

File :Y:\Forensics\14889-0451662.D
Operator : LJJ
Acquired : 13 Nov 2013 21:10 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: THJ
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 12

