

File :Y:\Forensics\10826-0448904.D
Operator : PDK
Acquired : 25 Jun 2013 22:20 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: JWH 147
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 14

