

File :C:\ISO\30JAN2017\21173-0497793.D
Operator : LauraN
Acquired : 30 Jan 2017 16:09 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: Nandrolone
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 5

