

File :Y:\Forensics\10913-0438977.D
Operator : PDK
Acquired : 5 Sep 2012 17:23 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: (±)-CP 47,497
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 5

