

File :C:\ISO\Data\06APR2015\17206-0467725.D
Operator : Laura N
Acquired : 7 Apr 2015 12:43 using AcqMethod CAY_DRUG.M
Instrument : ISO_GCMS1
Sample Name: 5-fluoro Cumyl pinaca
Misc Info :
Vial Number: 15

