

File :Y:\Forensics\10010117-0420042-4.D
Operator : PDK
Acquired : 20 Sep 2012 17:16 using AcqMethod CAY_DRUG_CB.M
Instrument : Instrument #1
Sample Name: CB-52
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 120C-50C/min-280C
Vial Number: 19

