

# 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

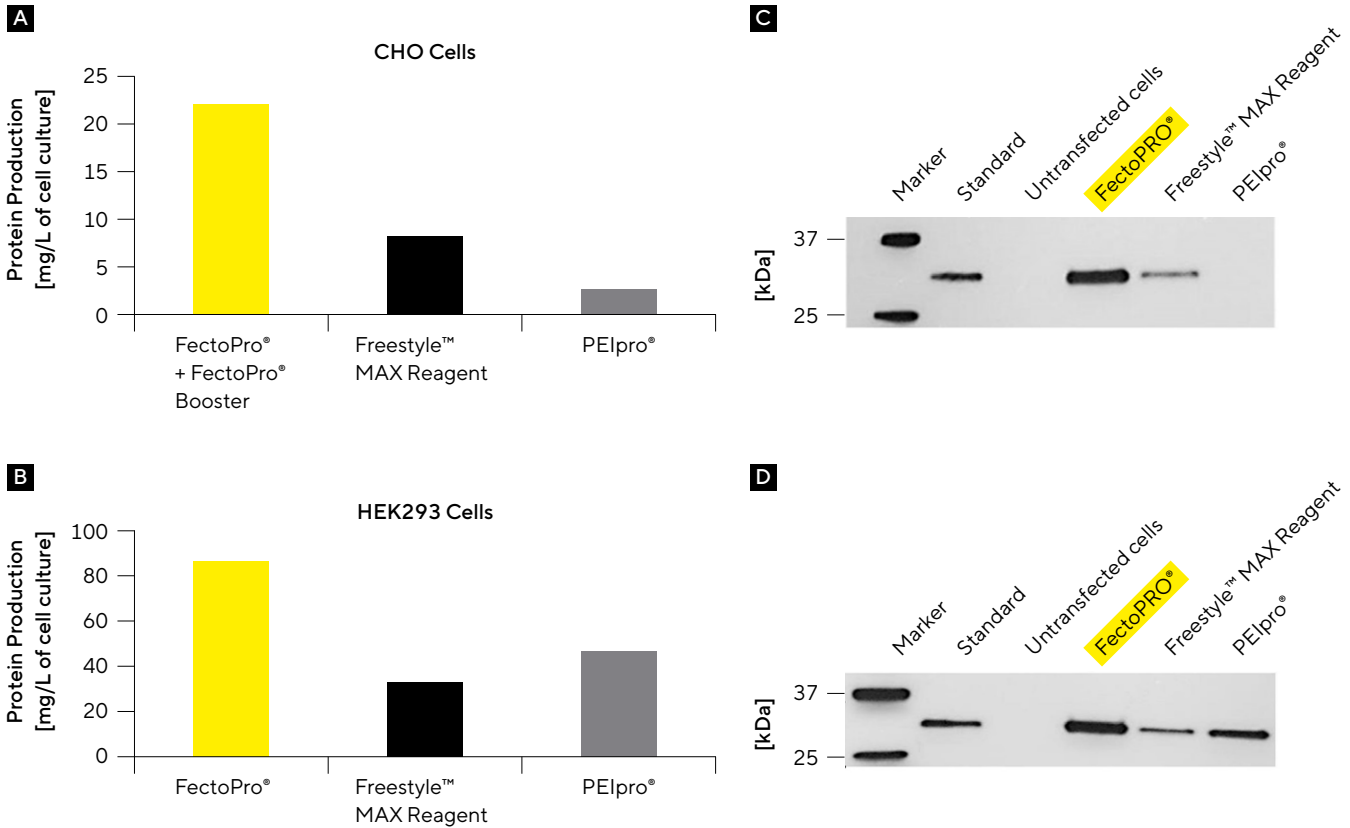
- Protein and monoclonal antibody (mAb) production
- Transient protein expression
- Academic research

## 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  $1 \times 10^6$  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

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

# Ordering Information

<b>Item</b>	<b>Description</b>	<b>Package   Volume   Quantity   Size</b>	<b>Order Number</b>
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	101000007
FectoPRO® 100 mL	Transfection reagent for protein and mAb production	100 mL bottle	101000019

**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](https://sartorius.com/transfection-reagents-plasmids)