

File :Y:\Forensics\15489-0455089-2.D
Operator : PDK
Acquired : 24 Mar 2014 16:00 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: 5-fluoro AMB
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 17

