

File : C:\ISO\Data\08MAR2016\18326-0476833.D
Operator : LauraN
Acquired : 8 Mar 2016 18:27 using AcqMethod CAY_DRUG_CATH.M
Instrument : ISO_GCMS1
Sample Name: 4'-chloro-alpha-Pyrrolidinovalerophenone (hydrochloride)
Misc Info : 30m x 0.32mm, Rtx-5MS, 1min@50C,30C/min to 300C
Vial Number: 12

