

GIBCOBRLLIFE  TECHNOLOGIES™**SEQUENCER_x Enhancer
Solution F****Cat. No. 12244-018**
reactions**Size: 2 x 100****Store at -20°C**Description:

SEQUENCER_x Enhancer Solution F is a novel reaction co-solvent (patent pending) which facilitate sequencing through problematic GC-rich regions, inverted repeats and/or hairpins. The component is added to the reaction mixture, and the mixture is incubated in a thermocycler as usual. Early terminations and/or low signal strength are typically improved within the troublesome region, permitting additional sequence data to be obtained. SEQUENCER_x Enhancer Solution F is fully compatible with BigDye terminator, dRhodamine dye terminator and rhodamine dye terminator chemistries and all sequencing platforms. Sufficient reagents are supplied based on a final 1X reaction volume of 20 µl.

<u>Component</u>	<u>Amount</u>
20X SEQUENCER _x Solution F	2 x 100 µl

Quality Control:

SEQUENCER_x Enhancer Solution F is evaluated with a GC rich (80% GC) template in a fluorescent sequencing assay using dRhodamine dye terminator chemistry.

Doc. Rev. 102700

This product is distributed for laboratory research use only. CAUTION: Not for diagnostic use. The safety and efficacy of this product in diagnostic or other clinical uses has not been established.

For technical questions about this product, call the Life Technologies TECH-LINE™ (800) 828-6686.

General Recommendations and Guidelines:

- Calculate the amount of Enhancer required based on sequencing reaction volume. The recommended final concentration in a sequencing reaction is 1X. For example, to a 20- μ l BigDye Terminator reaction, add 1 μ l of 20X SEQUENCER_x Enhancer Solution F. However, the optimal effective concentration may vary depending on the exact nature of the difficult region within the DNA template. A final concentration of 0.5X to 2X can be used.
- In place of an equal volume of the water addition, add the appropriate volume of Enhancer to either a master mix or an individual sequencing reaction. Mix well after Enhancer addition.
- There is no need to make any other changes to the composition of a standard sequencing reaction or to the thermocycling conditions.
- Resolution of problem templates may be negatively impacted if sequencing reactions are scaled down below 0.25X to 0.5X.
- For optimal results, do not combine more than one SEQUENCER_x Enhancer Solution in a single sequencing reaction.
- These reagents are not designed to solve sequencing problems associated with long (>30) poly A structures.

Limited Label License No. 12: The purchase of this product conveys to the buyer the non-transferable right to use the product and components of the product in research conducted by the buyer. The buyer cannot sell or otherwise transfer this product or its components to a third party and in particular, no rights are conveyed to the buyer to use the product or its components for commercial purposes. Commercial purposes means any activity for which a party receives consideration and may include, but is not limited to, (1) use of the product or its components in manufacturing, (2) use of the product or its components to provide a service, information or data, (3) use of the product or its components for diagnostics purposes, or (4) resell the product or its components, whether or not such product or its components are resold for use in research. If purchaser is not willing to accept the limitations of this limited use statement, Life Technologies, Inc. is willing to accept return of the product with a full refund. For information on purchasing a license to use this product for purposes other than research, contact the Director of Licensing, 9800 Medical Center Drive, Rockville, MD 20850. Phone (301) 610-8000. Fax (301) 610-8383.