

FREQUENTLY ASKED QUESTIONS FOR VECTOR NTI SOFTWARE

1. Are all of the Vector NTI Advance, Express and Express Designer software versions going to be discontinued at the same time? Yes, all three software versions (including all available catalog numbers) are going to be discontinued at the same time.
2. What is the last date for discontinuation of Vector NTI software sales? The last day for the sales of Vector NTI desktop software licenses or support contracts will be December 31, 2019. Until this day, existing and new customers will be able to purchase or renew their licenses. Customers will be able to obtain support until December 31, 2020.
3. Will the existing licenses and support contracts also stop with the discontinuation of Vector NTI? Yes, all existing NTI licenses types and support contracts will stop with the discontinuation of Vector NTI for all the customers on December 31, 2020.
4. What is the last date for discontinuation of the technical support? The last day to receive support will be December 31, 2020. After this date, users will no longer have access to technical support, replacement keys or reinstallation requests.
5. How long will the support be provided to existing customers after the official discontinuation (i.e. after December 31, 2020)? After December 31, 2020, users will no longer have access to any technical support, replacement keys or reinstallation requests.
6. What are the guidelines for customers who have purchased licenses up to the December 31, 2019? Customers will be able to use support services provided by the support contract till December 31, 2020. After December 31, 2020, access to replacement keys and support services no longer will be available. Current users are encouraged to export data from Vector NTI software using launched Vector NTI Data Export Tool and migrate to alternative software solutions (see Question 11 below) before the last day of support. No return of licenses and support process will be maintained if customer decides to return the license(s) or support contract(s) after discontinuation of sale of licenses and support contracts (December 31, 2019).
7. What are the guidelines for customers after the end of Vector NTI software support (December 31, 2020)? Vector NTI licenses and support contracts will no longer exist in our catalog. Thermo Fisher Scientific will no longer be able to assist customers with Vector NTI software - technical, license renewal or support issue. Free downloads of the software will no longer be available after December 31, 2020.
8. Will Vector NTI Data Export tool be available to export of the data (databases) before December 31st, 2020? Yes, Vector NTI data Export Tool will be available to export data (databases) before December 31, 2020.
9. What are the guidelines for customers regarding utilization of Vector NTI Export Data tool after December 31, 2020? If customers choose not to transition or export their data before December 31st, 2020, they do so at their own risk. No support regarding data export will be provided after December 31, 2020.
10. What option will exist for customers who own non-expiring and dynamic license as well as signed advanced support? The same answer as to question No.6.

11. What are the recommendations for alternate software that can be used after exporting files from Vector NTI? Is the software dependent on Vector NTI version used previously?
Vector NTI software supports the export of data into standard file formats like GenBank, GenPept, Fasta, Rebase, CSV and Text. The Data Export Tool can export the entire Vector NTI database into GenBank, GenPept, Rebase and CSV format. Any software that supports the above-mentioned file formats can be used as an alternative to Vector NTI. However, customers should perform their own analysis to determine which software will best suit their needs.
12. What software will be used to validate the data exported out from Vector NTI?
UGene, Geneous, Benchling, SnapGene
13. Can I export my ContigExpress and alignment projects through Vector NTI Data Export Tool? No. Unsupported data types which are not maintained by Data Export Tool are: Citations, Blast Analysis, Analysis Results, Contig Projects, Alignment Projects, Cloning Projects, Gene Synthesis Projects. the tool will not be exporting project specific information:
 - Name of the project.
 - Name and sequence information of fragments used in the project. e.g. the fragments used to create a cloning construct or a contig or an aligned sequence.
 - Input values used to execute the algorithm for the project. For e.g. the input values used to run 'cap3' algorithm in Contig Express or the input values used to run 'clustalw' algorithm in AlignX.
14. Will subfolder structure be maintained in the export process? Yes. It will be maintained as folders.
15. What other software are available that can read the parts circuits and devices SBOL files?
Data Export tool can only export out Part/Devices/Circuits data in XML format. It does not support SBOL format.
16. Does the new export tool maintain parent child relationships that are present in Vector NTI Advanced?
The parent-child relationship is maintained in the database of that particular software. Different software would maintain it in different formats and when the molecules is loaded, it displays this information. The Data Export tool will create multi record GenBank files, which include descendant and its parent(s) information. The tool will also create a .csv file which contains information like child sequence positions in the parent.
17. Will the fragments used in alignment or Contig assembly projects be exported? Will these have to be re-aligned or re-assembled in other compatible software after the export?
Yes, saved fragments will be exported as individual molecules. Contig/AlignX project settings would be lost. Exported molecules have to be re-aligned/re-assembled using selected software. This is because the project settings of Vector NTI are in a proprietary format. This format may not work with other software.
18. What are the formats for Enzymes and Oligos molecules compatible with other software to work with?
Enzymes are exported out in a Rebase format described here:
<http://rebase.neb.com/rebase/rebase.f7.html>
Oligos are exported out in csv format.