

File :Y:\Forensics\11864-0456525.D
Operator : PDK
Acquired : 15 Apr 2014 14:07 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: 4-hydroxy DMT
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 3

