

Automated RNA-seq Library Preparation on the Beckman Coulter Life Sciences Biomek i7 Hybrid Workstation Using a Shortened KAPA RNA HyperPrep Workflow

Amy Janiak¹, Amy Elias¹, Elisa Vega¹, Zach Smith², Nikita Dsouza¹, Alejandro Quiroz Zarate¹, Marsha McMakin¹, Rachel Kasinskas¹
¹Roche Diagnostics Corporation, Wilmington, MA
²Beckman Coulter Life Sciences, Indianapolis, IN

INTRODUCTION

As RNA sequencing (RNA-seq) has become more important to our understanding of disease and to the development of new pharmaceuticals, the need for faster, scalable, reproducible workflows has also grown. However, RNA-seq library preparation methods are complex and require high levels of precision, posing a challenge to the required scaling and reproducibility.

In this study, a two-fold approach was taken to overcome these challenges:

1. a new, shorter RNA-seq library prep workflow was developed using on-market kits without the need for additional reagents, and
2. this shorter workflow was automated on a liquid handler that is frequently used for NGS library preparation.

WORKFLOW MODIFICATIONS

From the original KAPA RNA HyperPrep Kit

- Shorter DNase digestion incubation (from 30 min to 15 min)
- Lower temperature for the 1st strand synthesis primer extension (20°C vs 25°C).
- Shorter 2nd strand synthesis incubation (42°C for 5 min, instead of 62°C for 30 min)
- Reduced adapter concentrations with lower input amounts, to minimize primer dimer formation
- Fewer 0.7X post-ligation cleanups (one, instead of two)

