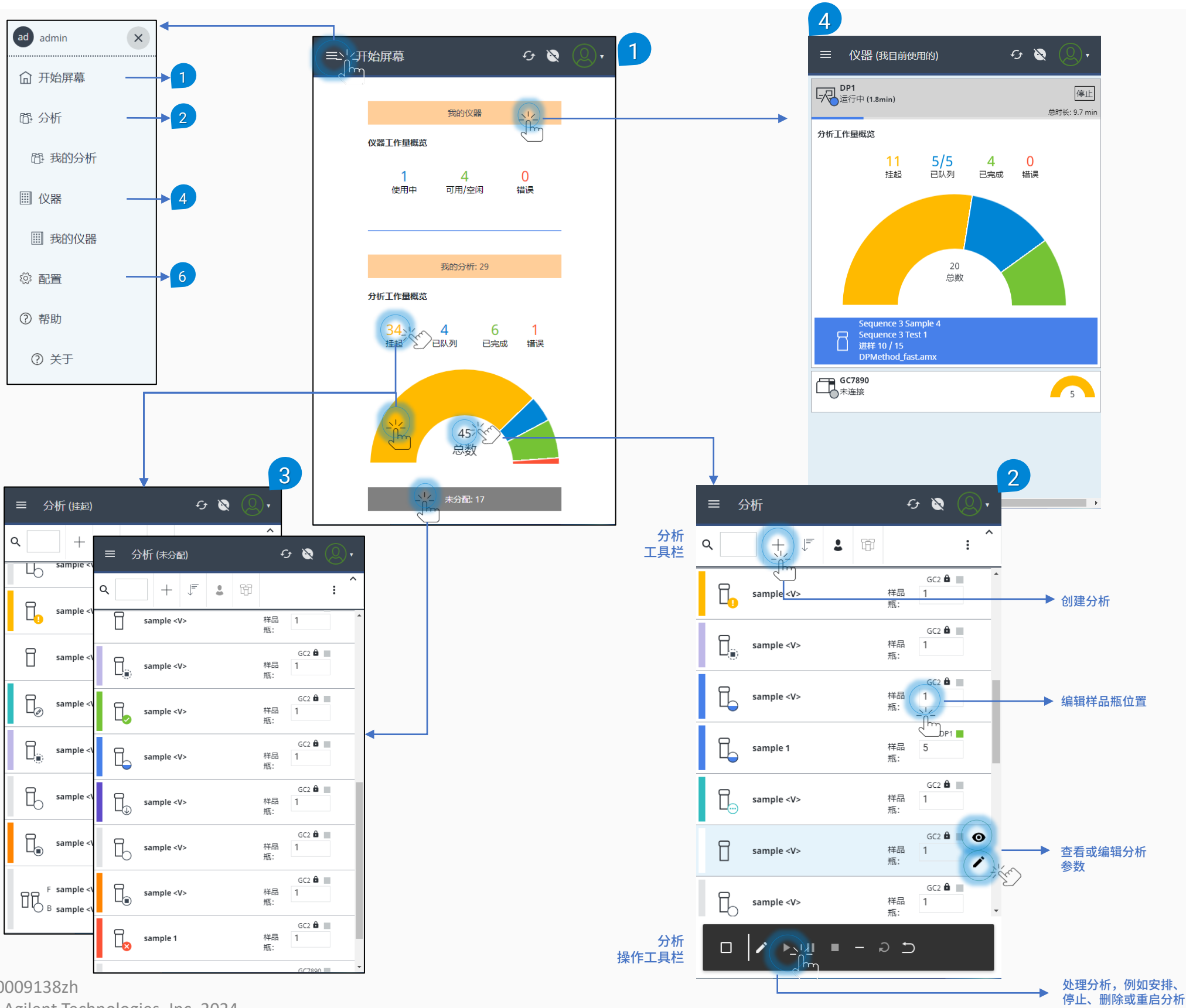


Sample Scheduler 入门指南



Sample Scheduler 入门指南



2 在分析视图中选择序列。展开窗口，直至“序列”视图按钮可用。

5 查看序列的详细信息。编辑单个序列行，或整个序列。

6 单击或点击配置按钮以查看和编辑运行队列设置。

