

File :C:\ISO\Data\17MAY2016\12036-0482560.D
Operator : LauraN
Acquired : 17 May 2016 15:20 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: AH-7921
Misc Info : 30m x 0.32mm, Rtx-5MS, 1min@50C,30C/min to 300C
Vial Number: 15

