

File :Z:\15969-0458217-8.D
Operator : Tom B
Acquired : 18 Sep 2014 11:57 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: 25C-NBF
Misc Info :
Vial Number: 2

