

File :C:\ISO\30JUN2017\15480-0508404.D
Operator : JianmeiL
Acquired : 30 Jun 2017 15:46 using AcqMethod CAY_DRUG_300DEGHOLD.M
Instrument : ISO_GCMS1
Sample Name: W-18
Misc Info : 30m x 0.32mm, Rtx5MS, 300 degC hold 40 minutes
Vial Number: 6

