

File : C:\ISO\Data\06APR2015\17614-0467677.D
Operator : Laura N
Acquired : 6 Apr 2015 15:42 using AcqMethod CAY_DRUG.M
Instrument : ISO_GCMS1
Sample Name: APP Chminaca
Misc Info :
Vial Number: 8

