

File :Y:\Forensics\14038-0449722.D
Operator : PDK
Acquired : 19 Aug 2013 11:26 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: AB-PINACA
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 1

