

File :Z:\9001933-0457892-2.D
Operator : Tom B
Acquired : 12 Sep 2014 13:15 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: N,N-Dimethylpentylone
Misc Info :
Vial Number: 10

