

File :Y:\Forensics\11946-0443535.D
Operator : PDK
Acquired : 19 Dec 2012 13:55 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: 25D-NBOMe
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 4

