

File :Z:\16201-0460016-2.D
Operator : Tom B
Acquired : 15 Sep 2014 12:05 using AcqMethod CAY_DRUG.M
Instrument : ISO_GCMS1
Sample Name: PX1
Misc Info :
Vial Number: 14

