

File :Y:\Forensics\15533-0455794.D
Operator : PDK
Acquired : 12 Apr 2014 3:39 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: EG-018
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 11

