



HEK media and feed portfolio

High performance batch, fed-batch and perfusion

CHARACTERISTICS

- 100% chemically-defined
- little/ no adaptation from other serum-free media
- stable growth at high viability in seed train culture
- high and robust performance

100% FREE OF

- serum and animal components
- hydrolysates and L-glutamine

AVAILABLE AS

- liquid and powder formulations
- research and production scale

MEDIA AND APPLICATIONS

Growth media suitable for various HEK cell lines. Designed as all-in-one solutions, from thawing, adaptation, growth and protein production- no need for usage of several media.

Medium	GF ¹	HT ²	Product information	Order no.
HEK GM	+	+	robust and effective base medium	851
HEK ViP NX	-	+	for growth, transfection/infection and production	892
HEK ViP NB	-	+	for growth, transfection/infection and production	891
HEK TF	+	+	for growth, transfection/infection and production	861

¹ GF: growth factor, ² HT: hypoxanthine/thymidine

FEEDS AND APPLICATIONS

Supplements for prolonged production of biologics in fed-batch processes. Contain highly concentrated nutrients to increase productivity of cultured cells.

HEK FS	with 40 g/L D-glucose; for all HEK cells	871
Basic Feed	with 20 g/L D-glucose; also for hybridoma	1092

📌 You could not find, what you were looking for?
Feel free to express wishes to info@xell.de

📌 For more information please visit
www.xell.ag/shop/cultur-media-and-feeds/

CONTACT US

Xell AG
Alte Verler Strasse 1
33689 Bielefeld
Germany

Research & Development
+49 (0)521 96989-200
+49 (0)521 96989-201 fax
info@xell.de
www.xell.de



Xell AG
Waldweg 21
33758 Schloss
Holte-Stukenbrock
Germany

Production & Analytics
+49 (0)5207 9597-200
+49 (0)5207 9597-201 fax
info@xell.de
www.xell.de