

File :C:\ISO\Data\29OCT2015\18173-0473513-GCMS.D
Operator : NathanL
Acquired : 29 Oct 2015 21:36 using AcqMethod CAY_DRUG.M
Instrument : ISO_GCMS1
Sample Name: Clonazepam
Misc Info : 30m x 0.32mm, Rtx-5MS, 1min@240C,30C/min to 300C
Vial Number: 7

