

File :Y:\17332-0465398-2.D
Operator : TomB
Acquired : 22 May 2015 17:21 using AcqMethod CAY_DRUG.M
Instrument : ISO_GCMS1
Sample Name: FUB-JWH 018
Misc Info :
Vial Number: 19

