

File :C:\ISO\Data\11NOV2015\17211-0465319_stune.D
Operator : LIMS import
Acquired : 18 Nov 2015 12:53 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: 17211-0465319
Misc Info : 30m x 0.32mm, Rtx-5MS, 1min@50C,30C/min to 300C
Vial Number: 5

