

File :Y:\Forensics\14106-0445071-2.D
Operator : PDK
Acquired : 21 Mar 2013 13:39 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: 3,4-dimethoxy-a-Pyrrolidinopentiophenone
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 9

