

File :Y:\Forensics\15388-0454860.D
Operator : PDK
Acquired : 23 Mar 2014 15:34 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: Mepirapim
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 5

