

File :Z:\15967-0458296-2.D
Operator : Tom B
Acquired : 19 Sep 2014 16:59 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: 25B-NBF
Misc Info :
Vial Number: 15

