

File :C:\ISO\Data\30MAR2016\18596-0479029.D
Operator : LauraN
Acquired : 30 Mar 2016 12:35 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: U-47700
Misc Info : 30m x 0.32mm, Rtx-5MS, 1min@50C,30C/min to 300C
Vial Number: 5

