

File :C:\ISO\12OCT2016\18365-0492127.D
Operator : LauraLS
Acquired : 12 Oct 2016 12:07 using AcqMethod CAY_DRUG.M
Instrument : ISO_GCMS1
Sample Name: ATHPINACA isomer 1
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@240C,30C/min to 300C
Vial Number: 4

