

FectoPRO®

Transfection Reagent for Protein and Monoclonal Antibody Production



Product Information

FectoPRO® is a transfection reagent specifically designed for enhanced transient gene expression in suspension CHO and HEK293 cells cultured in various serum-free media. FectoPRO®-mediated transfection processes are easily scalable, ensuring robust, reproducible protein production. The increased protein yields achieved with less plasmid DNA (pDNA) make FectoPRO® a cost-effective solution.

Features and Benefits

- **Productive:** High transfection efficiency and supports productivity
- **Cost-Effective:** Enhanced production yields and lower pDNA requirement
- **Versatility:** Compatible with various expression media and cell systems, adapted to high-density cell systems

Introduction

Relevant Applications

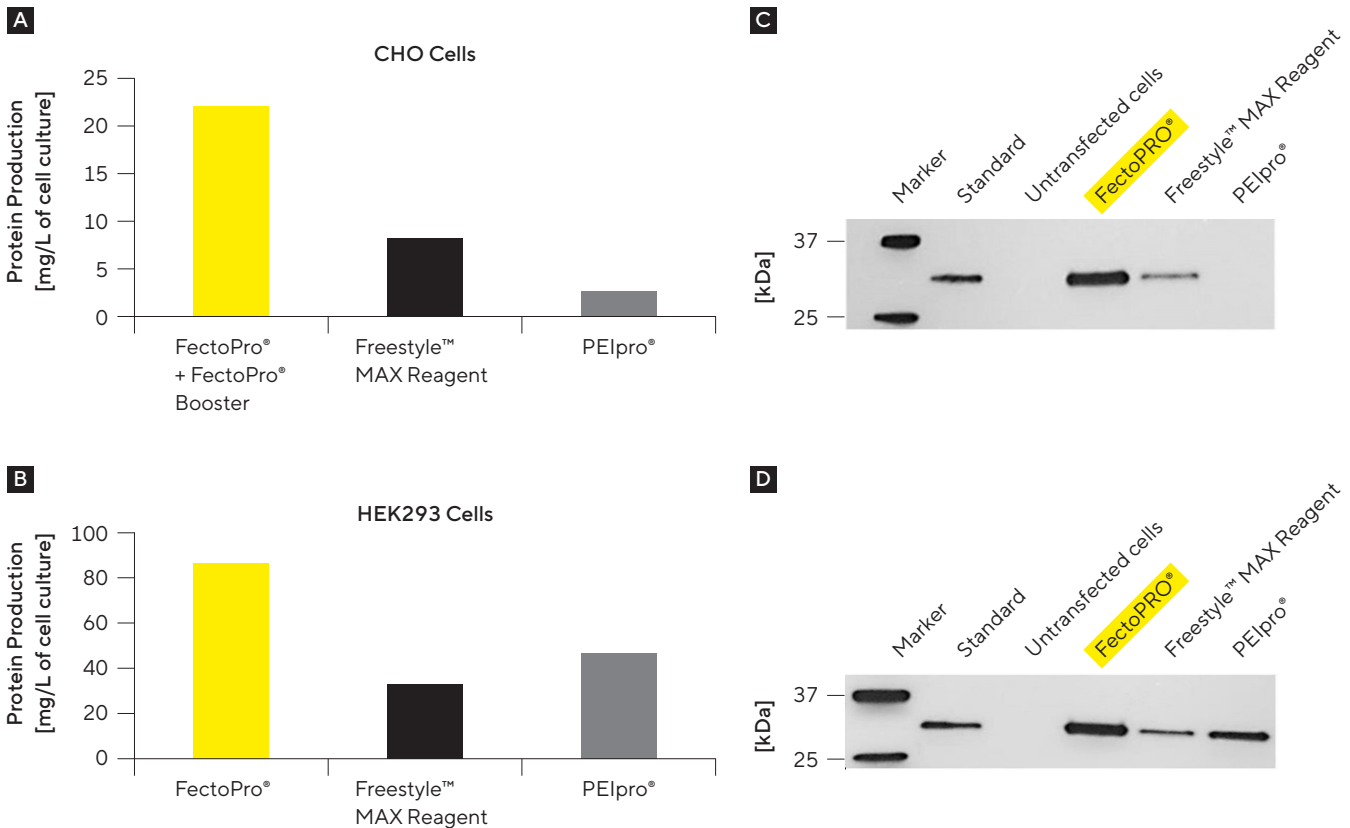
- Protein and monoclonal antibody (mAb) production
- Transient protein expression
- Academic research

Relevant Process Steps

- Upstream process
- Transfection

FectoPRO® gives superior transfection efficiency leading to higher protein production yields in both suspension CHO and HEK293 cells compared to other transfection reagents such as linear PEI.

Figure 1: FectoPRO® Gives Superior Production Yields in Suspension CHO and HEK293 Cells



Note. FreeStyle™ CHO-S and HEK293F cells were seeded at 1×10^6 cells/mL in 30 ml of FreeStyle™ CHO or 293 Expression Medium and transfected with FectoPRO® + FectoPRO® Booster ($0.4 \mu\text{g DNA/mL}$), FreeStyle™ MAX Reagent ($1.25 \mu\text{g DNA/mL}$) or PEIpro® ($1 \mu\text{g DNA/mL}$) following the recommended protocols. IgG3-Fc (34 kDa) production was assayed 72 hours (CHO cells) or 120 hours (HEK293 cells) after transfection by (A and B) Protein G affinity quantification (HPLC) and (C and D) Western blot. Standard corresponds to purified IgG3-Fc (34 kDa).

Technical Specifications

| | |
|--------------------------------|--|
| Molecule delivered | DNA |
| Cell types | Suspension CHO, HEK293, and derivatives |
| Number of transfections | 1 mL of FectoPRO® is sufficient to transfect 800 mL to more than 1 L of cell culture |
| Storage | 5 °C ± 3 |
| Provided with | FectoPRO® Booster |

Ordering Information

| Item | Description | Package Volume Quantity Size | Order Number |
|------------------|---|---|---------------------|
| FectoPRO® 1 mL | Transfection reagent for protein and mAb production | 1 mL bottle | 101000014 |
| FectoPRO® 10 mL | Transfection reagent for protein and mAb production | 10 mL bottle | 101000015 |
| FectoPRO® 100 mL | Transfection reagent for protein and mAb production | 100 mL bottle | 101000016 |

Germany

Sartorius Stedim Biotech GmbH
August-Spindler-Strasse 11
37079 Goettingen
Phone +49 551 308 0

USA

Sartorius Stedim North America Inc.
565 Johnson Avenue
Bohemia, NY 11716
Toll-Free +1 800 368 7178

France

Polyplus®—Now part of Sartorius
75 Rue Marguerite Perey
67400 Illkirch
Phone +33 390 406 180



For more information, visit

sartorius.com/transfection-reagents-plasmids