

File :Y:\Forensics\15937-0457042-6.D
Operator : PDK
Acquired : 14 Apr 2014 18:22 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: AKB48 4-fluorobenzyl analog
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 9

