

File :Z:\9001948-0458842-2.D
Operator : Tom B
Acquired : 24 Sep 2014 18:04 using AcqMethod CAY_DRUG.M
Instrument : ISO_GCMS1
Sample Name: PTI-1
Misc Info :
Vial Number: 1

