

File :Z:\16434-0461309.D
Operator : Tom B
Acquired : 19 Sep 2014 14:46 using AcqMethod CAY_DRUG.M
Instrument : ISO_GCMS1
Sample Name: PX 2
Misc Info :
Vial Number: 10

