

USER INSTRUCTIONS

GenExact™ ssDNA and GenWand™ dsDNA

Storage Conditions

- Keep the ssDNA/dsDNA tube(s) tightly sealed at -20 °C for long-term storage and avoid repeated freeze-thaw cycles.
- Avoid repeated freeze-thaw cycles after dissolving. If necessary, divide the stock solution into small aliquots.
- Stock solutions at high concentration (> 100ng/μL) can be stored at -20 °C for 6 months.

Tips for Re-suspension

- Centrifuge tube(s) at 12000 rpm for 3 minutes before opening to ensure samples are at the bottom of the tube(s).
- Use nuclease-free water or other desired buffer to dissolve the single-stranded DNA into the stock solution or working concentration. When dealing with mammalian cell lines, embryos, or primary cells, the buffer should be free of endotoxin and microorganisms.
- To open potential secondary structures, it is recommended to heat the ssDNA/dsDNA sample at 70°C for 5 minutes and store on ice before use.
- When using a working concentration of 1 μg/μl or higher, we recommend to pipet the solution multiply times to ensure all the ssDNA/dsDNA is dissolved evenly. Then, another round of centrifugation is recommended to ensure the samples are at the bottom of the tube(s).

Files Instructions

- Please view sequencing file to verify if the sequence is 100% accurate. If any mismatches are found, contact our customer service immediately. Note: If you have opted into GenScript's Green Program for paperless state-ments, then you will not receive printouts of contigs; you may login to your online account to download them.
- To view sequence chromatograms, login to your account via <https://www.genscript.com/customer/login> and click "My Orders" on the left navigation, then click Order ID & scroll down to view the chromatogram data and construct map. If you forgot your password, click "Forget your password?" on your login page and follow the instructions on the screen.
- Sequencing chromatogram (trace) data can be viewed by using Chromas (PC) or Editview (Mac). Sequence alignment files can be opened by DNASTAR® SeqMan.
- Construct maps (for gene synthesis, mutagenesis, and cloning service) can be opened by Adobe PDF Reader.