

File :Z:\9001950-0458839-2.D
Operator : Tom B
Acquired : 11 Sep 2014 16:38 using AcqMethod CAY_DRUG.M
Instrument : ISO_GCMS1
Sample Name: PTI-2
Misc Info :
Vial Number: 2

