

File :C:\ISO\Data\09OCT2015\9001543-0473746-R.D
Operator : LIMS import
Acquired : 9 Oct 2015 17:03 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: Heroin RM
Misc Info : 30m x 0.32mm, Rtx-5MS, 1min@50C,30C/min to 300C
Vial Number: 5

