

File :C:\ISO\02NOV2016\9001934-0491448.D
Operator : LauraN
Acquired : 2 Nov 2016 16:07 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: alpha-Pyrrolidinohexanophenone (hydrochloride)
Misc Info : 30m x 0.32mm, Rtx-5MS, 1min@50C,30C/min to 300C
Vial Number: 5

