

File :Y:\9002102-0461817.D
Operator :TomB
Acquired : 17 Jun 2015 18:02 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: 4-EAPB
Misc Info :
Vial Number: 19

