

File : C:\ISO\19JUN2019\23694-0550838-5.D  
Operator : JianmeiL  
Acquired : 19 Jun 2019 12:35 using AcqMethod CAY\_DRUG\_CATH.M  
Instrument : ISO\_GCMS1  
Sample Name: Vortioxetine (hydrobromide)  
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C  
Vial Number: 12

