

File :Y:\Forensics\15166-0453912.D
Operator : LauraN
Acquired : 30 Jan 2014 14:21 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: 4-fluoro-a-Pyrrolidinopentiophenone□hydrochlo
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 14

