

File :C:\ISO\Data\11AUG2015\17283-0471433.D
Operator : LauraN
Acquired : 11 Aug 2015 14:44 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: ALÉPH-4
Misc Info : 30m x 0.32mm, Rtx-5MS, 1min@50C,30C/min to 300C
Vial Number: 4

