

File :C:\ISO\22AUG2018\20859-0537082.D

Operator : JianmeiL

Acquired : 22 Aug 2018 11:52 using AcqMethod CAY_DRUG_CATH.M

Instrument : ISO_GCMS1

Sample Name: Tetrahydrofuran fentanyl HCl

Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C

Vial Number: 7

