

File :C:\ISO\27JUN2017\15609-0507371.D
Operator : JianmeiL
Acquired : 27 Jun 2017 15:35 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: 4-Methyl-alpha-ethylaminopentiophenone (HCl)
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 8

