

File :Z:\15155-0459249-3.D
Operator : Tom B
Acquired : 23 Sep 2014 16:14 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: FAB-144
Misc Info :
Vial Number: 7

