



*XenoTech Now Offers Over 1,100 Cell Lines to Bridge the Gap Between Research and Discovery*

**JCRB (Japanese Collection of Research Biosources)**

XenoTech has teamed up with the National Institute of Biomedical Innovation to bring you access to roughly 1,100 cell lines, helping you to bridge the gap between research and discovery. XenoTech is the exclusive US provider of JCRB's comprehensive catalog of cell lines.

The Japanese Collection of Research Biosources was founded by the ministry of Health, Labor and Welfare in 1984. They provide one of the most comprehensive cell banks in the world to users across the globe.

Types of cell lines offered include:

- General cells
- Luciferase-expressing cancer cells
- Mouse homozygous mutant ES cells
- Immortalized Mesenchymal stem cells
- Genetically-modified cells

To view the complete list of cell lines offered by XenoTech, visit [www.xenotech.com/products/jcrb-cell-bank](http://www.xenotech.com/products/jcrb-cell-bank)

*Ordering Process Made Simple*

XenoTech is committed to making the ordering process for JCRB cell lines seamless. All aspects of your order will be handled by XenoTech including invoicing and importation. Simply place your order for the desired cell lines and XenoTech takes care of the rest.

XenoTech imports JCRB cell lines on a bi-weekly basis. Upon placement of your order, XenoTech will notify you of your estimated product delivery date and invoice you at the time of shipment.

Contact a customer service representative at **913.438.7450** or email us at [info@xenotechllc.com](mailto:info@xenotechllc.com)

Cell Line Pricing		
Cell Type	Non-Profit Organization	Profit Organization
General cells	\$740	\$790
Luciferase-exp cancer cells	\$800	\$850
Mouse homozygous mutant ES cells	\$800	\$850
Immortalized Mesenchymal stem cells	\$740	\$790
Genetically-modified cells	\$800	\$850

All pricing is inclusive of shipping & handling charges

*"We received the cell lines in great condition. I can't tell you how wonderful and smooth it was to receive materials from XenoTech! We are so happy that you are distributing the JCRB lines in the US."*

**- Associate Director, Life Science Company**