

File :Z:\15518-0459009.D
Operator : Tom B
Acquired : 12 Sep 2014 17:17 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: Diphenidine
Misc Info :
Vial Number: 4

