial synthesis is rather superior than that obtained for genes provided by classical PCR with no formation of all the mutations or Single Nucleotide Polymorphisms (SNP), which may cause troubles during PCR or other applications.

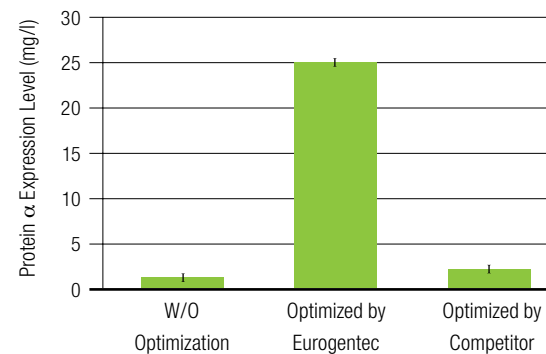
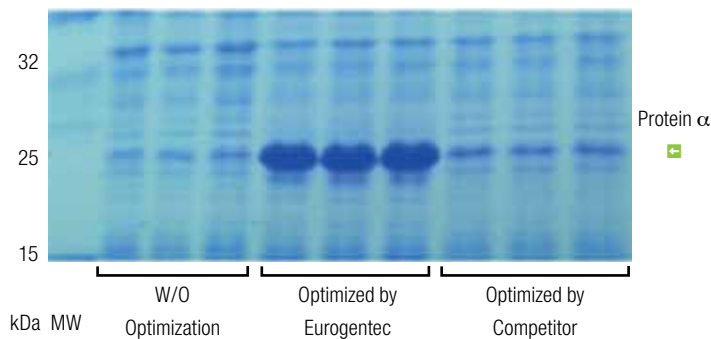
Custom Genes bring you both optimization and flexibility. Any synthesized genes can be cloned in any vector to lead to protein expression 100 fold higher than classical PCR amplifications.

With Custom Genes synthesize any natural, mutant or virtual sequences adapted to your needs!

Case Study

Increase your protein expression level in *E. coli* !

Eurogentec's gene optimization delivered up to 20 times higher protein expression levels compared to non-optimized native gene sequence.



For example the expression level of protein α was 13 times more than that from Competitor's optimization method.

Specifications

- ▣ Any Gene in Any Vector
- ▣ Highly complex genes
- ▣ Up to 50 000 bp
- ▣ Gene optimization
- ▣ 100 % Guaranteed Sequence
- ▣ Fast Turnaround Times
- ▣ Trusted Quality

Process Overview

