ProDye™ Terminator Sequencing System

Sanger 测序完美契合您的工作流程

ProDye[™] Terminator Sequencing System 是一种经过改进的荧光 Sanger 测序方法,该方法采用包含在预混液中的专有热稳定 DNA 聚合酶。ProDye[™] Terminator Sequencing System 兼容多种 DNA 模板,可与 Spectrum Compact CE System 以及 Applied Biosystems[®] 3130、3130xl、3500、3500xL、3730、3730xl 和 SeqStudio[®] Genetic Analyzers 配合使用。ProDye[™] Terminator Sequencing System 使用 dTMR、dCXR、dRSixG 和 dROneTen 终止物染料,其与 BigDye[®] Terminator V3.1 Cycle Sequencing Kit 中的染料相同。

ProDye[™] Terminator Sequencing System 与各种热循环仪兼容,其具有的简易循环方案设计用于获取最大的读取长度。测序反 应的所有组分均包含在即用型预混合形式的包装内,只需加入模板及其专用引物。包括了额外的 5X 测序缓冲液和无核酸酶水。

- 易于实现的相同方案、工作流程和校准
- 设计适用于广泛的测序应用
- 降低成本且提高生产力

与多种样品和运行条件兼容

使用"标准"或"快速"测序方案对质粒 DNA 和其他样品进行测序。



图 1. 采用 ProDye[™] Terminator Sequencing System 对 NIST 质粒 ERCC-00083 进行循环测序,采用 Spectrum Compact CE System 的"标 准"测序方案进行检测获得的图谱。采用 ProDye[™] Terminator Sequencing System 对 NIST 质粒 ERCC-00083 进行循环测序。测序反应用乙醇 / EDTA 沉淀法纯化,并在 Hi-Di[™] 甲酰胺中变性,然后在 Spectrum Compact CE System 上使用 T_Seq_36_Std 方案(图 A)和 T_Seq_36_Fast 方案(图 B)进行检测。

测序分析读取长度

将 ProDye™ Terminator Sequencing System 与 Spectrum Compact CE System 相结合,在更短的时间内获得更多的碱基。



图 2. 采用 ProDye™ Terminator Sequencing System 对 pGEM[®]-3Zf(+) 质粒对照组进行循环测序,并采用 Spectrum Compact CE System 的 T_ Seq_36_Std 方案(图 A)和 T_Seq_36_Fast 方案(图 B)进行检测获得的图谱。

您的样品、我们的测序试剂盒,适用于任何仪器

● 🗿 🎊 👐 🛶 🍳 🍳 🤋 🖕

View Report Help

将 ProDye™ Terminator Sequencing System 无缝整合到熟悉的工作流程中。



准备反应 模板 DNA 纯化 准确定量 **约 30 分钟** **热循环** 混合 2.5X 反应预混液、 模板、引物和水 准备进行热循环 **约 2-2.5 小时** **反应纯化** 去除多余的染料标记 的终止物 **5 分钟 -1.5 小时**

毛细管电泳 分离 DNA 片段 检测荧光信号 "Basecaller"软件自动 分析样品 30 分钟 -3 小时 / 每次注射 数据分析

为进一步分析生成的电 泳图(原始和经处理) 和测序信息

数小时内

订购信息

产品	规格	目录号
ProDye [™] Terminator Sequencing System	24 reactions	CR4324
	200 reactions	CR4302
	1,000 reactions	CR4310
ProDye™ 4C Matrix Standard	5 preps	CR4500
ProDye™ Sanger Sequencing Standard, 4 × 4 孔	16 wells	CR4604
ProDye™ Sanger Sequencing Standard, 2 × 96 孔	192 wells	CR4696
ProDye™ 5X Sequencing Buffer	12 ml	CR1011

仅供研究使用。不得用于诊断。

了解更多信息请登录: www.promega.com/ProDye

Promega (Beijing) Biotech Co., Ltd 普洛麦格(北京)生物技术有限公司

地址:北京市东城区北三环东路 36 号环球贸易中心 B 座 907-909 电话: 010-58256268 网址:www.promega.com 微网站:wechat.promega.com.cn 技术支持电话: 400 810 8133(手机拨打) 技术支持邮箱: chinatechserv@promega.com



欢迎关注 Promega 官方微信