

# SARTORIUS

## HEK293 Media Sample Kit

CFV3FB4000



# Cell Culture Media and Applications

Growth and transfection media suitable for various HEK cell lines.

Designed as all-in-one solutions, from thawing, growth, transfection and production; no need for usage of several media. On the next page you can find further information to decide which medium best suits your needs.

Medium (1 L Bottle)	Growth Factor	Characteristics/Comparison	Order Number*
HEKGM	with	<ul style="list-style-type: none"><li>Stable/Robust</li><li>High nutrient level</li></ul>	851-0001
HEKViP NX	without	<ul style="list-style-type: none"><li>Highest nutrient level</li></ul>	892-0001
HEKViP NB	without	<ul style="list-style-type: none"><li>Medium nutrient level</li><li>Leanest formulation (lowest number of components)</li></ul>	891-0001
HEKTF	with	<ul style="list-style-type: none"><li>Broad application</li></ul>	861-0001

\* These order numbers are for ordering individual media bottles

## Characteristics

- 100% chemically-defined
- Little/No adaptation from other media
- Stable growth in seed train culture
- High and robust performance
- For suspension cultures
- CFV3FB4000 for research use only
- Available for further manufacturing use upon request

## 100% free of

- Serum and animal components
- Hydrolysates and L-glutamine

## Available as

- Liquid and powder formulations
- Research and production scale

## Missing information?

Please send an email to [support@xell.de](mailto:support@xell.de)

## Download associated documents here

[www.xell.ag/shop/document-search](http://www.xell.ag/shop/document-search)

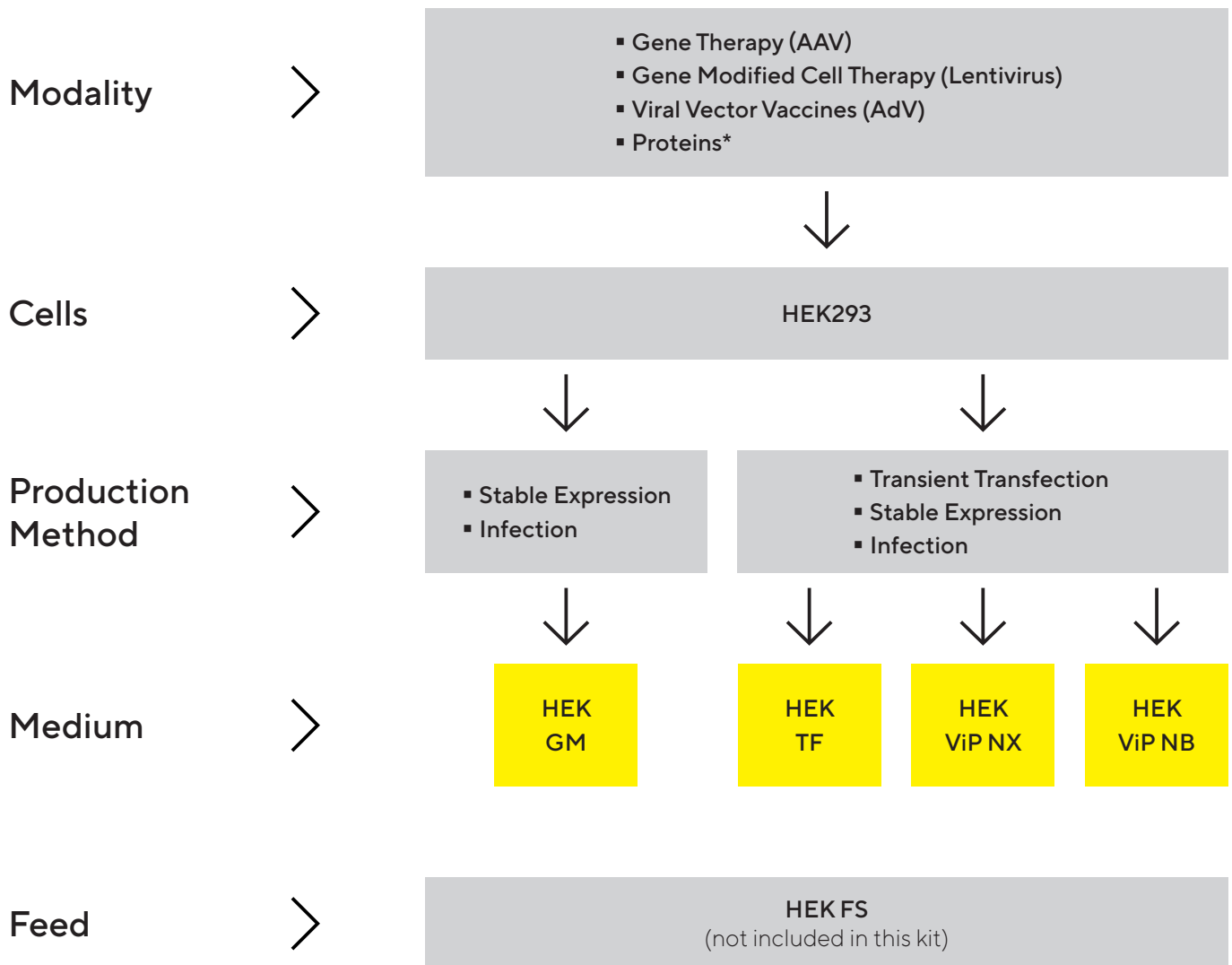
## For more information, please visit

[www.xell.ag/shop/culture-media-and-feeds](http://www.xell.ag/shop/culture-media-and-feeds)

Scan to sign up for your free  
HEK293 Media Sample Kit



# Simplifying Your Decision



\* HEK293 media can be used for mAb or protein production

## **Germany**

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

Sartorius Xell GmbH  
Waldweg 21  
33758 Schloss Holte-Stukenbrock  
Phone +49 521 96989 200

## **USA**

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

 **For more information, visit**

[www.sartorius.com](http://www.sartorius.com)

[www.xell.de](http://www.xell.de) | [info@xell.de](mailto:info@xell.de)

Specifications subject to change without notice.

© 2022 Copyright Sartorius Stedim Biotech GmbH, August-Spindler-Strasse 11, 37079 Goettingen, Germany

Status March 2022