

JCRB Cell Line Ordering Stepwise Procedure

From XenoTech Website:

1. Click on “Place your JCRB Cell Line Order” at bottom of page.

[Place Your JCRB Cell Line Order](#)

2. Review the “[How to order](#)” page on JCRB Website. Confirm that Pricing Information, Distributorship information, Cancellation Policy, and Shipping schedule are all acceptable within your project needs.
3. Click “Cell search top”, located in the top left hand corner of the side navigation panel.
4. If you know which cell line you are trying to order, use the “Quick Search” or “Advanced Search” applications to locate the cell line of interest. If you want to browse for cell lines, use the “[Alphabetical List](#)” or “[Animal/Tissue List](#)”. XenoTech has also created a [sortable list in excel](#) to ease browsing.
5. Identify the cell line(s) you wish to order and click “Order” to the left of the description to add the respective item to your shopping cart. To add more items to your cart, click “Additional order” and proceed to Step 6 of this document. To proceed to Checkout, select “Order form(SSL)” and proceed to Step 8 of this document. To remove an item from your order, click “Cancel” to the right-hand side of the description.
6. Add more items: You have selected “Additional order” and are now viewing what you currently have in your cart. Select the green plus sign next to “Search Box” to the upper left-hand side of your item descriptions for a drop down version of the “Quick search” application you used in Step 4. At any time, you may select “View Cart” on the upper right-hand side of the item descriptions to view your selected items and proceed to checkout.
7. Repeat Steps 4-6 as necessary.
8. You are ready for checkout and have selected “Order form(SSL)”. Please review the ordering process, cancellation policy and the agreement. If you agree to terms, select “Order form for outside Japan(SSL)”.
9. Confirm your items are correct. If you have an account, select “User login” and enter your credentials; if you do not have an account, select “User registration”.
10. Follow onscreen instructions until the order is placed. Note: Payment is NOT required at time of order placement. XenoTech will follow-up to collect payment.

After Order Submission:

You will be prompted to sign “Form A”. Please sign and send back to JCRB promptly. This will expedite the shipping process.

If you have any questions, please contact a XenoTech representative at info@xenotechllc.com.