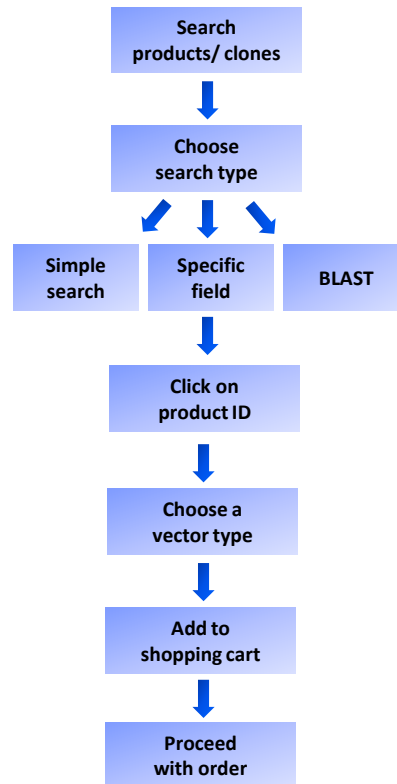


**Follow these simple steps to find your favorite clones from GeneCopoeia.**

1. Go to “Search products/clones” webpage.
2. Choose any of three options provided to start the search
  - Use “Simple Search.” Type any keyword.  
(Description, Nucleotide accession, Entrez gene ID, Symbols/Aliases or Catalog/product ID)
  - Search by a particular field
  - Search by BLAST
3. The search engine gives options based on the search query entered. Click on the column “Product ID.”
4. The webpage with all the information available for that gene with the chosen product ID comes up.
  - Choose among 80 different ready-for-expression vector types including mammalian, bacterial, lentiviral, insect and yeast.
  - Gateway shuttle system and ORFeome Collaboration clones are also available.
  - Related products used with the clone of interest is also listed for quick access.
5. To learn more about the specific gene or to view the gene annotation page, click on “Product ID” again.
6. To view price and proceed with the order, select the appropriate expression vector type.



## Search Clones