

File :C:\ISO\03OCT2017\BSB\16616-0515924.D
Operator :JianmeiL
Acquired : 3 Oct 2017 20:30 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: MAB-CHMINACA
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 6

