

File :C:\ISO\03OCT2016\16966-0491071.D
Operator : SteveA
Acquired : 3 Oct 2016 18:06 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: MDMA-FUBINACA
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 4

