

File :C:\ISO\07AUG2018 skb\18705-0537068.D
Operator : SBouhall
Acquired : 7 Aug 2018 17:02 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: Furanyl fentanyl (hydrochloride)
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 7

