

ProDye™ Sanger Sequencing Standard

Instructions for Use of Products CR4604 and CR4696.

Quick Protocol

The ProDye™ Sanger Sequencing Standard can be used for a performance check on various capillary electrophoresis (CE) platforms as well as for spectral calibration on some platforms.

Using the Sequencing Standard

Materials to Be Supplied by the User

- MicroAmp® Optical 8-Tube Strip, 0.2ml (Applied Biosystems Cat.# 4316567), and Strip Septa Mat, 8-Well (Cat.# CE2308), for use on the Spectrum Compact CE System
- or**
- MicroAmp® Optical 96-Well Plate (Applied Biosystems Cat.# N8010560) and Septa (Applied Biosystems Cat.# 4412614)
- Hi-Di™ formamide (Applied Biosystems Cat.# 4401457)

The table below lists the sequencing standard product and quantity recommended for use with each capillary electrophoresis instrument.

Capillary Electrophoresis Instrument	Cat.#	Number of Tubes
Spectrum Compact CE System	CR4604	1
Applied Biosystems® 3500 Genetic Analyzer	CR4604	2
Applied Biosystems® 3500xL Genetic Analyzer	CR4696	1
Applied Biosystems® 3730 Genetic Analyzer	CR4696	1
Applied Biosystems® 3730x/ Genetic Analyzer	CR4696	1
Applied Biosystems® 3130 Genetic Analyzer	CR4604	1
Applied Biosystems® 3130x/ Genetic Analyzer	CR4604	4
	CR4696	1
Applied Biosystems SeqStudio® Genetic Analyzer	CR4604	1

1. Thaw the ProDye™ Sanger Sequencing Standard at room temperature. Vortex and centrifuge it briefly in a microcentrifuge to collect the material at the bottom of the tube.
2. To each tube of sequencing standard, add the volumes of Hi-Di™ formamide listed in the table below:

Product	Hi-Di™ formamide volume per tube
CR4604	33µl
CR4696	750µl

3. Vortex and centrifuge briefly in a microcentrifuge to collect the material at the bottom of the tube.
4. Aliquot 10µl of the standard to each desired well of the 8-well strip tube or 96-well plate.
5. Cap the wells with strip or plate septa. Centrifuge briefly in a centrifuge that can accommodate your samples to collect the material at the bottom of the wells and to remove any air bubbles from the samples.
6. Follow the instructions specific to the instrument you are using to begin your performance check or spectral calibration run.

Additional protocol information in Technical Manual #TM672, available online at: www.promega.com

For information on how to operate the Applied Biosystems® Genetic Analyzers, refer to the manufacturer's User Guide.

ProDye is a trademark of Promega Corporation.

Applied Biosystems is a registered trademark of Applied Biosystems, LLC. MicroAmp and SeqStudio are registered trademarks and Hi-Di is a trademark of Thermo Fisher Scientific, Inc.

