

File :Y:\Forensics\11113-0440571.D
Operator : LJJ
Acquired : 14 Nov 2013 15:34 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: 2C-I (hydrochloride)
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 2

