

File : C:\ISO\05DEC2018\22665-0545063.D
Operator : JianmeiL
Acquired : 5 Dec 2018 18:30 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: Bromazolam
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 15

