

File :Y:\Forensics\11928-0442099.D
Operator : PDK
Acquired : 14 Nov 2012 13:36 using AcqMethod CAY_DRUG_CATH.M
Instrument : Instrument #1
Sample Name: UR-144 Degradant
Misc Info : 30mx0.32mm, 0.5u Rtx-5MS, 50C-30C/min-300C
Vial Number: 2

