

File :C:\ISO\28NOV2017\9002414-0517879R.D
Operator : JianmeiL
Acquired : 29 Nov 2017 11:24 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: 4-fluoro IPV (HCl)
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 4

