

File :C:\ISO\Data\01MAR2016 cust inq\14039-0451007.D  
Operator : LauraN  
Acquired : 1 Mar 2016 11:41 using AcqMethod CAY\_DRUG\_CATH.M  
Instrument : ISO\_GCMS1  
Sample Name: AB-FUBINACA  
Misc Info : 30m x 0.32mm, Rtx-5MS, 1min@50C,30C/min to 300C  
Vial Number: 5

