SARTUR

HEK293 Media Sample Kit

CFV3FB4000 | Chemically Defined Cell Culture Media Designed for HEK293 Cell Lines



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	Stable/RobustHigh nutrient level	851-0001
HEK ViP NX	without	 Highest nutrient level 	892-0001
HEK ViP NB	without	 Medium nutrient level Leanest formulation (lowest number of components) 	891-0001
HEKTF	with	 Broad application 	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

100% free of

- Serum and animal components
- Hydrolysates and L-glutamine

Available as

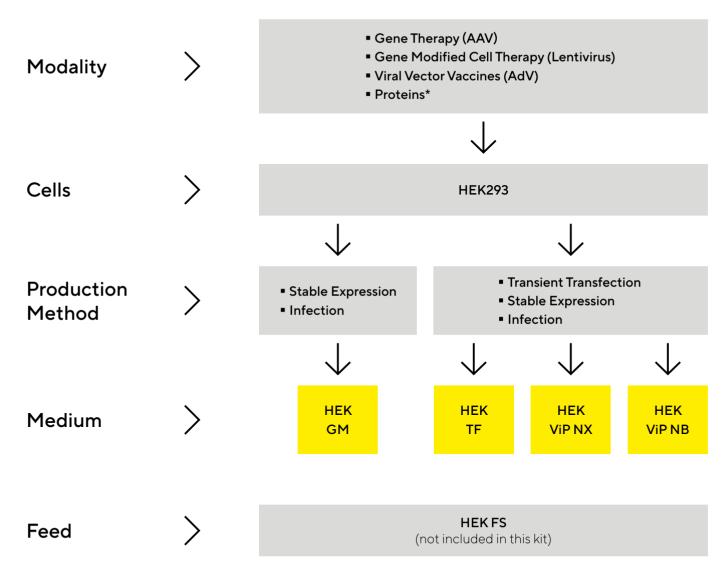
- Liquid and powder formulations
- Research and production scale

Scan to sign up for your free HEK293 Media Sample Kit



For more information, please visit www.sartorius.com/cell-culture-media

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

For more information, visit www.sartorius.com

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

Specifications subject to change without notice.

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