

File :Y:\Forensics\12065-0440921.D
Operator : PDK
Acquired : 17 Sep 2012 16:53 using AcqMethod CAY_DRUG.M
Instrument : Instrument #1
Sample Name: AKB48 N-(5-fluoropentyl) analog
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 240C-30C/min-300C
Vial Number: 12

