

File :Y:\Forensics\9001069-0446419.D
Operator : PDK
Acquired : 30 Apr 2013 15:21 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: 4-Methylethcathinone
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 11

