

File :C:\ISO\Data\06APR2015\17210-0467623.D
Operator : Laura N
Acquired : 6 Apr 2015 16:16 using AcqMethod CAY_DRUG.M
Instrument : ISO_GCMS1
Sample Name: Cumyl Thpinaca
Misc Info :
Vial Number: 9

