

File :C:\ISO\Data\23MAR2015\17394-0467004.D
Operator : Laura
Acquired : 23 Mar 2015 15:10 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: 3-Chloromethcathinone HCl
Misc Info :
Vial Number: 4

