

File :Y:\Forensics\14293-0448608.D
Operator : PDK
Acquired : 10 Sep 2013 16:41 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: ADBICA
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 8

