

File :Y:\Forensics\15064-0457195.D
Operator : PDK
Acquired : 29 May 2014 16:42 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: 4-methoxy PV9
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 5

