

File :Y:\Forensics\14528-0455137-2.D
Operator : PDK
Acquired : 24 Mar 2014 20:44 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: 30C-NBOMe
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 3

