

File :Y:\Forensics\15971-0457890-2.D
Operator : PDK
Acquired : 30 May 2014 17:33 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: MMB2201
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 3

