

File :C:\ISO\Data\13JAN2016\16965-0471309.D

Operator : LauraN

Acquired : 14 Jan 2016 10:10 using AcqMethod CAY\_DRUG\_CATH.M

Instrument : ISO\_GCMS1

Sample Name: MDMA-CHMICA

Misc Info : 30m x 0.32mm, Rtx-5MS, 1min@50C,30C/min to 300C

Vial Number: 4

