

File :C:\ISO\15JUL2017\19410-0508010r.D
Operator : JianmeiL
Acquired : 15 Jul 2017 20:26 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: Carfentanil
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 3

