

# BD Gentest High Viability CryoHepatocyte Recovery Kit

## Features

- Improves cryopreserved hepatocyte cell viability, yield, and plating efficiency
- Saves researchers time with an optimized single centrifugation recovery step versus traditional two-step processes
- Provides recovery medium for two purifications; up to five vials of cryopreserved hepatocytes can be processed per purification
- Supports human plateable cryopreserved hepatocyte recovery and plating for induction assays or other assays requiring plated cells such as metabolism, transporter, and toxicity studies

## Viability Improvement

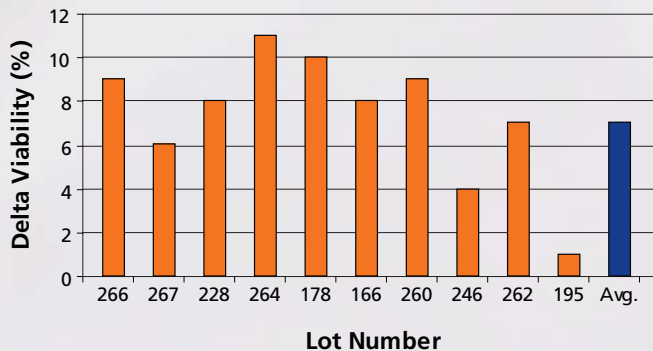


Figure 1. Viability Improvement versus Conventional Recovery Medium. Post-thaw viability increased 7% on average for a representative 10 BD Gentest inducible cryopreserved hepatocyte lots when the BD Gentest high viability cryohepatocyte recovery medium was compared to conventional recovery medium.

## Recovery Improvement

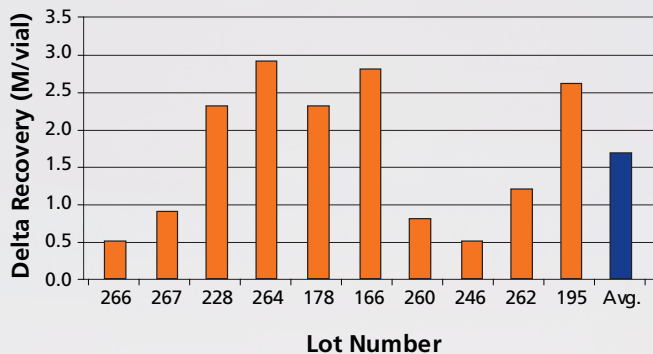


Figure 2. Recovery Improvement versus Conventional Recovery Medium. Post-thaw recovery increased 1.7 million cells/vial on average (25% increase) for a representative 10 BD Gentest inducible cryopreserved hepatocyte lots when the BD Gentest high viability cryohepatocyte recovery medium was compared to conventional recovery medium.

The BD Gentest™ high viability cryohepatocyte recovery kit provides a high-viability, high-yield recovery medium optimized for a single centrifugation step as well as an optimized plating medium for high plating efficiency.

## Increases post-thaw viability and yield

Tests on 25 lots of BD Gentest inducible-qualified cryopreserved hepatocytes showed the BD Gentest high viability cryohepatocyte recovery kit achieved  $\geq 80\%$  post-thaw viability for all lots. Recovery of viable cells also increased by an average of 1.5 million cells/vial (24% increase) when compared to other recovery and thawing medium.

## Single recovery centrifugation step saves time versus traditional two-step processes

The non-cytotoxic nature of the recovery medium allows researchers to use a one-step centrifugation step and move directly to plating hepatocytes. The need for a second wash and centrifugation step is eliminated.

## Convenient format

The BD Gentest high viability cryohepatocyte recovery kit provides two tubes of post-thaw recovery medium and two tubes of plating medium for high throughput screening or pooled product applications.

Visit [bdbiosciences.com/admetox](http://bdbiosciences.com/admetox) for more information.



# BD Gentest High Viability CryoHepatocyte Recovery Kit

## Plating Efficiency and Cell Health

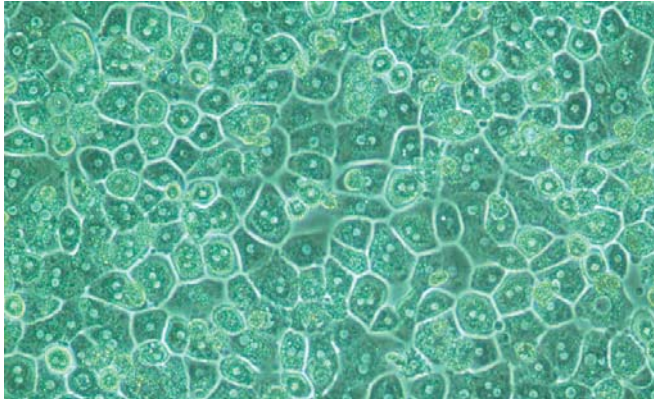


Figure 3. Plating Efficiency versus Conventional Recovery Medium.

The BD Gentest high viability cryohepatocyte recovery kit maximizes plating efficiency when compared to conventional plating medium. BD Gentest inducible cryohepatocytes, Lot 281, demonstrated 95% cell confluency after 24-hour post-thaw. Confluency determined by visual inspection. Magnification 200X.

## Related Products

- BD Gentest™ inducible-, platable metabolism-, and platable transporter-qualified human cryohepatocytes
- BD™ hepatocyte culture media and supplements
- BD BioCoat™ Hepatocyte Differentiation Environment
- BD extracellular matrices including BD Matrigel™ matrix
- BD BioCoat collagen-coated plates, dishes and flasks
- BD Falcon™ cell cultureware
- BD Falcon pipets
- BD Gentest ADME/Tox chemicals including substrates, inhibitors, and metabolites
- BD Gentest<sup>SM</sup> contract research services including induction and metabolic stability assays using hepatocytes

## Multiple Vial Post-thaw Viability and Recovery

Up to 5 vials of hepatocytes can be processed with one tube of medium without loss of viability and recovery.

No. vials of cryopreserved hepatocytes used per tube of recovery medium	Lot 251		Lot 247	
	Viability (%)	Recovery (M/Vial)	Viability (%)	Recovery (M/Vial)
1	84.2	6.7	80.8	11.8
2	84.5	7.4	79.0	12.3
3	83.9	6.8	81.7	12.5
4	81.3	6.6	76.7	12.3
5	81.3	6.8	80.3	13.3
6	78.4	6.5	75.5	12.4

Post-thaw viability and recovery using up to 5 vials of cryohepatocytes per tube of recovery medium demonstrates the equivalent viability and recovery as individually thawed tubes of cryohepatocytes. Greater than 5 vials of cryopreserved hepatocytes per tube of recovery medium demonstrated a 5% drop in viability.

## Order Information

BD Gentest High Viability CryoHepatocyte Recovery Kit

Description	Qty.	Cat.No.
<b>Platable human cryopreserved hepatocyte recovery and plating medium, kit contains:</b> • 2X tubes (45 mL each) recovery media • 2X tubes (45 mL each) plating media	1 kit	454534

BD Gentest High Viability Recovery Medium

Description	Qty.	Cat.No.
<b>BD Gentest High Viability Recovery Medium</b>	1 tube @ 45 mL	454560

BD Gentest CryoHepatocyte Plating Medium

Description	Qty.	Cat.No.
<b>BD Gentest CryoHepatocyte Plating Medium</b>	1 tube @ 45 mL	454561

BD Gentest Human Plateable Cryopreserved Hepatocytes

Description	Qty.	Cat.No.
BD Gentest Plateable Transporter-Qualified Human CryoHepatocytes	≥5 million cells	454541
BD Gentest Plateable Metabolism-Qualified Human CryoHepatocytes	≥5 million cells	454543
BD Gentest Inducible-Qualified Human CryoHepatocytes	2-5 million cells	454550
	≥5 million cells	454551

\*Please note that all orders must be lot specific. Each lot number corresponds with a specific donor. For the most recent listing of available donor information, please visit [bdbiosciences.com/support/resources/adme/index.jsp#admetools](http://bdbiosciences.com/support/resources/adme/index.jsp#admetools)

U.S. Orders  
tel: 877.232.8995  
e-mail:  
BDBCcustomerService@bd.com

Purchase Orders should be made out to:  
BD Biosciences  
2350 Qume Drive  
San Jose, CA 95131

## Important Ordering Information

**HepatoLink** is the BD Biosciences immediate response system providing researchers with instant notification when BD Gentest hepatocyte products are available. For your convenience, select a preferred contact method—e-mail, phone, and/or SMS text alerts. Upon registration, you will be sent donor characterization data when a new lot becomes available. Register online at

[www.bdbiosciences.com/admetox](http://www.bdbiosciences.com/admetox)

Place your order directly through BD's online system. If you have any questions regarding BD's program or your participation, please contact us at 978.901.7260 or e-mail [hepatocytes@bd.com](mailto:hepatocytes@bd.com).



**BD Biosciences**  
Two Oak Park  
Bedford, MA 01730  
tel: 877.232.8995  
fax: 800.325.9637  
[bdbiosciences.com](http://bdbiosciences.com)