

File :C:\ISO\Data\09JUN2015\17299-0468145.D
Operator : LauraN
Acquired : 9 Jun 2015 18:50 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: 5-methoxy EiPT
Misc Info : 30mx0.32mm,0.5uRtx-5MS,1min@50C-30C/min to 300C
Vial Number: 12

