

File :Y:\Forensics\15434-0455353-2.D
Operator : PDK
Acquired : 24 Mar 2014 11:27 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: AB-CHMINACA
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 1

