

File :C:\ISO\21NOV2017-non ISO\18030-0517875.D
Operator : JianmeiL
Acquired : 22 Nov 2017 10:42 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: 3',4' trimethylene-alpha-pyrrolidinohexanophenone (hydrochloride)
Misc Info : 30m x 0.32mm, Rtx5MS, 1min@50C,30C/min to 300C
Vial Number: 3

