

File :Z:\16436-0461224-3.D
Operator : Tom B
Acquired : 18 Sep 2014 18:56 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: 4-Chloromethcathinone
Misc Info :
Vial Number: 4

