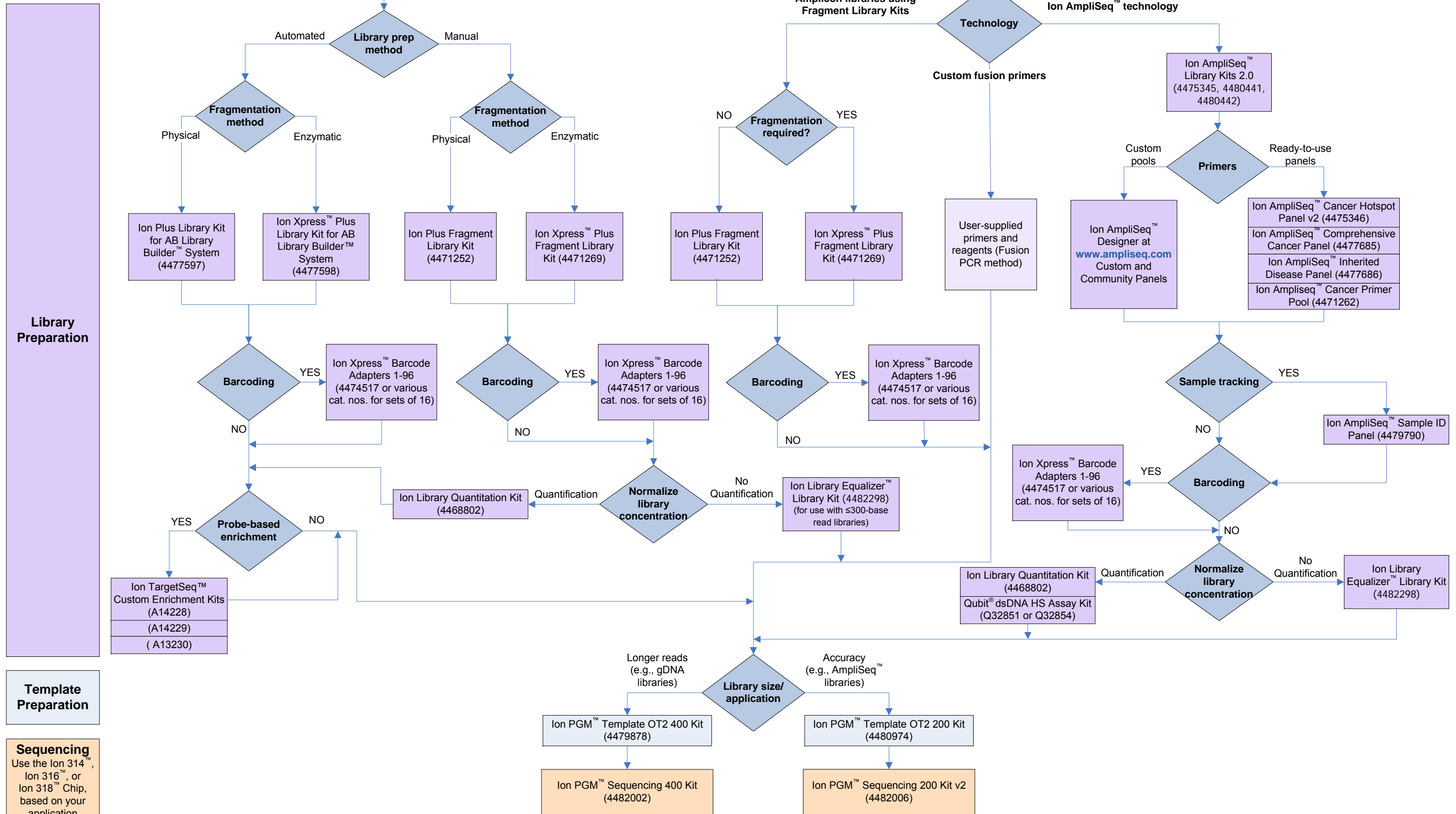


Decision Tree for DNA Sequencing on the Ion PGM™ System



Library Preparation

Template Preparation

Sequencing
Use the Ion 314™, Ion 316™, or Ion 318™ Chip, based on your application

Other applications: Visit the [Applications page](#) on the Ion Community.
Protocols: Visit the [Protocols page](#) on the Ion Community.