

File :Z:\16200-0459976-7.D
Operator : Tom B
Acquired : 12 Sep 2014 18:06 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: MDMB-CHMINACA
Misc Info :
Vial Number: 6

