

File :C:\ISO\Data\11MAR2015\9002195-0466680.D

Operator : Laura

Acquired : 11 Mar 2015 12:11 using AcqMethod CAY\_DRUG\_CATH.M

Instrument : ISO\_GCMS1

Sample Name: 9002195-0466680

Misc Info :

Vial Number: 18

