

File :C:\ISO\Data\09OCT2015\15963-0469770.D
Operator : LIMS import
Acquired : 9 Oct 2015 17:22 using AcqMethod CAY_DRUG.M
Instrument : ISO_GCMS1
Sample Name: 5-fluoro PCN RM
Misc Info : 30m x 0.32mm, Rtx-5MS, 1min@240C,30C/min to 300C
Vial Number: 11

