

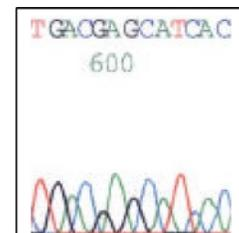
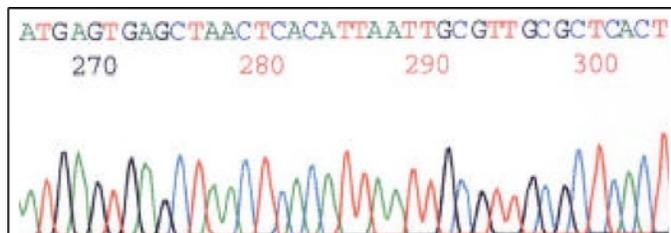
HalfBD DNA Sequencing Reagents

HalfBD® is a specially formulated DNA sequencing reagent, compatible with all PE Applied Biosystem BigDye systems. Its simple approach allows you to save up to 90% of your sequencing costs, by enabling you to reduce the volume of BigDye mix used in each sequencing reaction. Its unique formula maintains your read length, accuracy and signal strength.

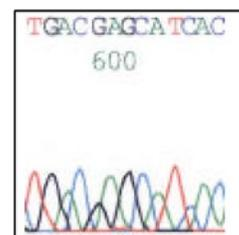
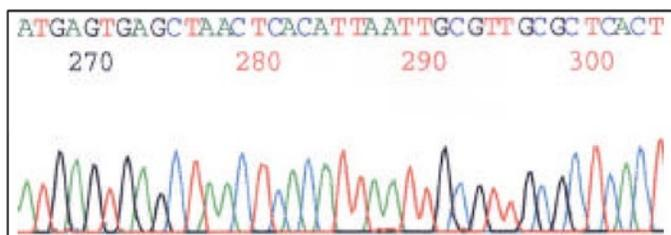
Key features

- For use on all ABI platforms including the 3700 capillary DNA sequencer.
- HalfBD has special enhancers maintain the quality of your sequence data while allowing you to use less BigDye mix per reaction.
- Simply use HalfBD to replace some of the BigDye mix in your sequencing reaction. Mix HalfBD with your BigDye mix at a ratio of 1:1 and sequence as normal.
- With good quality template DNA a ratio of 3:1 or even 7:1 can be used, while still achieving excellent quality sequence data.
- Fully compatible with BigDye version 3.
- Unique stabilizers improve the freeze/thaw characteristics of your BigDye.

3 μ L HalfBD + 1 μ L ABI BigDye mix



0 μ L HalfBD + 4 μ L ABI BigDye mix



Ordering information

Catalog number	Number of half-volume reactions*	Volume
K1102	400	0.8 mL
K1104	800	16 mL
K1110	2000	4.0 mL
K1115	10,000	20 mL

* This assumes 2 μ L HalfBD and 2 μ L BigDye per reaction.

Contact Us

Phone: +1-800-635-5577
 Web: www.moleculardevices.com
 Email: info@moldev.com

Check our website for a current listing of worldwide distributors.

Regional Offices

USA and Canada +1-800-635-5577
 Brazil +55-11-3616-6607
 China (Beijing) +86-10-6410-8669
 China (Shanghai) +86-21-3372-1088
 Germany 00800-665-32860

Japan (Osaka) +81-6-7174-8831
 Japan (Tokyo) +81-3-6362-5260
 South Korea +82-2-3471-9531
 United Kingdom +44-118-944-8000

